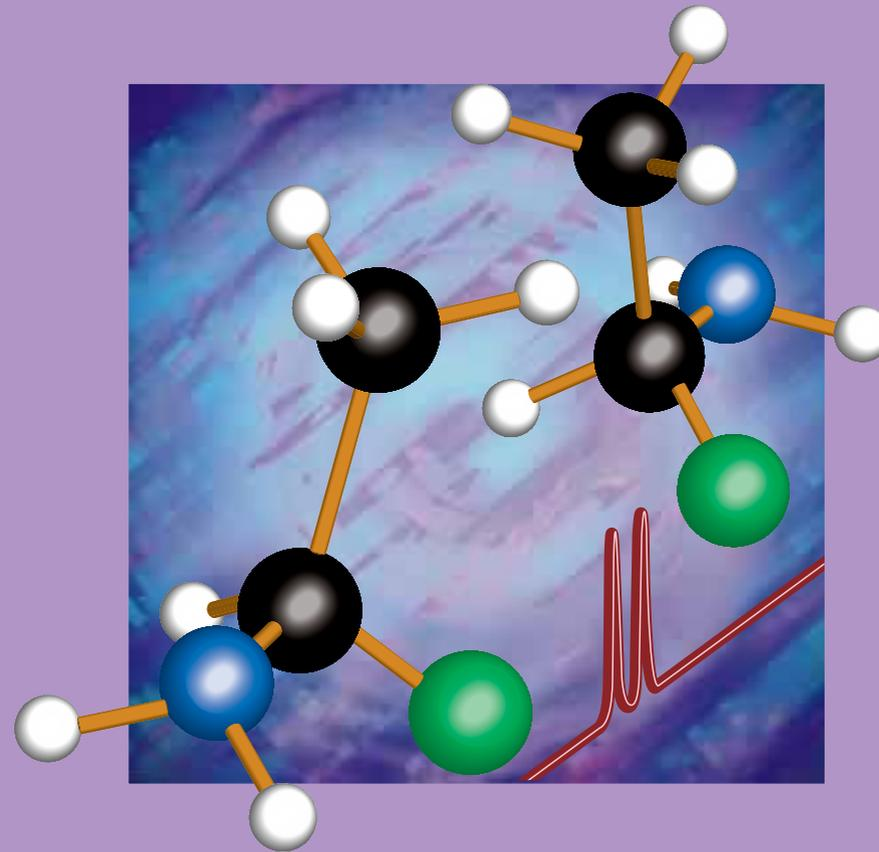


An Introduction to Chiral Analysis by Capillary Electrophoresis



BIO-RAD

An Introduction to Chiral Analysis by Capillary Electrophoresis

Salvatore Fanali

**Istituto di Cromatografia del Consiglio Nazionale
delle Ricerche
Area della Ricerche di Roma
P.O. Box 10 00016 Monterotondo Scalo (Roma) Italy**

Table of Contents

| | |
|---|-----------|
| About the Author | ii |
| Section 1 Introduction..... | 1 |
| 1.1 Abbreviations Used in the Text..... | 1 |
| 1.2 Introduction to Chiral Analysis | 2 |
| Section 2 Chirality and Optical Isomers Separation | 3 |
| Section 3 Separation of Diastereoisomers by Indirect Methods | 4 |
| Section 4 Direct Methods for Chiral Separation in Capillary Electrophoresis | 8 |
| 4.1 Inclusion Complexation | 12 |
| <i>Cyclodextrins and their Derivatives in Capillary Zone Electrophoresis</i> | <i>12</i> |
| <i>Cyclodextrins and their Derivatives in Capillary Gel Electrophoresis, Isotachopheresis, and Electrochromatography.....</i> | <i>25</i> |
| <i>Chiral Crown-Ether</i> | <i>26</i> |
| 4.2 Ligand Exchange | 27 |
| 4.3 Chiral Micelles..... | 28 |
| 4.4 Affinity Interactions..... | 30 |
| <i>Proteins as Chiral Selectors</i> | <i>30</i> |
| <i>Interactions with Saccharides (Linear Polymers) and Antibiotics</i> | <i>33</i> |
| 4.5 Combination of Chiral/Chiral or Chiral/Achiral Selectors ... | 33 |
| Section 5 Optimization of the Method..... | 35 |
| Section 6 Quantitation | 39 |
| Section 7 Applications..... | 40 |
| Section 8 Conclusions | 42 |
| Section 9 References..... | 43 |
| Acknowledgments | 50 |

About the Author

Dr. Salvatore Fanali is a researcher at the Istituto di Cromatografia del Consiglio Nazionale delle Ricerche in Montelibretti (Rome) Italy. After experience in high performance paper electrophoresis and paper chromatography, his research concerned the study of capillary isotachopheresis in the pharmaceutical field. Since 1988 his research has focused on electromigration methods, mainly applied to the separation of enantiomers. He is author or co-author of seventy publications, two booklets, and chapters in two books. He is a member of the advisory editorial boards of *Bollettino Chimico Farmaceutico*, *Chromatographia*, *Electrophoresis*, *Journal of Capillary Electrophoresis*, *Journal of Chromatography*, and *Journal of High Resolution Chromatography*.

Section 1 Introduction

1.1 Abbreviations Used in the Text

| | |
|----------------------|--|
| APA | aryl propionic acid |
| BGE | background electrolyte |
| CD | cyclodextrin |
| CGE | capillary gel electrophoresis |
| CM- β -CD | carboxymethylated- β -cyclodextrin |
| CM- β -CD pol | carboxymethylated- β -cyclodextrin polymer |
| CMC | critical micellar concentration |
| Crown-ether | 18-crown-6-ether tetra-carboxylic acid |
| CSP | chiral stationary phase |
| CZE | capillary zone electrophoresis |
| DAT | (+)-O,O'-diacetyl-L-tartaric anhydride |
| DBT | (+)-O,O'-dibenzoyl-L-tartaric anhydride |
| di-OMe- β -CD | 2,6-di-O-methyl- β -cyclodextrin |
| EHC | electro-chromatography |
| eof | electro-osmotic flow |
| GC | gas chromatography |
| HEC | hydroxyethylcellulose |
| HP- β -CD | hydroxypropyl- β -cyclodextrin |
| HPLC | high performance liquid chromatography |
| HPMC | hydroxypropylmethylcellulose |
| ITP | isotachopheresis |
| LE | leading electrolyte |
| LEC | ligand exchange chromatography |
| MEKC | micellar electrokinetic chromatography |
| NDA | naphthalene-2,3-dicarboxyaldehyde |
| NSAIDs | non-steroidal anti-inflammatory drugs |
| PhL | phenyl lactic acid |
| PVA | polyvinyl alcohol |
| PVP | polyvinyl pyrrolidone |
| SBE- β -CD | sulfobutyl-ether(IV)- β -cyclodextrin |
| SDAla | sodium N-dodecanoyl-L-alaninate |
| SDS | sodium dodecyl sulfate |
| SDVal | sodium N-dodecanoyl-L-valinate |
| STC | sodium taurocholate |
| STDC | sodium taurodeoxycholate |
| TLC | thin layer chromatography |
| tri-OMe- β -CD | 2,3,6-tri-O-methyl- β -cyclodextrin |

1.2 Introduction to Chiral Analysis

The separation of enantiomers is widely studied in analytical chemistry, especially in the pharmaceutical and biological fields, since chiral drugs are administered either as enantiomers or as racemic mixtures. Very often two enantiomers of the same racemic drug possess different pharmacological effects. For example, S(-)-propranolol is considerably more active than its enantiomers. The anesthetic ketamine is administered as racemate, and the S(+)-ketamine form is more potent than the R(-) form. In addition, the R(-) form may cause post-operative effects. Because of the side effects that could be caused by the presence of an undesirable component in a racemic drug, the current tendency in the pharmaceutical industry is to prepare drugs with one enantiomer only. However, the production of such drugs through stereoselective reaction or preparative enantiomeric separation can provide impure material. Thus rapid, sensitive analytical methods of high resolving power are required to control the synthetic chiral process of the pharmaceutical.

Analytical methods which have been used for the separation of chiral compounds include high performance liquid chromatography (HPLC),¹⁻³ gas chromatography (GC),^{4,5} thin layer chromatography (TLC),⁶ and recently, capillary electrophoresis (CE).⁷⁻¹⁵ CE is especially useful for the analysis of different classes of compounds, including organic and inorganic ions, peptides, proteins, saccharides, drugs, optical isomers, and others. In CE the separation will take place if the analytes, under the influence of an applied electric field, move toward the detector with different velocities (different effective mobilities). Thin capillaries, with internal diameters of 10–100 μm , allow application of relatively high electric fields (200–1,000 V/cm) permitting analytes to be separated in short times with high resolution and high efficiency, because the Joule heat, the main cause of zone dispersion, is easily controlled.

The different separation modes available in CE, namely zone electrophoresis (CZE), micellar electrokinetic chromatography (MEKC), gel electrophoresis (CGE), and isotachopheresis (ITP), combined with the high resolving power and high efficiency, make the CE technique competitive with others.¹⁶ CE offers additional advantages over other techniques such as HPLC, including

- A relatively small volume of sample and buffer (nl and μl , respectively) is required
- Expensive chiral columns can be avoided because the chiral selector can be easily added to the BGE

- The separation is highly reproducible because the buffer with the chiral selector is replenished after each run

The use of CE for analysis of chiral compounds is well documented and, considering the enormous increase in publications, it can be expected that in the near future chiral analysis by CE will be widely applied to pharmaceutical, clinical, and environmental samples.

Published reviews,¹⁷⁻²⁴ and the special issue edited by us²⁵ dealing with the separation of optical isomers by CE, are available for further reading. The article by Fanali *et al.* provides a list of enantiomeric compounds which have been separated using capillary electrophoresis.²⁴

Section 2 Chirality and Optical Isomers Separation

When four different ligands are bound to a tetravalent carbon, an asymmetric molecule is generated in which the carbon is the asymmetric center. As shown in Figure 1, two optical isomers can be generated due to the different spatial orientations of the ligands around the chiral center.

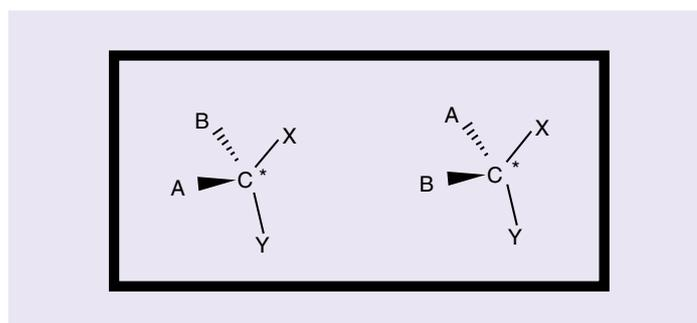


Fig. 1. Mirror image of two optical isomers.

Enantiomers are two stereoisomers which exhibit non-superimposable mirror images. Diastereomers generally possess at least two asymmetric centers (one of them has the same configuration), and are not mirror images. The most common chiral center is represented by tetrahedral carbon, although other atoms, like nitrogen, sulfur, and phosphorous, can be found in stereoisomers. Compounds possessing at least two enantiomers are chiral compounds.

The main property of stereoisomers is represented by the rotation of polarized light in different directions, counter-clockwise (levo-rotatory) and clockwise (dextro-rotatory). They are l(-)- and d(+)-isomers, respectively. The optical rotation of two enantiomers (in a racemic mixture) is of the same magnitude, while in the case of diastereomers the two parameters may not be the same due to different physico-chemical properties. Unlike enantiomers, diastereomers may have different melting points, boiling points, and solubility.

Although the d, or (+), and l, or (-), symbols show a very important physical property of the molecule, they unfortunately do not give any information concerning the spatial arrangement of the chiral center. The Fisher convention, widely employed for sugars and amino acids, makes use of D and L symbols and is based on the comparison of the substituents of the chiral center of the compound under investigation with that of (+)-glyceraldehyde. However, the D and L convention can be confused with d and l terminology, and thus the currently recognized convention is the Cahn-Ingold-Prelog.²⁶ Here the priority of the ligand to the asymmetric center based on the atomic number is controlled, and the group with lower priority is positioned far away from the observation point. Then the priority of the other substituents is examined. If it decreases clockwise, the R configuration is assigned, otherwise the S configuration is given.

Section 3

Separation of Diastereoisomers by Indirect Methods

Because two diastereoisomers of the same compound have different physico-chemical properties, they can, in principle, be separated from each other by CE using a non-chiral electrophoretic system. As an example, consider the chemical structure of nor-ephedrine and ψ -nor-ephedrine (Figure 2). The two compounds possess the same molecular weight and the same substituent groups at the two asymmetric carbons, and the configuration of only one of the asymmetric centers is the same; thus they are diastereoisomers. Each diastereoisomer exists as a pair of enantiomers (d- and l-norephedrine and d- and l- ψ -norephedrine, respectively).

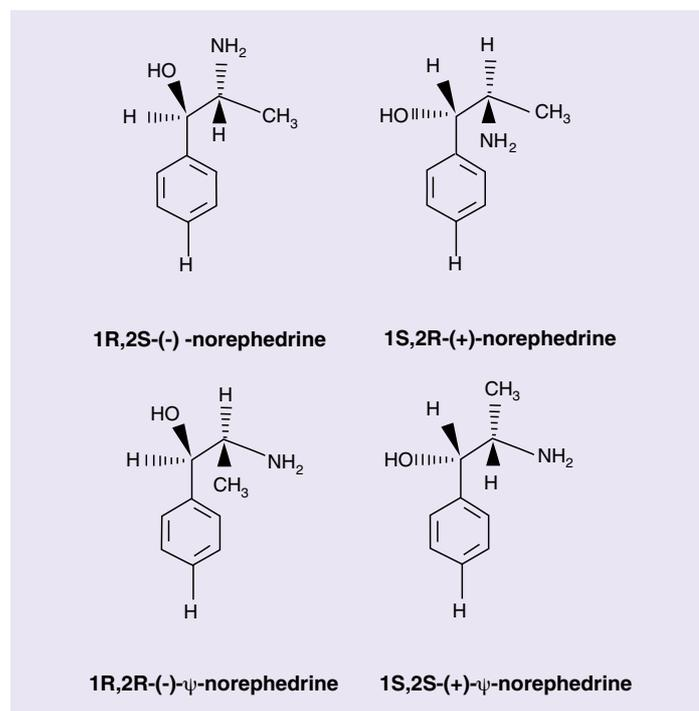


Fig. 2. Chemical structure of nor-ephedrine and ψ -nor-ephedrine.

When a mixture of the two compounds is injected for an electrophoretic run using a non-chiral background electrolyte (phosphate buffer at pH 2.5, see Figure 3), two peaks, corresponding to the two diastereoisomers, are obtained; each peak represents a couple of enantiomers, but the electrophoretic system is unable to separate them.

The above principle can be advantageously used for the resolution of a pair of enantiomers. Thus it is necessary to use a chiral environment that will modify the physico-chemical properties of the two analytes, transforming enantiomers into diastereoisomers. By adjusting the selectivity of the separation, the resolution can be improved.

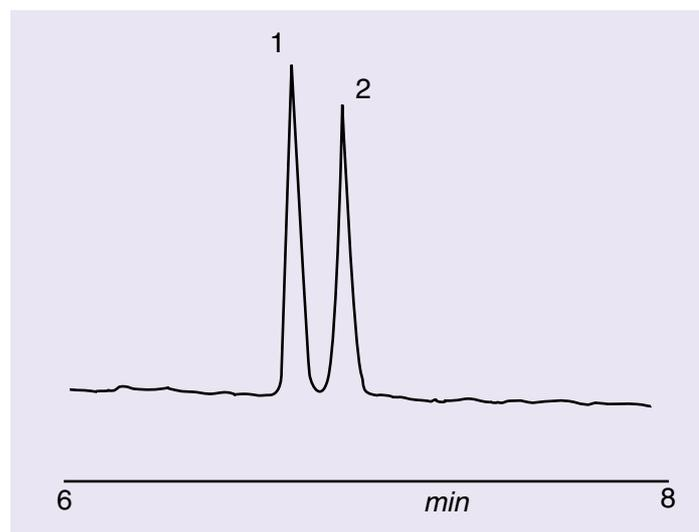


Fig. 3. Electropherogram of the separation of 1) ψ -norephedrine and 2) norephedrine. Apparatus, BioFocus[®] 3000 system; capillary, (coated) 35 cm x 0.05 mm ID; background electrolyte, 50 mM phosphate buffer, pH 2.5; applied voltage 15 kV; injection, electrokinetic 7 kV, 7 s of 2×10^{-6} M of each compound; carousel and capillary temperature 25 $^{\circ}$ C; detection at 206 nm.

The indirect separation method is based on the reaction, before the analysis, of a racemic mixture with a chiral reagent (R or S), producing a mixture of two diastereoisomers that can be resolved using a non-chiral electrophoretic system. The product of the reaction is a mixture of two stable compounds where relatively strong bonds (covalent) are involved in the process.

The derivatization of α -amino- or α -hydroxyacids with (+)-O,O'-diacetyl-L-tartaric anhydride (DAT) or (+)-O,O'-dibenzoyl-L-tartaric anhydride (DBT) can serve as an example. Figure 4 shows the chemical structure of the two diastereoisomers formed between DBT and L- and D-amino acids. The configuration of the two carbons of the tartrate is the same, and that of the amino acid chain is different.

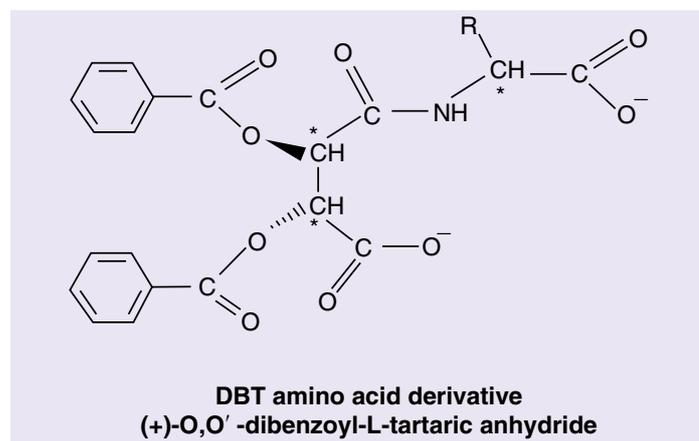


Fig. 4. Chemical structure of the two diastereoisomers formed in the reaction between (+)-O,O'-dibenzoyl-L-tartaric anhydride (DBT) and L-, D-amino acid.

The separation of such derivatives was obtained by CE using a BGE at pH 5.8 containing polyvinyl pyrrolidone (PVP) as a physical network.²⁷ The effect of the concentration of PVP added to the BGE for the separation of mandelic acid and tryptophan after derivatization with DBT²⁸ is shown in Figure 5.

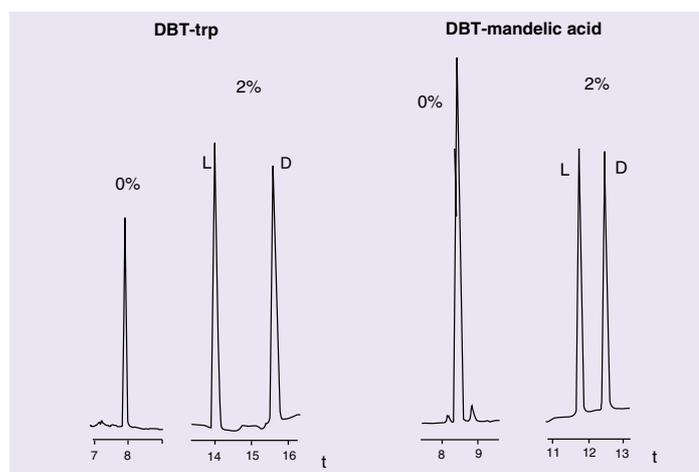


Fig. 5. Effect of the concentration of polyvinylpyrrolidone (PVP) on the separation of diastereoisomers of DPT-D,L-mandelic acid and DBT-D,L-tryptophan. Capillary, (coated) 56 cm x 0.1 mm ID; background electrolyte, 25 mM phosphate buffer, pH 5.8; applied voltage 12 kV; injection, hydrodynamic (5 s at a height distance of 10 cm) of 10^{-4} M of sample; detection at 233 nm (modified from reference ²⁸).

Although the indirect separation method using the appropriate chiral reagent (R or S) can be used advantageously when an inversion of migration order is required, the method presents several drawbacks. It is time consuming, it requires the use of very pure chiral reagent, the enantiomers must contain reacting groups (hydroxyl, amino), and the kinetics of the reaction can lead to different peak areas for the two diastereomers formed. Probably due to these and other drawbacks of the indirect separation method, the direct separation method is becoming very popular for enantiomeric separation by CE.

Section 4

Direct Method for Chiral Separation in Capillary Electrophoresis

In the direct separation method the chiral selector can be added to the BGE, bound to the capillary wall, or included in a gel matrix. It interacts with the two enantiomers during the electrophoretic process, forming labile diastereoisomeric complexes. Relatively weak bonds are involved, *e.g.*, hydrogen, π - π , or hydrophobic.

The separation of two enantiomers can take place only if the two diastereoisomers formed possess different stability constants, causing the two analytes to move with different velocities. The effective mobility of the most complexed enantiomer is lower than that of its isomer (this is not true if the chiral selector is negatively charged). For example, Figure 6 shows the separation of the four optical isomers analyzed in Figure 3, using the same background electrolyte supplemented with a chiral selector (2,6-di-O-methyl- β -CD) that allows the separation not only of the diastereoisomers but also of the enantiomers present in the mixture.

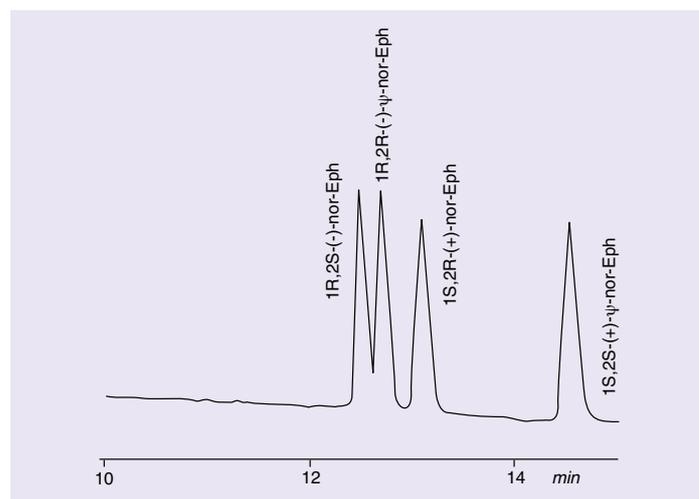


Fig. 6. Electropherogram of the separation of the four optical isomers of norephedrine and ψ -norephedrine. Apparatus, BioFocus 3000 system; background electrolyte, 50 mM phosphate buffer, pH 2.5, and 20 mM of 2,6-di-O-methyl- β -cyclodextrin. For other experimental conditions, see Figure 3.

The direct separation of enantiomers by CE is easier to perform than the indirect separation. This is demonstrated by the increasing number of publications reported in the literature (see references 24 and 29). The direct method is less time consuming since derivatization and purification are not required, a wide number of chiral selectors are commercially available, and small amounts of chiral selectors can be used. Furthermore the chiral purity of the system is not of paramount importance. In fact, the presence of impure chiral selectors will only cause a reduction of resolution of the two enantiomers in the direct method. In the indirect method, the presence of impure chiral selectors will produce four diastereoisomers (more difficult to separate). This is demonstrated in Figure 7, where a mixture of (+)-Co(en)₃³⁺ and (-)-Co(en)₃³⁺ is separated using sodium L(+)-tartrate; when the BGE contained the same chiral selector with 10% D(-)-tartrate, the resolution of the two enantiomers was reduced, while use of 100% D(-)-tartrate allowed baseline resolution with inversion of migration order.

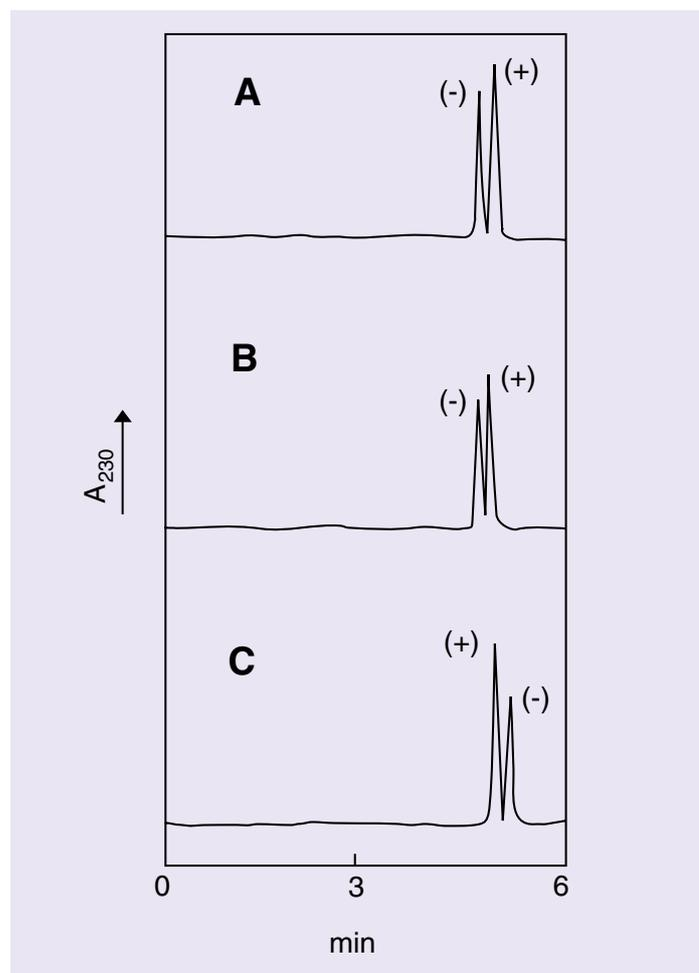


Fig. 7. Electropherograms of the separation of (+/-)-Co(en₃)³⁺. Apparatus, BioFocus 3000 system; capillary, (coated) 24 cm x 0.05 mm ID; background electrolyte, 50 mM sodium tartrate, pH 5.1 [A= 100% L(+), B= 90% L(+) and 10% D(-), C= 100% D(-)]; applied voltage 5 kV, 9.4 A; injection, 5 psi*s 35 mg racemic Co-complex + 9 mg (+)-antipode in 100 ml; detection at 230 nm.

Table 1 summarizes the main chiral selectors used in CE for the direct separation of enantiomers, as well as the different CE modes used and the main mechanisms involved in the separation process.

Table 1. Chiral Inclusion-Complexing Agents Used In Capillary Electrophoresis.

| Compound | CE Type | Mechanism | References |
|--|---------------------|-----------------|----------------|
| Uncharged Compounds | | | |
| α -cyclodextrin (α -CD) | CZE, ITP, CGE, MEKC | inclusion | 14, 30–32 |
| β -cyclodextrin (β -CD) | CZE, ITP, CGE, MEKC | inclusion | 15, 33–35 |
| γ -cyclodextrin (γ -CD) | CZE, ITP | inclusion | 14, 36–38 |
| glycosylated- α -cyclodextrin (G- α -CD) | CZE | inclusion | 39 |
| heptakis-2,6-di-O-methyl- β -cyclodextrin (di-OMe- β -CD) | CZE, ITP | inclusion | 13, 15, 30, 40 |
| heptakis-2,3,6-tri-O-methyl- β -cyclodextrin (tri-OMe- β -CD) | CZE, ECH | inclusion | 40–43 |
| hydroxyethyl- β -cyclodextrin (HE- β -CD) | CZE | inclusion | 41 |
| hydroxypropyl- β -cyclodextrin (HP- β -CD) | CZE | inclusion | 33, 39, 41, 44 |
| allyl derivatized- β -CD (Al- β -CD) | CGE | inclusion | 45 |
| Uncharged- β -cyclodextrin polymer | CZE | inclusion | 46 |
| Chargeable Compounds | | | |
| 6A-methylamino- β -cyclodextrin (6A-NH- β -CD) | CZE | inclusion | 47 |
| 6A,6D-dimethylamino- β -cyclodextrin (6A,D-di-NH- β -CD) | CZE | inclusion | 47 |
| 4-sulfobutyl-ether- β -cyclodextrin (SBE- β -CD) | CZE | inclusion | 33, 48–52 |
| mono-(6- β -aminoethylamino-6-deoxy)- β -cyclodextrin (β -CDen) | CZE | inclusion | 53 |
| carboxymethyl- β -cyclodextrin (CM- β -CD) | CZE | inclusion | 9, 54 |
| carboxymethyl- β -cyclodextrin polymer | CZE | inclusion | 55 |
| 18-crown-6-ether tetracarboxylic acid (18-crown ether) | CZE | inclusion | 32, 56 |
| copper(II)/L-proline or L-hydroxyproline or aspartame | CZE, MEKC | ligand exchange | 57–59 |
| maltodextrins | CZE | affinity | 60 |
| bile salts | MEKC | affinity-MEKC | 7, 61, 62 |
| proteins | CZE | affinity | 63–65 |
| antibiotics | CZE | inclusion | 66, 67 |
| L-tartrate | CZE | outer-sphere | 68 |

4.1 Inclusion Complexation

The resolution of enantiomers has been obtained in CE through inclusion-complexation mechanisms using either cyclodextrins or crown-ether derivatives. Here the two analytes (guest compounds) fit the cavity of the chiral selector (host compound), forming host-guest or inclusion-complexes. The separation of the two enantiomers can take place only if the two diastereomeric complexes, formed during the electrophoretic process, possess different stability constants. The chiral separation is obtained due to the formation of secondary bonds between the substituent groups on the chiral center of the analytes and those of the chiral selectors positioned outside the cavity (hydroxyl or modified hydroxyl and carboxylic groups for cyclodextrins and crown-ethers, respectively). To clarify the recognition mechanism of chiral resolution using inclusion-complexation, it is necessary to describe the structure and the main properties of the chiral selectors involved in the electrophoretic process.

Cyclodextrins and their Derivatives in Capillary Zone Electrophoresis

Cyclodextrins (CD) are neutral and natural cyclic oligosaccharides composed of several glucopyranose units. They are chiral due to the presence of asymmetric carbons on the glucose units. In spite of the fact that cyclodextrins containing 6–12 D(+)-glucopyranose units are reported in the literature, only 6, 7, and 8 unit-formed molecules, named with the Greek alphabet letters α , β , and γ , respectively, are in frequent use in analytical chemistry.^{69,70}

The shape of cyclodextrins is similar to that of a truncated cone with a cavity of different dimensions depending on the CD type (the number of the glucose units and the substituent groups). Their cavity is relatively hydrophobic and able to accept guest compounds of different types, particularly those with non-polar groups. The outside is relatively hydrophilic due to the presence of hydroxyl groups (primary and secondary).

Figure 8 shows the chemical structure of α -cyclodextrin, while Table 2 gives the main properties of the most commonly used CDs.

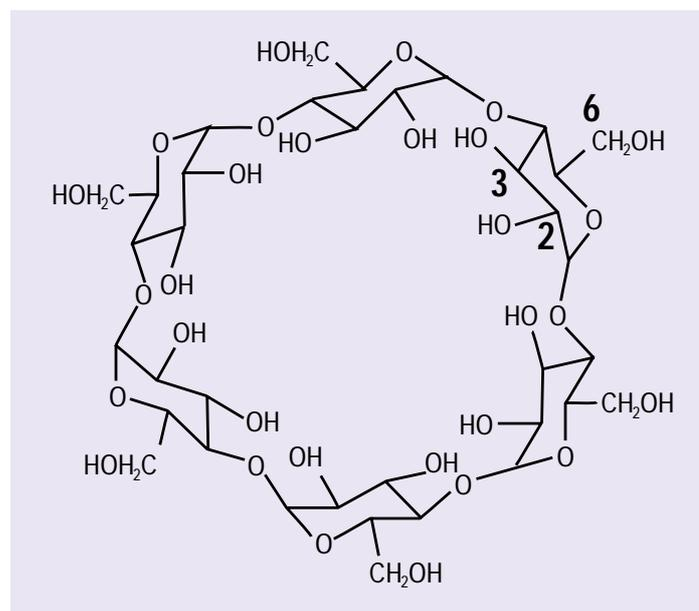


Fig. 8. Chemical structure of α -cyclodextrin.

Table 2. The Main Properties of Native Cyclodextrins ^{69,71}

| Cyclodextrin type | α | β | ψ |
|--|-----------|-----------|-----------|
| Number of glucopyranose | 6 | 7 | 8 |
| Molecular weight | 973 | 1,135 | 1,297 |
| Internal diameter (nm) | 0.47–0.52 | 0.60–0.64 | 0.75–0.83 |
| Depth (nm) | 0.79–0.80 | 0.79–0.80 | 0.79–0.80 |
| Specific rotation [α] _{25 D} | 150.5 | 162.5 | 177.4 |
| Melting point (K) | 551 | 572 | 540 |
| Solubility in water g/100 ml 25 °C | 14.50 | 1.85 | 23.20 |

The solubility of β -CD is relatively low when compared to that of α and γ ; this could be a problem when using the chiral selector as an additive to the background electrolyte. When higher concentrations of β -CD must be used, water-methanol, -ethanol, or -urea mixtures can be successfully used. In fact the solubility of β -CD in aqueous solutions of 4 and 8 M urea increases up to 0.089 and 0.226 M, respectively.⁷²

The hydroxyl groups present on the two entrances of CD at position 2, 3, and 6 of each glucopyranose can be modified, by chemical reactions, to produce CD derivatives with different properties than those of the parent ones. By manipulating the reaction type, the operational reaction conditions, and the ratio of the reagents, the modification can be obtained in only one or more than one hydroxyl group. Reaction with hydroxyl groups alters the properties of the CD, allowing

- improved solubility of the CD
- formation of different bonds with analytes that can improve the inclusion-complexation
- analysis of uncharged optical isomers

For example, the solubility of 2,6-di-O-methyl- β -CD and of the negatively charged β -cyclodextrin polymer in water were found to be 57 g/100 ml⁷³ and >40 g/100 ml, respectively, relatively high in comparison to that of β -CD (1.85 g/100 ml). Furthermore, the cavity of di-OMe- β -CD is deeper than its parent due to the presence of methoxy groups at the entrances, and thus is more hydrophobic. Considering other derivatives like carboxymethyl, methyl amino, phosphate, and sulfate, it is easy to recognize the advantages in analytical chemistry, especially in CE. In fact, such types of CDs can be charged to allow the optimization of the separation, *e.g.*, migration in the opposite direction of the analyte, or introduction of charge to neutral compounds through complexation (for a comprehensive review of CD derivatives, see reference 74).

Inclusion complexation and stereoselectivity are influenced by several experimental parameters, such as CD type and concentration, applied voltage, capillary temperature and length, ionic strength, organic solvent, electro-osmotic flow, and polymeric additives. Some effects of these parameters are discussed in this section, with the aim of describing the appropriate experimental conditions when enantiomers must be separated using cyclodextrins.

CD Type and Concentration

The first requirement for inclusion complexation is fitting the analyte into the CD cavity, and thus CD selection must be made considering the shape of the analyte. As previously discussed, the internal diameters of native CDs increase by increasing the number of glucose units $\gamma > \beta > \alpha$. Thus, selecting the appropriate CD is related to the shape and dimensions of the analyte. As an example,

consider the chemical structure of terbutaline and nicergoline (see Figure 9a and 9b). Terbutaline is currently used in the treatment of asthma diseases, and nicergoline is used as a vasodilator in acute myocardial infarction with diastolic hypertension. Terbutaline contains only one aromatic ring and one asymmetric carbon at the α position, while nicergoline is formed by more than one aromatic ring and possesses several chiral centers.

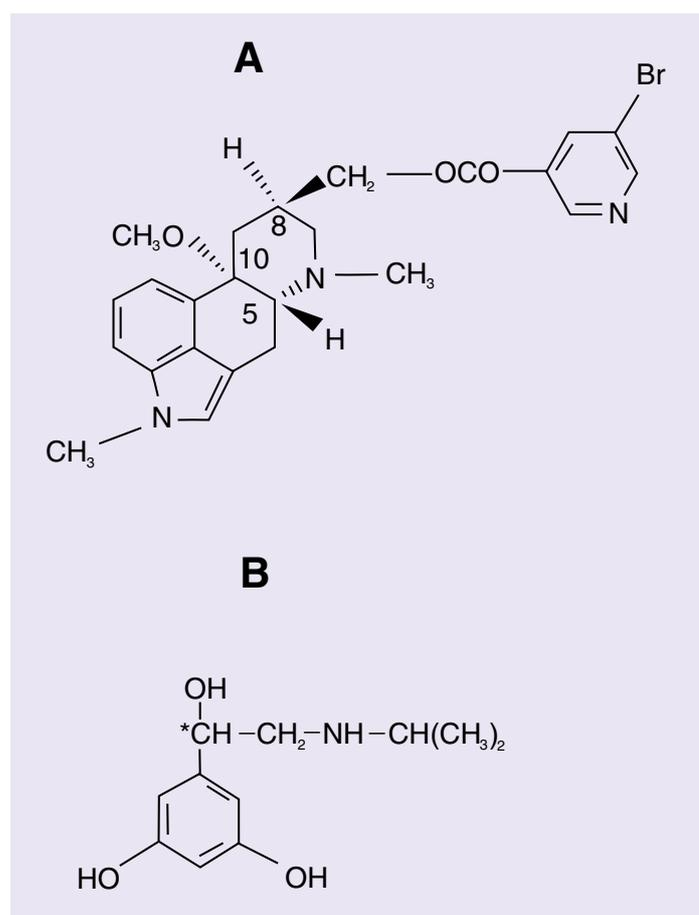


Fig. 9. Chemical structure of (A) nicergoline and (B) terbutaline.

Figures 10a, 10b, and 10c show the electropherograms of the separation of terbutaline and nicergoline into their enantiomers by capillary zone electrophoresis when α , β , and γ -CD were separately added to a BGE at pH 2.5.

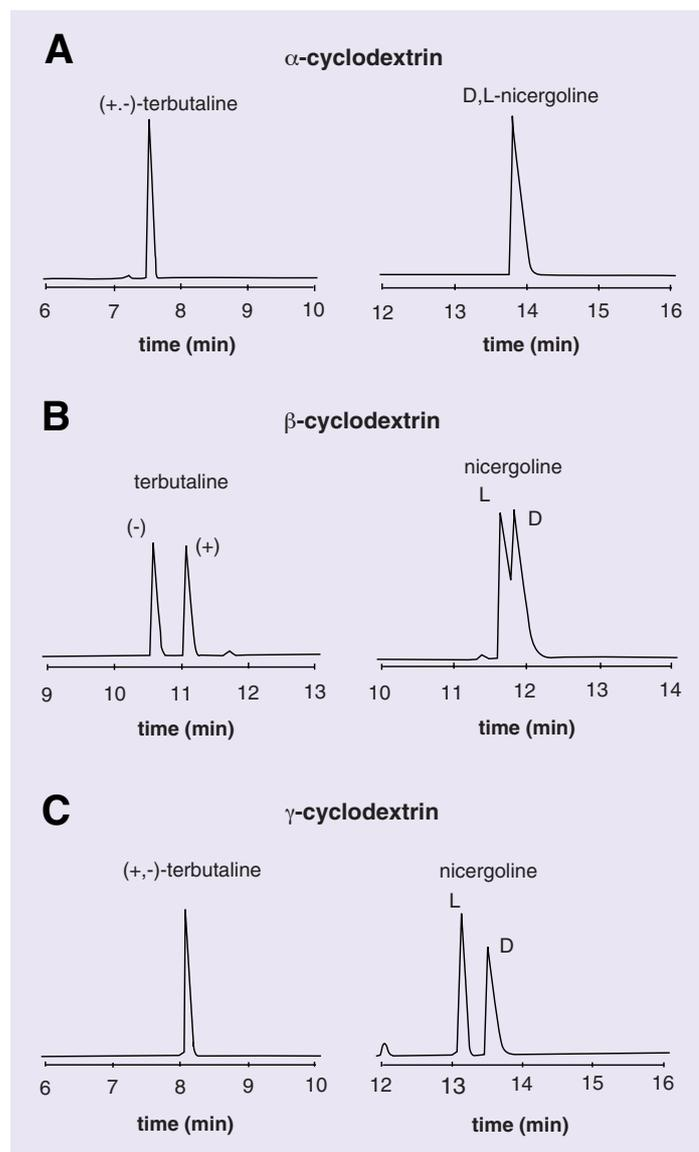


Fig. 10. Electropherograms of the enantiomeric separation of racemic terbutaline and nicergoline using different types of cyclodextrins as chiral selectors by CE. Apparatus, BioFocus 3000 system; capillary, (coated) 35 cm x 0.05 mm ID; background electrolyte, 0.1 M phosphate buffer +a) 30 mM α -CD, b) 20 mM β -CD and c) 30 mM γ -CD; applied voltage 15 kV; injection, electrokinetic 7 kV, 7 s of 5×10^{-5} M of racemic terbutaline or nicergoline.

Clearly α -CD was not able to separate both racemic mixtures into their enantiomers because its cavity is too small to guest the two compounds. β -CD allowed the resolution of terbutaline, while nicergoline was partly resolved, and γ -CD was not able to guest terbutaline (the molecule is too small) but was very effective for nicergoline.

In a previous study it has been shown that for the enantiomeric separation of terbutaline, either β -CD and dimethylated- β -CD can be successfully used in CE but the latter forms the strongest inclusion complexes and is more stereoselective than the former. The complexation of terbutaline increased with increasing concentration of both β -CD and its derivative, while the resolution showed a maximum at 15 mM and 5 mM for β -CD and dimethylated- β -CD, respectively.¹⁵ Figure 11 shows the effect of the concentration of β -CD and di-O-methyl- β -CD on the resolution of (+) and (-)-terbutaline.

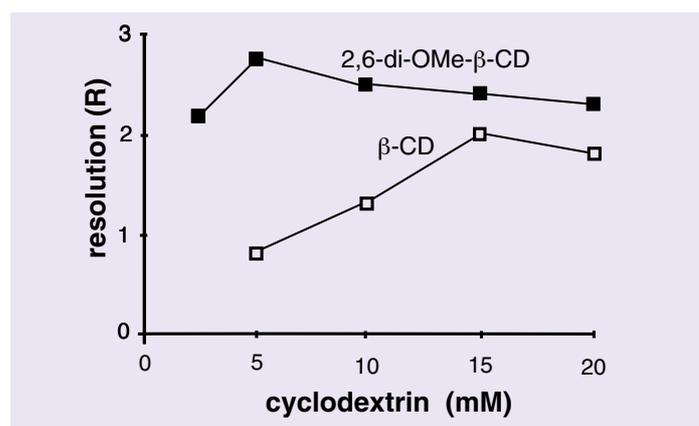


Fig. 11. Effect of the concentration of β -cyclodextrin and di-O-methyl- β -cyclodextrin on the resolution of racemic terbutaline into their enantiomers.

It is important to consider the degree of substitution (DS) of the cyclodextrin used. The effect of DS was demonstrated for the chiral separation of some hydroxyacids using different β -CDs modified with methyl amino⁴⁷ or hydroxypropyl groups⁷⁵ where the presence of more than one substituent can strongly affect the chiral recognition. The CD type can also influence the migration order of the two enantiomers; this effect was observed for the separation of dansyl-norvaline enantiomers using trimethylated- β -CD and β -CD.⁷⁶ Other authors have discussed the influence of the concentration of CD in CZE, indicating that the concentration of the CD can play a very important role for chiral resolution by CE, and thus this parameter should be

carefully controlled in order to find the optimal experimental conditions.^{8,34,38} Wren and Rowe discussed the optimum concentration of CD for enantiomeric separation and found that it is dependent upon the formation constants of the diastereoisomers formed during the electrophoresis (equation 1).⁷⁷

$$(1) \quad [C_{CD}] = \frac{1}{[K_{(+)} K_{(-)}]^{1/2}}$$

where C_{CD} is the optimum concentration of the CD used, and K is the formation constants of the two enantiomers (+) and (-). The mathematical model was verified for the enantiomeric separation of propranolol using β -CD or its dimethylated derivative.

Analyte shape

The shape of the analytes is another important parameter to be considered for inclusion-complexation with cyclodextrin. This effect was studied for the resolution of several tryptophan derivatives (methyl, hydroxyl) where the position of the substituent group on the indole ring strongly influenced the stability of the complex formed, and thus the chiral recognition.³⁰ The shape of the enantiomeric pairs can be modified, for example, by derivatizing with hydrophobic achiral compounds that can be included into the CD cavity. Amino acids as dansyl-, naphthalene-2,3-dicarboxaldehyde- (NDA), and naphthylamide derivatives, and monosaccharides derivatized as naphthalene sulfonate, were resolved into their enantiomers using cyclodextrins.⁷⁸⁻⁸¹

Charge of CD

The charge of the cyclodextrin can play an important role in the resolution mechanism when the electrophoretic separation of enantiomers has to be carried out. In this case the electrostatic interaction with the analytes, the movement of the chiral selector in the opposite direction of the two enantiomers, and the possibility for separating uncharged compounds (inducing the charge) are the main advantages for using such chiral selectors.

The use of charged cyclodextrins was first demonstrated by Terabe,⁵³ who separated several racemic dansyl-amino acids using a positively charged β -cyclodextrin. In our laboratory we used a methylamino- β -cyclodextrin for the enantiomeric resolution of several α -hydroxyacids⁴⁷ and a chargeable β -cyclodextrin polymer (carboxymethylated- β -CD polymer, CM- β -CD pol)⁵⁵ for the separation of racemic basic compounds into their enantiomers.

The inversion of migration order of studied enantiomers was demonstrated^{55,82} using a charged chiral selector in the presence of electro-osmotic flow; this approach can be advantageously employed when the enantiomeric impurity migrates behind the main isomer. In this case, it was necessary to increase the injected amount of the sample in order to detect the impurity (< 1%), and a loss of resolution occurred.

Electro-osmotic flow, concentration of the chargeable cyclodextrin, and the pH of the BGE have to be controlled to optimize the separation method.

When carboxymethylated- β -CD polymer is used in CE, the weakly acidic carboxylic groups can be charged or uncharged, depending upon the pH of the BGE. At low pH, 2–3, the carboxylic groups of the CD are protonated, the chiral selector behaves as a quasi stationary phase, and the groups form hydrogen bonds with analytes. At pH >3.5, the CD is charged due to the dissociation of the carboxylic substituents and migrates with its own mobility, while forming inclusion complexes and allowing ion-pair interactions with analytes.

Recently a new modified β -cyclodextrin (sulfobutyl-ether(IV)- β -cyclodextrin, SBE- β -CD) has been synthesized and characterized by CE with indirect UV detection.⁸³ This modified β -CD has been used in CE for the enantiomeric separation of several classes of compounds (positively charged, negatively charged, or uncharged). As shown in Figure 12, the SBE- β -CD contains four modified primary hydroxyl groups (position 6 of four glucose units) with a butyl chain and sulfonic groups. Due to its chemical properties, the modified CD is negatively charged at any commonly used pH in CE. These features allow its use in a charged mode over a wide pH (2–11) range.

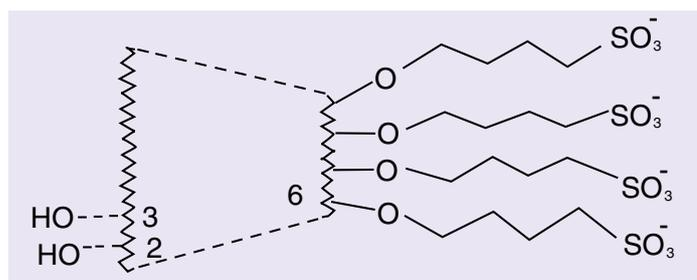


Fig. 12. Chemical structure of sulfobutyl-ether(IV)- β -cyclodextrin (SBE- β -CD).

The chiral selector was used at pH 10, close to the pK of several ephedrine, to obtain good enantiomeric resolution.⁴⁸ Operation at pH 2.5 allowed the enantiomeric separation of several basic compounds of pharmaceutical interest, including thalidomide, dimethindene, and mefloquine.⁴⁹ Recently we showed the usefulness of this new chiral selector for the enantiomeric separation of several positively and negatively charged compounds of pharmaceutical interest as well as uncharged phenyl-alcohols and dansyl-amino acids.⁵² In the case of acidic compounds, ion-pair interactions are not involved in the chiral resolution mechanism, but electrostatic repulsion can be responsible for the different stability constants of the two diastereoisomers formed during the electrophoretic process. Figure 13 shows the electropherograms of the enantiomeric separation of warfarin, bupivacaine, 2-phenyl-2-butanol and dansyl-phenylalanine.

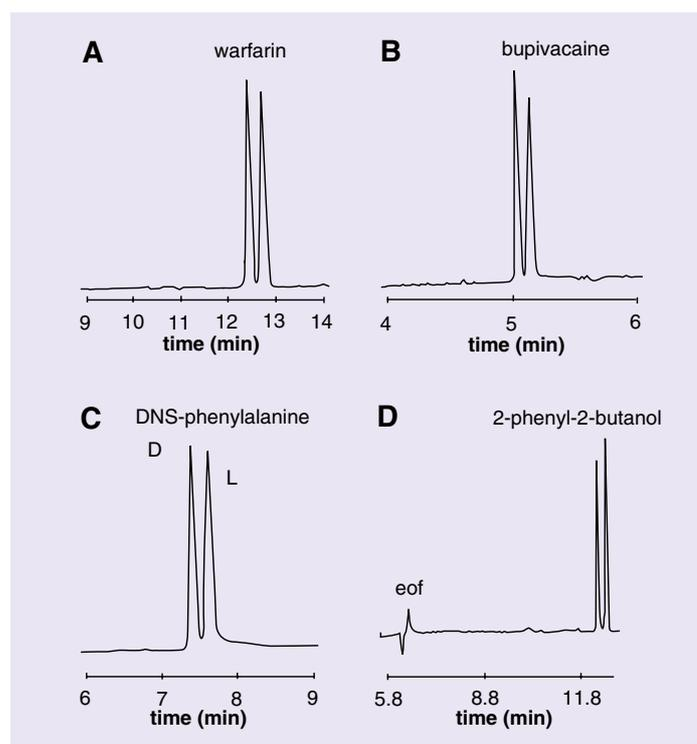


Fig. 13. Electropherograms of the enantiomeric separation of warfarin, bupivacaine, 2-dansyl-phenylalanine, and phenyl-2-butanol. Experimental conditions: apparatus, BioFocus 3000 system; capillary, 50 cm x 0.05 mm ID; applied voltage, 15 kV. Background electrolyte: A and B, 50 mM phosphate buffer (pH 6) and SBE- β -CD 6 and 10 mg/ml, respectively; C, 50 mM Tris/HCl, pH 8, with 3 mg/ml of SBE- β -CD; D, 50 mM borate buffer, pH 9, and 20 mg/ml SBE- β -CD; injection, 5 psi*s 10^{-4} M of each racemic compound.

Composition of the Background Electrolyte

Chiral resolution can be strongly influenced by the composition of the BGE, and thus the selection of the appropriate buffer system, its concentration, its ionic strength, its pH, the presence of polymeric additives, and the content of the organic additive should be carefully considered. Decreasing the ionic strength of the BGE generally causes a reduction of migration time and resolution; peak tailing has also been observed.³⁴ The negative effect related to the low ionic strength of the buffer is due to electromigration dispersion that can be avoided using a BGE at a concentration 10^2 times higher than that of the sample. The peak shape can be controlled by selecting a co-ion with similar electrophoretic mobility to that of the analyte.

The addition of an organic solvent to the BGE can produce a negative effect on the binding constant of the inclusion-complex with cyclodextrins due to the competition between the organic additive and the analyte. The affinity of the analytes for the organic additive must be considered. However, organic solvents can improve the selectivity of the enantiomeric separation by differential influence on the binding constants of the two enantiomers. Propranolol enantiomers were not separated even using 40 mM of β -CD (urea was added to the BGE in order to increase the solubility of the CD), while separation was obtained when 30% (v/v) methanol was employed as a component of the buffer.¹⁵ Recently we demonstrated that the use of methanol can greatly improve the stereoselectivity of the separation of racemic non-steroidal anti-inflammatory drugs (NSAIDs). Racemic flurbiprofen was baseline resolved into its enantiomers at pH 5 using 30 mM of 2,3,6-tri-O-methyl- β -cyclodextrin, but the drug was poorly resolved when the CD concentration was lowered to 5 mM. The addition of methanol to the BGE containing 5 mM of chiral selector caused an increase in migration time as well as an increase in resolution, allowing baseline enantiomeric separation when the buffer contained 20% organic additive. The effect of methanol on the enantiomeric separation of propranolol and flurbiprofen is shown in Figure 14.

A theoretical model of the effect of organic solvent in the BGE when cyclodextrins are used as chiral selector in CE has been discussed by Wren.³⁴ When the CD concentration was at or below the optimum value (maximum resolution), the addition of organic solvent (methanol or acetonitrile) caused a reduction of resolution. This effect was due to the change in the formation constants of the inclusion-complexes, modifying the optimum CD concentration.

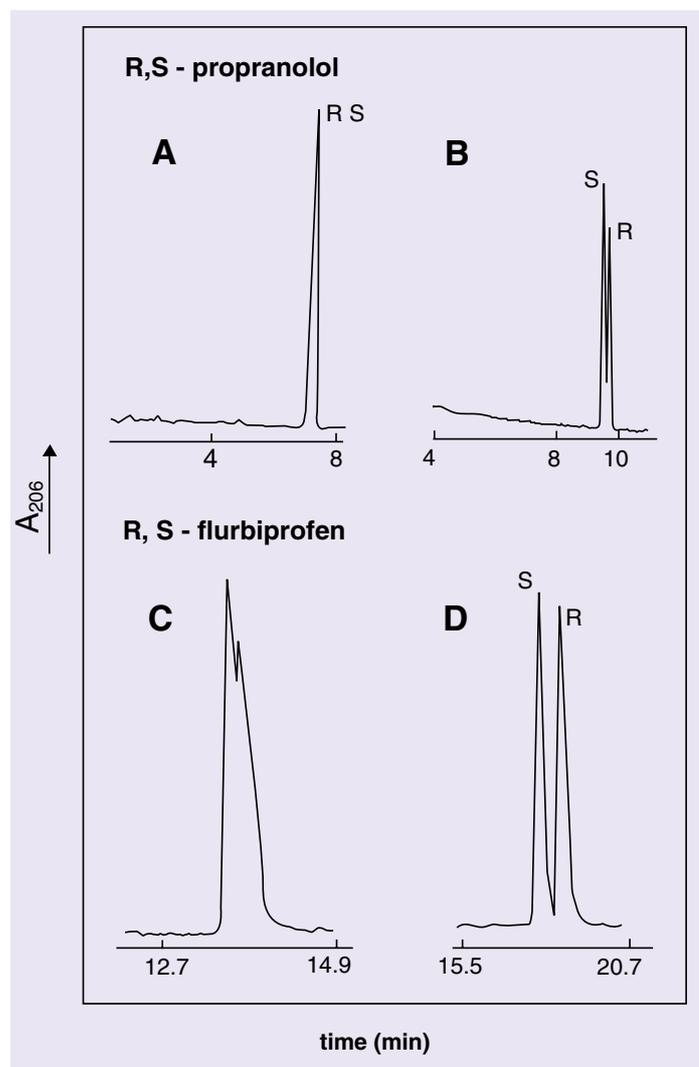


Fig. 14. Effect of methanol added to the BGE on the resolution of R,S-propranolol and R,S-flurbiprofen. A, (without methanol) and **B,** (with 30% methanol); apparatus HPE 100 system, capillary, 20 cm x 0.025 mm ID; background electrolyte, 0.1 M phosphate pH 2.5, 4 M urea and 40 mM β -CD. **C,** (without methanol) and **D,** (with 20 % methanol); apparatus BioFocus 3000 system; capillary, (coated) 36 cm x 0.05 mm ID; 0.1 M MES, pH 5, and 5 mM tri-OMe- β -CD; applied voltage 20 kV.

The influence of the pH of BGE was discussed by several authors and a theoretical model has also been described.⁸⁵

Strong electro-osmotic flow is not recommended for good enantiomeric separation by CE, because the time spent by the analyte for inclusion-complexation equilibrium is reduced. A coated capillary of short length can be used to improve the resolution or to perform the analysis in a reasonable time.

For many applications in the separation of enantiomers of basic compounds, acidic pH values have been used. At low pH the electro-osmotic flow is reduced, increasing the resolution.²⁹ Similar effects can be obtained by adding polymers such as hydroxyethylcellulose (HEC), hydroxypropylmethylcellulose (HPMC), or polyvinylalcohol (PVA) to the BGE, or by using capillaries coated internally, for example with polyacrylamide. The addition of polymers will reduce the eof and increase the viscosity. In all cases, the adsorption of analytes is also avoided or reduced.

The pH of the BGE not only influences the electro-osmotic flow but can also affect the charge of the analytes, and thus their effective mobility. When the analysis of weakly acidic compounds must be carried out in the presence of electro-osmotic flow and a native CD is used, both CD and analytes (uncharged and charged, respectively) are transported to the detector by the electro-osmotic flow. The migration time of the analyte in the presence of the CD is shorter than in its absence due to inclusion-complexation. Increasing the pH will cause an increase in migration time due to increased analyte charge, and this phenomenon can help the resolution. Of course, for the optimum experimental conditions, the magnitude of the electro-osmotic flow must be considered.

Capillary Temperature

Efficient temperature control is recommended to prevent loss in resolution when cyclodextrins are used as chiral selectors for enantiomeric separations. This can be easily obtained using a BioFocus electrophoresis system, which provides a system for circulating liquid coolant into the cartridge containing the capillary.

Temperature increases cause a decrease in buffer viscosity, and thus a decrease in migration time. Furthermore, a change in the capillary temperature can strongly influence the stability of the inclusion complex formed between analyte and cyclodextrin (generally an increase in temperature causes a decrease in the binding constant).⁸⁶ This effect is shown in Figure 15, where the separation of norepinephrine, epinephrine, and isoproterenol enantiomers was obtained using a dimethylated- β -CD.⁸⁷

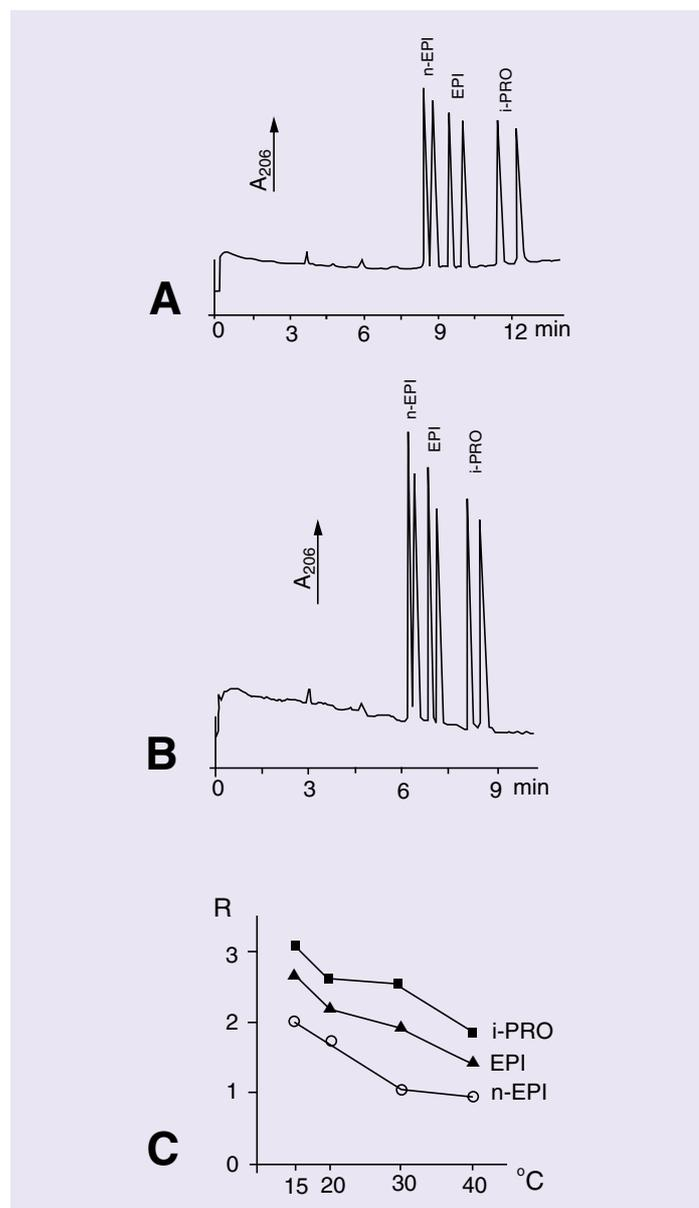


Fig. 15. Influence of the capillary temperature on the resolution of racemic norepinephrine (n-EPI), epinephrine (EPI), and isoproterenol (i-PRO) into their enantiomers. Apparatus, BioFocus 3000 system; capillary, (coated) 30 cm x 0.05 mm ID; background electrolyte, 50 mM phosphate buffer, pH 2.5, and 20 mM of di-OMe- β -cyclodextrin; applied voltage 10 kV; column temperature, **A**) 20 °C, **B**) 40 °C. **C**) resolution (R) vs. column temperature (modified from reference 87).

Cyclodextrins or their Derivatives in Capillary Gel Electrophoresis, Isotachopheresis, and Electrochromatography

Cyclodextrins or their derivatives have also been used as chiral selectors for enantiomeric separations with other capillary electrophoretic techniques, although there are fewer publications on this than on CZE and micellar electrokinetic chromatography. The use of capillary gel electrophoresis (CGE) for the separation of enantiomers was first described by Guttman et al.⁸ A capillary, filled with crosslinked polyacrylamide gel and also containing unbound α - or β - or γ -cyclodextrin, was used for the enantiomeric separation of several dansyl-amino acids. The electro-osmotic flow was suppressed by coating the capillary wall, and the separations were performed at pH 8.3. The CD type and its concentration, as well as the column temperature, influenced the migration time and the resolution of the amino acid derivatives studied. The addition of 10% methanol improved the separation of enantiomers. After this pioneering work, gel filled capillaries were not employed for chiral separations in CE until 1992, when Cruzado⁴⁵ used a modified β -CD (allyl carbamoylated- β -CD) copolymerized with acrylamide for the enantiomeric separation of several dansyl-amino acids and homatropine. The gel-modified cyclodextrin presented several drawbacks, including short lifetime and poor reproducibility.

The feasibility of enantiomeric separation using a chiral stationary phase (CSP) in CE has been shown by Mayer and Schurig.⁴³ The CSP was bound to the capillary wall after static thermal coating with Chirasil-Dex containing monokis-6-O-octamethylene-per-methyl- β or γ -cyclodextrin. The coating stability was tested by gas chromatography and by CE.

The separation of several racemic mixtures into their enantiomers was obtained with a borate/phosphate buffer at pH 7 in presence of electro-osmotic flow. Recently, the enantiomeric separation of both neutral and charged analytes, *e.g.*, benzoin, hexobarbital, and several amino acid derivatives, has been achieved using a β -CD bonded chiral stationary phase (CSP with 5 μ m particle diameter employed in HPLC).⁸⁸ The type of BGE influenced the direction of the electro-osmotic flow. The results obtained suggested perspectives on the use of a wide number of chiral phases.

Enantiomeric separations have also been performed by capillary isotachopheresis (ITP) employing cyclodextrins or their derivatives as chiral additives to the leading electrolyte (LE). The first paper by Smolkova-Keulemansova's group⁸⁹ dealt with the separation of some

pseudo ephedrine alkaloid enantiomers using β -CD or di-OMe- β -CD as a chiral additive to the LE. In this study, the CD type and concentration, and the pH of the LE, influenced the efficiency of the separation. After this pioneering study the same group showed the enantiomeric resolution of some phenothiazines with β - or γ -CD and discussed the importance of the counter-ion type on the ITP separation.⁹⁰ Ketotifen and its intermediate enantiomers⁹¹ have also been separated by this group using ITP. Furthermore the possibility of using a coupled column system, in which the two capillaries contained two different cyclodextrins, for a two-step separation has been demonstrated. The authors called the system two-dimensional separation. Unfortunately, this technique did not become popular, even though it offered several advantages over CZE, including a concentrating effect, analysis of compounds in traces, and the possibility of working with larger volumes of sample. Considering the advantages, we are convinced that in the near future ITP will be widely used, at least in combination with CZE, in order to improve the sensitivity of the separation method.

Chiral Crown-Ether

Another class of compounds used for enantiomeric resolution by the inclusion-complexation mechanism is represented by crown-ethers. These macromolecules, discovered in 1967 by Pederson,⁹³ are able to form inclusion complexes with several inorganic and organic ions in which the guest compound fits the cavity, forming weak bonds (ion-dipole) with the etheroatoms (O, S) of the crown. Here the inclusion-complexation is completely different than that of CDs; the hydrophilic part of the analyte is included, while with CDs it is the hydrophobic part. Although the inclusion of the amino group is necessary for enantiomeric separation, other interactions are required in order to perform chiral recognition. Figure 16 shows the chemical structure of the chiral 18-crown-6-ether tetracarboxylic acid used in capillary zone electrophoresis.

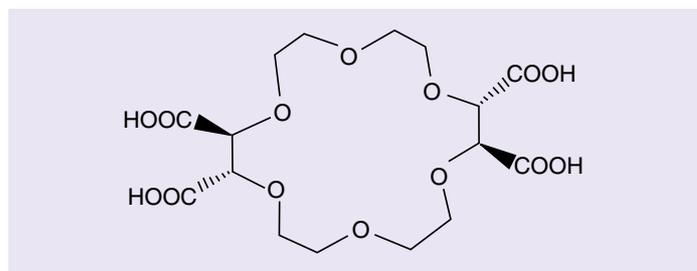


Fig. 16. Chemical structure of 18-Crown-6-ether tetracarboxylic acid.

Two different chiral resolution mechanisms have been proposed by Kuhn⁹⁴ in which the four carboxylic groups of the crown were involved in the chiral recognition process, either forming electrostatic interactions or causing a steric barrier effect with the included molecule.

The chemical structure of the analyte plays a very important role for chiral recognition with the modified crown-ether; in fact the presence of a primary amino group is fundamental for inclusion-complexation. For the barrier mechanism, nonpolar substituents on the asymmetric center have a strong influence. For example, naphthylethylamine was resolved, while phenylethylamine was not at all. Enantiomers of aliphatic amino acids with branched chains were better separated than those with linear chains (valine > norvaline). Finally the distance of the amino group from the chiral center has to be considered. Better resolution is obtained with short distances.

Other parameters which must be controlled for the optimization of the chiral separation using chiral crown ether by CZE include the pH of the BGE, the organic additive, and the buffer composition (K^+ , NH_4^+ , Na^+ ions should be avoided because they can compete with analytes in the inclusion process while Tris seems to be the most effective cation).⁹⁴ Applications using 18-crown-6-ether tetra carboxylic acid as chiral selector in capillary zone electrophoresis cover a wide range of amino derivatives, including racemic amino acids,³² amines and peptides,⁵⁶ and amino alcohols and sympathomimetic drugs.⁹⁵

4.2 Ligand Exchange

The ligand-exchange mechanism, first introduced by Davankov in chromatography,⁹⁶ was successfully applied in CE for the enantiomeric separation of dansyl amino acids^{57,58} and α -hydroxy acids.⁵⁹ The separation is based on the interaction between a metal complex with a chiral ligand, added to the BGE, and the two enantiomers to be resolved. L- or D-proline, L- or D-hydroxyproline and aspartame are the ligands which have been employed in CE.

Figure 17 shows the separation of D and L-phenyllactic (3-PhL) acid when Cu(II)-L-hydroxyproline complexes were used as the chiral selector. The two analytes migrate as anions, toward the anode in the coated capillary, while the positively charged metal chelate migrates in the opposite direction. The 3-PhL forms a ternary complex with Cu(II), causing a reduction of the effective mobility of both enantiomers; the two complexes possess different stability constants and thus are separated by the end of the electrophoretic run. The optimization

of the separation method requires the control of several parameters, namely the metal and chiral ligand type, their ratio and concentration, the pH, the ionic strength, the organic additive and the column temperature.

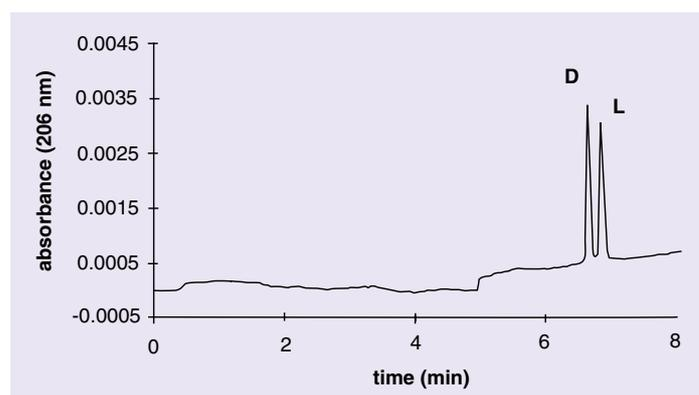


Fig. 17. Electropherograms of the separation of D- and L-phenyllactic acid using a ligand-exchange mechanism by CE. Apparatus, BioFocus 3000 system; capillary, (coated) 35 cm x 0.05 mm ID; background electrolyte, 0.02 M phosphate buffer, pH 4.4, and copper(II) acetate (2 mM)/ L-hydroxyproline (4 mM); applied voltage 15 kV, 15 A; injection 10 psi*s of 10^{-4} M of racemic 3-phenyllactic acid; carousel and capillary temperature, 25 °C.

One drawback that can be expected when ligand exchange is used in CE is the UV absorption of the metal complex, which reduces the sensitivity of the method. This drawback can be negated by using different approaches, *e.g.*, by filling only part of the capillary or by using a different detector such as laser induced fluorescence⁵⁷ or conductivity.

4.3 Chiral Micelles

The use of micelles in capillary electrophoresis was first introduced by Terabe *et al.* for the separation of neutral compounds. They called this technique micellar electrokinetic chromatography (MEKC).⁹⁷ Surfactants such as sodium dodecyl sulfate (SDS) were added to the aqueous solution in order to improve the selectivity of the separation of neutral analytes. With neutral-alkaline solutions, the electro-osmotic flow is relatively high and the bulk solution (analytes and micelles) moves toward the cathode. The migration of the negatively-charged micelles is retarded relative to neutral compounds. The analyte interacts with the two phases (aqueous and micellar) and thus its

migration is also retarded. The extent of the retardation is a function of the percentage of analyte distributed into the micellar phase.

Chiral recognition can be achieved in MEKC by including a chiral selector in the micellar system, either bound to the micelles or added to the electrolyte.

Sodium N-dodecanoyl-L-valinate (SDVal) and sodium N-dodecanoyl-L-alaninate (SDAla) were applied as chiral micelle selectors for the separation of several amino acid derivatives, namely 3,5-dinitrobenzoyl, 4-nitrobenzoyl and benzoyl-O-isopropyl esters.⁹⁸ SDVal was also used for the enantiomeric separation of phenylthiohydantoin amino acids by Terabe's group. Addition of methanol to the SDVal micellar solution as an organic modifier improved resolution, and the addition of urea significantly improved peak shapes.^{99,100}

Other very interesting chiral micelles are represented by bile salts, introduced for the first time in MEKC by Terabe *et al.*⁷ These surfactants, which possess a steroid nucleus and a side chain carrying a carboxylic group that can be conjugated with taurine, have both hydrophobic and hydrophilic properties. When taurine is conjugated, due to the presence of sulfonate groups, the chiral selector is charged even at low pH. Figure 18 shows the chemical structure of the main bile acid salts used in CE.

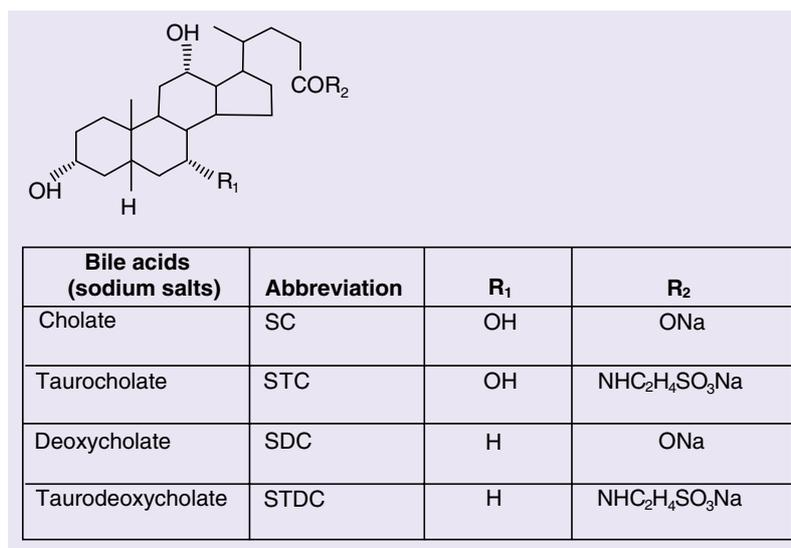


Fig. 18. Chemical structure of bile acid salts.

Sodium taurocholate (STC) and taurodeoxycholate have been used successfully as chiral selectors for the separation of several dansyl amino acids into their enantiomers.⁷ The reduction of the electro-osmotic flow by decreasing the pH to 3 produced good separations not obtainable under other experimental conditions. The same class of chiral surfactants proved to be useful chiral selectors for the separation of several racemic drugs into their enantiomers, *e.g.*, diltiazem and trimetoquinol,^{61,101} fonoldopam, 4-hydroxymephenytoin,¹⁰² laudanosine,¹⁰¹ tetrahydropapaveroline,¹⁰³ naphthol and binaphthyl derivatives.⁶² Digitonin,¹⁰⁰ glycyrrhizic acid, and β -escin¹⁰⁴ are natural surfactants which have also been used for chiral separations at a relatively low pH, and also combined with SDS in mixed micelle systems.

A very promising approach in MEKC is the use of chiral micelle polymers. Poly(sodium N-undecylenyl-L-valinate) has been shown to be a good chiral selector for the separation of 1-1'-bi-2-naphthol and laudanosine into their enantiomers. The advantages of such polymers over the monomers are their enhanced stability and rigidity. Furthermore, since the polymericelles have no CMC, they can be used at any concentration. Resolution was influenced by the polymer concentration and the pH. At low pH the charged polymerized micelles have a compact conformation, while at higher pH they lose it.¹⁰⁵

The addition of cyclodextrins to the micellar phase (CD-modified MEKC) enables the enantiomeric separation of uncharged compounds which could also be analyzed with charged cyclodextrins or with electrochromatography.

4.4 Affinity Interactions

Proteins as Chiral Selectors

Proteins are natural biopolymers with helical conformation able to interact selectively with a wide number of compounds of small size, such as pharmaceuticals. They have been used successfully as chiral selectors in capillary electrophoresis. The enantiomeric separation is based on stereoselective interactions between the protein and the two analytes, with the formation of two labile diastereoisomeric complexes during the electrophoretic run. Tanaka *et al.*⁶⁴ discussed a theoretical model for enantiomeric separation with proteins where the electrophoretic mobility of the free analyte (μ_s) and of the protein (μ_p), and the concentration of the protein (P), are correlated by the following equation:

$$(2) \quad \mu_{\text{app}} = \mu_{\text{eo}} + \frac{\mu_s + \mu_p K[P]}{1 + K[P]}$$

where μ_{app} , μ_{eo} , and K are the apparent mobilities of the analyte, electro-osmotic flow, and binding constant, respectively.

Calculating $\Delta\mu_{\text{app}} = \mu_{\text{app1}} - \mu_{\text{app2}}$, the following equation provides the optimum concentration of protein for the best chiral resolution (see equation 3)

$$(3) \quad [P]_{\text{op}} = (K_1 K_2)^{-1/2}$$

The theoretical model is similar to that discussed by Wren, where an uncharged chiral selector was used.⁷⁷ The Tanaka model was verified for the enantiomeric separation of several acidic compounds using avidin as a chiral selector. Avidin, a basic protein with a pI of 10.0–10.5, is positively charged at acidic/neutral pH and migrates in a direction opposite that of the analytes.

When proteins are used as chiral selectors in CE, several drawbacks have to be expected. The protein may adsorb on the capillary wall, the sensitivity in the low UV region can be reduced due to strong absorption of the chiral selector, and the protein may not be stable under the operating conditions. Coated capillaries were used in order to reduce the adsorption of the protein to the wall and to reduce the electroosmotic flow.

Figure 19 shows schematically the separation principle of affinity EKC for analysis of an acidic compound with avidin at acidic pH.

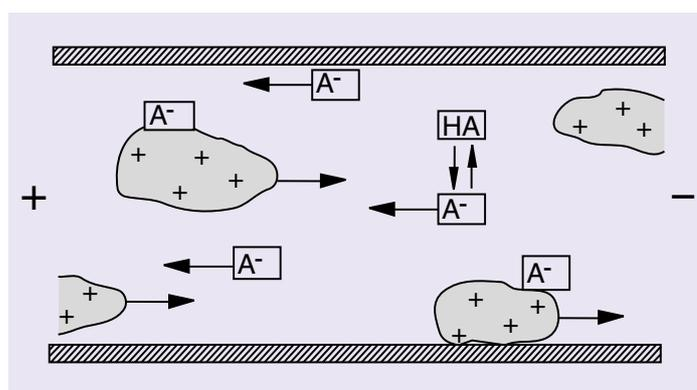


Fig. 19. Electrophoretic separation scheme for anions using avidin as a chiral selector (with permission from reference 48).

The separated enantiomeric compounds were vanilmandelic acid, warfarin, ibuprofen, ketoprofen, flurbiprofen, and the two diastereoisomers of folinic acid. Several parameters were investigated, *e.g.*, the concentration of the protein, the pH of the BGE, the content of the organic solvent and the capillary temperature. As shown in Figure 20, increasing the concentration of avidin from 0 to 25 μM gave better resolution of the enantiomeric pair of vanilmandelic acid at pH 4.

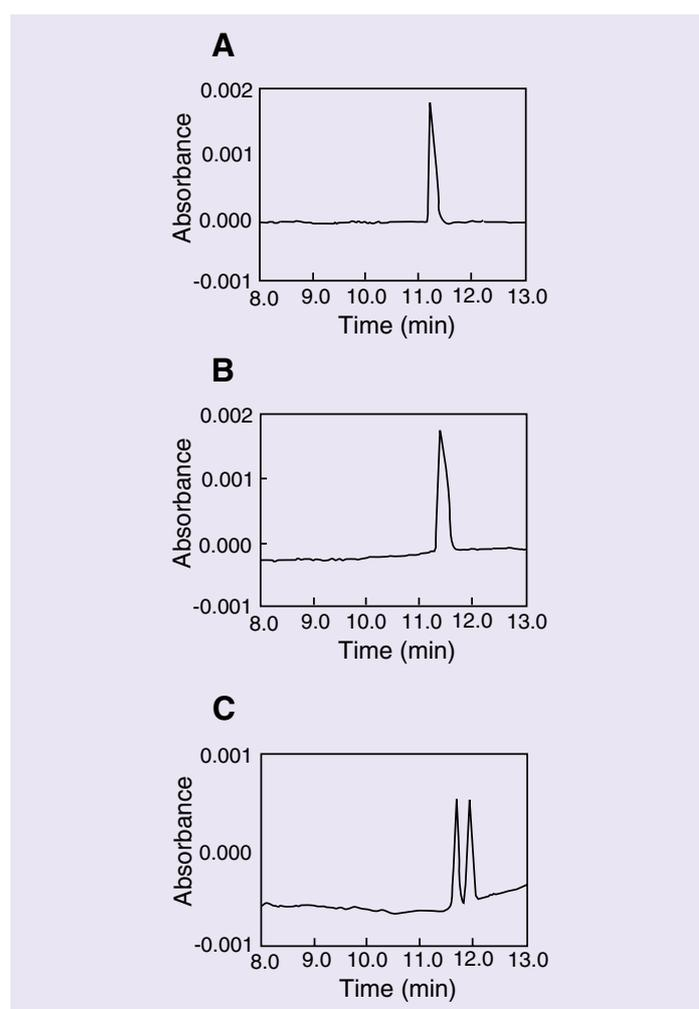


Fig. 20. Effect of the concentration of avidin on the enantiomeric separation of racemic vanilmandelic acid. Apparatus, BioFocus 3000 system; capillary, (coated) 36 cm x 0.05 mm ID; background electrolyte, 0.05 M phosphate buffer, pH 4, and avidin **A)** 0 M; **B)** 10 M and **C)** 25 M; applied voltage -12 kV; detection at 240 nm; injection, 1 psi*2s (sample 0.4 mg/ml five fold dilution) (with permission from reference 48).

It is worthwhile to mention here an interesting approach used by Valtcheva et al., employing the enzyme cellulase for the enantiomeric separation of several pharmaceutical compounds such as propranolol, pindolol, metoprolol, and labetalol.¹⁰ The chiral selector was negatively charged at the operating pH (pH=5.1), and the analytes migrated in the direction opposite to the enzyme. To optimize the method, the detector window was chiral selector-free, improving the detection limit and allowing use of a relatively high concentration of buffer (0.4 M) supplemented with 2-propanol.

Interactions with Saccharides (Linear Polymers) and Antibiotics

Linear polymers of α -(1-4)-linked D-glucose, termed maltodextrins, have proven to be good chiral selectors in CZE for the enantiomeric separation of non-steroidal anti-inflammatory drugs 2-APA NSAIDs (flurbiprofen, ibuprofen, carprofen, suprofen, and indoprofen) as well as coumarinic anticoagulant drugs.^{60,106}

An anionic biopolymer, heparin, was studied for the enantiomeric separation of several drugs including antimalarials and antihistamines. Heparin is a di-, tetra-, or hexasaccharide composed of uronic acid and glucosamine with a helical conformation in aqueous solutions and a molecular weight of 10–30,000. The enantiomeric separations were obtained at pH 4 and 5 and the chiral recognition was influenced by the size of the analyte and by the electrostatic interactions.¹⁰⁷

Another interesting class of compounds that can be used in CE as chiral selectors for enantiomeric separations is antibiotics. Vancomycin, a glycopeptide, was effective for the separation of enantiomeric dipeptides D-Ala-D-Ala/L-Ala-L-Ala and their acetyl derivatives¹⁰⁸ and for the separation of racemic derivatized tri- and tetrapeptides;¹⁰ in both studies the binding constants have been calculated. A recent paper demonstrated the applicability of the macrocyclic antibiotic rifamycin b for the separation of chiral compounds by affinity electrophoresis.¹¹⁰ Indirect UV detection was used because the antibiotic had strong absorption at low wavelengths.

4.5 Combination of Chiral/Chiral or Chiral/Achiral Selectors

Improvement in the selectivity of an enantiomeric separation can be obtained in CE by employing mixed selector systems. This approach consists of the use of a mixture of either chiral/achiral or chiral/chiral compounds. In a wide number of chiral separations, the combination of cyclodextrins with micelles was effective for the resolution of several uncharged compounds. Figure 21 shows schematically the electrophoretic separation when CD-MEKC is employed.

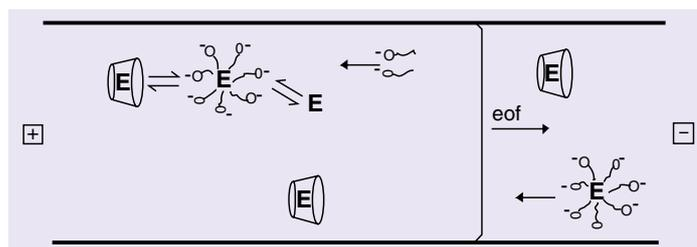


Fig. 21. Scheme of the electrophoretic separation when CD-modified MEKC is used.



The system is composed of an ionic micelle (usually SDS) and an aqueous phase (buffer + cyclodextrin). The analytes move with the eof, and are distributed between the micellar and the aqueous phase. The stereoselectivity is due to the interactions with CDs. The higher the interaction of the analyte with the chiral selector, the lower the migration time.

The selectivity can be modified by controlling several parameters, including surfactant type and concentration, CD type and concentration, organic additive, buffer type and concentration, and pH.

CD-MEKC allows the separation of enantiomers of amino acids, as dansyl-⁷⁸ or naphthalene-2,3-dicarboxaldehyde (NDA) derivatives,¹¹¹ using β - and/or γ -CD, as well as compounds of pharmaceutical interest, such as mephenytoin and its hydroxy derivative in urine.¹¹²

Figure 22 shows the electropherogram of an extracted undeglucuronidated urine sample from an extensive metabolizer (phenotyping). The presence of S-4-hydroxymephenytoin and R-mephenytoin was confirmed by spiking the sample mixture with the racemic

compounds (Figure 22 A and B) and comparing the UV spectra of the pure enantiomers with the separated peaks (Figure 22 D and E).

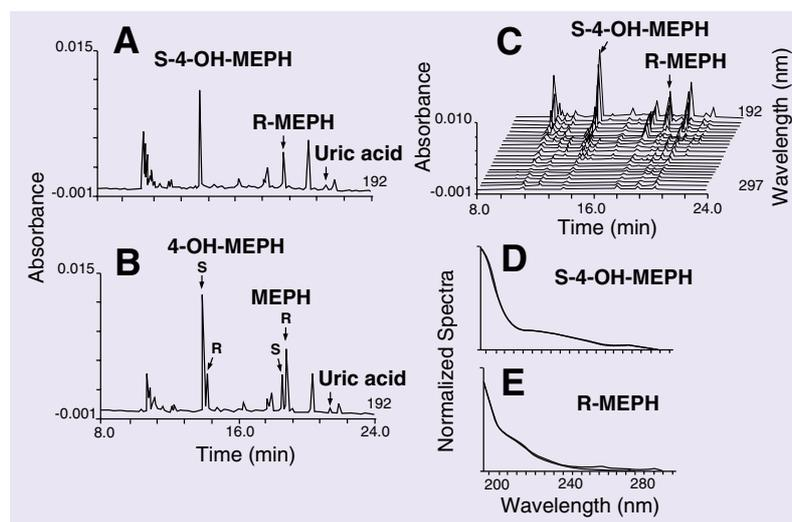
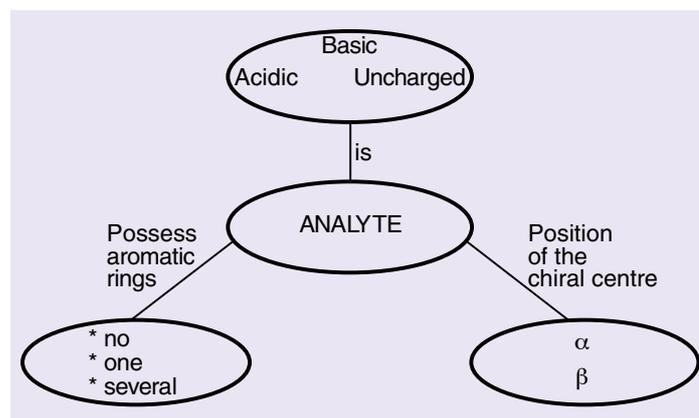


Fig. 22. Electrophoretic separation and characterization of S-4-hydroxymephenytoin (S-4-OH-MEPH) and R-mephenytoin (R-MEPH). Undeglucuronidated urine sample from an extensive metabolizer. Single and multiwavelength detection (**A** and **C**, respectively); in **B** the sample from **A** was spiked with the racemic compounds, while in **D** and **E** the spectra of the two detected compounds in the sample are compared with those of the pure enantiomers. (with permission from reference 112).

Another example of combining two chiral selectors, namely β -CD and L-tartrate, has been shown by us for the separation of several stereoisomers of cobalt(III) complexes with ethylenediamine and amino acids like Gly, Pro, and Phe. The L(+)-tartrate alone did not allow the resolution of the enantiomers of $\text{Co}[(\text{en})_2\text{D-Phe}]^{2+}$ and $\text{Co}[(\text{en})_2\text{L-Phe}]^{2+}$ and the addition of β -CD provided satisfactory results. In this separation both inclusion- and outer-sphere complexation are involved in the separation mechanism.⁶⁸

Section 5 Optimization of the Method

When a racemic mixture has to be separated into its enantiomers, it is necessary to consider several parameters in order to select the optimum electrophoretic experimental conditions. First the chemical structure has to be inspected to determine the following information:



When basic compounds must be analyzed (this case represents a wide number of pharmaceutical compounds), an acidic pH of the BGE is selected and the separation can be performed either in a coated or uncoated capillary. The recommended buffers are phosphate or citrate at pH 2.5 and 3.5, respectively (50–100 mM). In some cases a pH one unit lower than the pK_a of the enantiomers allows the migration of analytes as cations, and a pH higher than 2.5–3.5 could be considered.

The use of a pH >4–5 can be helpful because it enables the enantiomeric separation and allows sufficient electro-osmotic flow for the analysis of achiral impurities in the sample. Moreover the electro-osmotic flow should be controlled because it reduces the migration time and thus minimizes the time spent by the analytes in contact with the chiral selector.

Increasing the BGE concentration generally improves the resolution of enantiomers, but increases the migration times due to ion-pairing effects and reduction of electro-osmotic flow. Furthermore, peak-tailing and electromigration dispersion can be avoided. An upper limit for buffer concentration is dictated by the increase in current and resulting Joule heating which reduces the efficiency of the separation. Use of a zwitterionic buffer can be considered as a good solution to this problem.

When the enantiomeric separation of negatively charged compounds must be performed, the general rules for basic compounds must be considered. Selecting a pH for the BGE in the range 4.5–8 will charge (negatively) the two enantiomers and produce a sufficient electro-osmotic flow for movement of the analytes toward the cathode. The use of a coated capillary permits operation in a reversed polarity mode, and inversion of migration can be obtained. This

approach can be advantageously used when a minor component in a mixture of enantiomers has to be quantified.

The separation of uncharged enantiomers does not represent a problem in CE, because different electrophoretic mechanisms and chiral selector types are available. With uncharged analytes, a relatively strong electro-osmotic flow is necessary to perform the analysis within a reasonable time, and a charged chiral selector should be employed. Carboxymethylated or ethylated- β -cyclodextrin, sulfobutyl-ether- β -cyclodextrin, or a charged β -cyclodextrin polymer could be effective for chiral recognition. By inducing a negative charge, a selective increase in the migration times allows separation of the enantiomers.

Another approach for the enantiomeric separation of uncharged compounds is the use of surfactants with chiral selectors, *e.g.*, SDS-cyclodextrins. The partition of the analytes into the micellar phase causes a selective migration of neutral compounds, while the chiral selector allows chiral resolution. A buffer at a pH in the range 6–9 containing 25–100 mM of SDS supplemented with variable amounts of cyclodextrins is recommended. The effective mobility should be controlled by changing the composition of the buffer, to select the appropriate experimental conditions. The electro-osmotic flow should be controlled, and in some cases the addition of organic solvents can help the separation of enantiomers. If the uncharged compounds cannot be resolved into their enantiomers, different chiral selectors should be employed, *e.g.*, bile salts.

The capillary temperature should also be controlled when performing enantiomeric separations because, in general, an increase in temperature will result in a reduction of resolution. This parameter influences both the viscosity of the BGE (reduction of migration time) and mass transfer kinetics of analytes and chiral selectors. Thus preliminary studies should be carried out at 20–25 °C, and then a lower temperature should be used to optimize the chiral separation.

The capillary length can also influence the resolution of two enantiomers, and thus its selection should be done considering parameters such as applied voltage and capillary type. The use of longer capillaries can increase the resolution as well as the analysis time. When coated capillaries are used and the electro-osmotic flow is reduced or eliminated, good enantiomeric separations can be obtained even with capillaries of 20 cm length. For coated capillaries, lengths in the range of 20–35 cm are recommended, while with uncoated capillaries 50–70 cm lengths should be tried.

Selection of the Chiral Selector

The most simple chiral selectors that can be used for enantiomeric separation by CE are represented by native cyclodextrins because they are commercially available, not expensive, and several applications have been published.

If the analyzed compound does not contain aromatic rings, α -cyclodextrin or its derivatives can be selected. The same choice could be made if the analyte contains only one aromatic ring, unless substituents in ortho or meta position are present. Such substituents will cause steric hindrance effects, responsible for poor inclusion complexation.

When the compounds possess in their structure two aromatic rings (*e.g.* condensed), β -CD has to be selected, while for more than two aromatic groups γ -CD could be the appropriate chiral selector.

If successful enantiomeric separation cannot be achieved using native cyclodextrins, the wide number of modified ones (*e.g.*, dimethylated, trimethylated, hydroxypropylated) should be considered. Charged cyclodextrins may also be investigated, selecting the appropriate electrophoretic mechanism and experimental parameters.

As a general suggestion, run the compound initially in the absence of CDs and calculate the effective mobility. Then perform the run with the same BGE, containing a relatively low amount of CD (2.5–5 mM) in order to verify a decrease in effective mobility. If complexation is observed but insufficient resolution is obtained, increase the concentration of the chiral selector until satisfactory separation is achieved. In the case of β -CD, concentration screening is limited by its solubility (< 20 mM); solubility can be increased by addition of urea (4–8 M) or methanol or ethanol (< 30% v/v) to the BGE. Alternatively, modified cyclodextrins can be selected to improve the solubility or to introduce different secondary stereoselective bonds.

When compounds containing primary amines in their chemical structures, (*e.g.*, amino acids, peptides, drugs) have to be separated into their enantiomers, 18-crown-6-ether tetracarboxylic acid can be used advantageously. In this case attention should be paid to the selection of the BGE (pH and cation type, the latter should not compete with analytes in the inclusion-complex mechanism).

Bile salts with planar structures, namely sodium cholate, taurocholate, deoxycholate, and taurodeoxycholate, are recommended for chiral compounds with rigid skeletons. Optimization of the enantiomeric separation can be performed by modifying those parameters

involved in micellar systems, *e.g.*, buffer type and concentration, concentration of the bile salt, pH, and organic solvent.

For the enantiomeric separation of amino acids or hydroxy carboxylic acids, cyclodextrins are the simplest chiral selector to be used; however a ligand exchange electrophoretic system can also be considered. Copper(II)-amino acid or aspartame complexes could be selected, but unfortunately the detection limit is reduced due to the absorption of the chiral BGE.

Section 6 Quantitation

Quantitative analysis of chiral compounds is a very interesting and important topic that, although it has been shown to be possible by CE, unfortunately up to now is not yet well established. However, considering the good results obtained by several authors, the applicability of CE will be extended.

When a racemic mixture of a certain compound is analyzed by CE using an appropriate chiral selector, it should be separated into two peaks whose area ratio should be 1.00. The electrophoretic analysis of a racemic mixture of epinephrine¹⁴ showed two peaks with two different areas (49% of (-)-epinephrine and 51% of its antipode). In order to explain the observed differences, we have to consider that during the electrophoretic process the formation of two diastereoisomers takes place, and due to the different association constants, a different UV absorbance co-efficient could be expected.¹⁴ Furthermore, the analyte residence time in the detection path is different, the second enantiomer spending a longer time than the first one because it possesses a lower velocity.¹¹³ In order to solve this problem, it has been suggested that the measured areas be normalized to migration time,^{14,113} and/or that an internal standard (I.S.) be used.¹⁴

It has been shown that CE can be successfully used for quantitation of chiral impurities of drugs with good precision and accuracy. Quantitative analysis of trimetoquinol enantiomers by MEKC has been carried out with STDC as a chiral selector. The calibration graph was linear for concentrations of the R-antipode in the range 2.5–15% and less than 1% of this compound could be detected.¹⁰¹

Peterson and Trowbridge performed quantitative determination of (-)-epinephrine/(+)-epinephrine with and without an internal standard (pseudophephrine). The correlation coefficient and the RSD were found to be 0.9998 and 1.3%, respectively using the I.S., while

without I.S. they found 0.9992 and 3.3%, respectively. The authors analyzed a pharmaceutical formulation with different ages (5, 10, and 29 months) and found the presence of (+)-epinephrine in the range 1.2–2.3% (this antipode should not be present at all).

The analysis of the antiviral compound 2'-deoxy-3'-thiacytidine (BCH189), a nucleoside analog containing 0.3% of the undesired (-)-antipode, was done using a BGE at pH 2.5 and 50 mM of di-OMe- β -CD with a detection limit < 0.1%.¹¹⁴

A capillary electrophoretic method has been validated for the determination of the chiral purity of fluparoxan. Detector linearity was quite good in the studied range 1.5–125% of a target concentration (1.25 mg/ml) and the limit of quantitation and detection were found to be 1% and 0.3%, respectively. The concentration of β -CD used for the method was relatively high (150 mM) and thus urea and isopropanol were added to the BGE.¹¹⁵

Section 7 Applications

Chiral capillary electrophoresis has been applied, until now, mainly for enantiomeric separation of compounds of pharmaceutical interest. Resolution mechanisms, binding constants, and parameters affecting chiral analysis have been widely discussed. Several applications dealing with biological fluids are also available, as well as those dealing with quantitative analysis. The former topic is only at the beginning, and we are convinced that in the near future will be extensively developed.

For simplicity and to avoid repetition of some information given previously, Table 2 presents selected applications for enantiomeric separation of compounds of pharmaceutical interest.

Table 2. Selected Applications on the Enantiomeric Separation of Drugs

| Compound | Chiral selector | CE type | References |
|---|---|-------------|------------|
| adrenaline | di-OMe- β -CD | CZE | 14, 116 |
| atenolol | di-OMe- β -CD tri-OMe- β -CD | CZE | 41, 42 |
| basic drugs (terbutaline, isoproterenol, metaproterenol, synephrine, metanephrine, salbutamol, epinephrine, norphenylephrine, ephedrine, ψ -ephedrine, octopamine,) | macrocyclic antibiotic (rifamycin B) | CZE | 67 |
| basic drugs (nor-phenylephrine, ketamine, nor-ephedrine, ephedrine, epinephrine, terbutaline, propranolol) | negative β -CD polymer | CZE | 55 |
| basic drugs (atenolol, bupivacaine, ephedrine, homatropine, ketamine, metoprolol, N-methylpseudoephedrine, norephedrine, octopamine, pindolol, terbutaline) | (S)-N- dodecanoylvaline (S)-N-dodecoxy- carbonylvaline | MEKC | 117 |
| bupivacaine | di-OMe- β -CD | CZE/MEKC | 40 |
| carvediol | di-OMe- β -CD | CZE | 40 |
| chloramphenicol | di-OMe- β -CD | CZE | 38 |
| chlorpheniramine | β -CD | CZE or MEKC | 118, 119 |
| cicletanine | γ -CD | MEKC | 120 |
| clenbuterol | β -CD | CZE | 34 |
| clenbuterol | SBE- β -CD | CZE | 49 |
| cyclophosphamide | glycoprotein | affinity | 121 |
| diltiazem | STDC | MEKC | 103 |
| dimethindene | SBE- β -CD | CZE | 49 |
| ephedrine | di-OMe- β -CD | CZE | 13 |
| epinephrine | di-OMe- β -CD | CZE | 13, 116 |
| ergot alkaloids (lisuride, meluol, nicergoline,terguride) | γ -CD | CZE | 36 |
| herbicides (phenoxy acids) | α -, β -, γ -, di-OMe- β -CD | MEKC | 122 |
| hexobarbital | Chirasil- β -CD | EHC | 123 |
| mefloquine | SBE- β -CD | CZE | 49 |
| mephenytoin hydroxy metabolite | SDS- β -CD β -CD-STDC | MEKC | 102, 112 |
| methotrexate | vancomycin | CZE | 66 |
| mianserine | SBE- β -CD | CZE | 49 |

(continued on the next page)

Table 2. (continued)

| Compound | Chiral selector | CE type | References |
|---|-----------------------|---------------|------------|
| naproxen | HP- β -CD | CZE | 124 |
| NSAIDs (APA) (flurbiprofen, indoprofen, carprofen, ketoprofen, suprofen, fenoprofen, ibuprofen, naproxen) | vancomycin | CZE | 66 |
| NSAIDs (APA) (flurbiprofen, ketoprofen, fenoprofen, ibuprofen, suprofen) | tri-OMe- β -CD | CZE | 125 |
| NSAIDs (APA) (ibuprofen, cicloprofen, flurbiprofen, carprofen) | Chirasil- β -CD | EHC | 123 |
| NSAIDs (APA) (warfarin, ibuprofen, ketoprofen, flurbiprofen) | avidin | affinity | 64 |
| octopamine | di-OMe- β -CD | CZE | 87 |
| quinagolide | crown-ether | CZE | 94 |
| racemethorphan, racemorphan | SDS- β -CD | MEKC | 126 |
| salbutamol (and its chiral and achiral impurities) | di-OMe- β -CD | CZE | 127 |
| thalidomide | SBE- β -CD | CZE | 33, 49 |
| trimetoquinol and its isomers | dextran sulfate | affinity MEKC | 128 |
| trimetoquinol | STDC | MEKC | 103 |
| warfarin | BSA | affinity | 63 |
| warfarin | SD-Val | MEKC | 129 |
| warfarin | SBE- β -CD | CZE | 52 |
| warfarin | di-OMe- β -CD | CZE | 130 |

Section 8 Conclusions

The enantiomeric separation of different classes of compounds can be obtained easily and rapidly by CE techniques, with good reproducibility and at a relatively low cost. Direct and indirect separation methods can be successfully applied, but the former seems to be the most popular due to its simplicity.

Cyclodextrins have become the most popular chiral selectors in CE, since they are stable and quite soluble in the buffers commonly

used in CE, and because the faster equilibration of the CD-solute complex provides high efficiency and good peak symmetry. The hydroxyl groups of the CD (primary and secondary) can be easily modified, using chemical reactions, to obtain CDs with different properties, thus extending the application range by allowing different electrophoretic separation mechanisms. However, several other chiral selectors are available, *e.g.*, chiral crown ethers for compounds containing amino groups, proteins, chiral surfactants, saccharides, and recently, antibiotics. It has been shown that the enantiomeric resolution is strongly influenced by several parameters, including chiral selector type and concentration, composition of the BGE (ionic strength, ion type and concentration, pH, organic solvent), polymeric additives in the BGE, applied voltage, and capillary temperature.

The advantages of CE over other separation methods, *e.g.* HPLC, include higher efficiency, shorter analysis time, and lower costs (only a few μL of chiral buffer are required). Poor detection limits and limited preparative capability are the main disadvantages not yet resolved; in fact the loading of relatively high amounts of sample is not possible in CE. Some solutions to this problem have been shown, *e.g.*, the combination of ITP and CE.

Section 9 References

1. Debowski, J., Sybilska, D. and Jurczak, J., Resolution of some chiral mandelic acid derivatives into enantiomers by reversed-phase high performance liquid chromatography via α - and β -cyclodextrin inclusion complexes, *J. Chromatogr.*, **282**, 83 (1983).
2. Ward, T. J. and Armstrong, D. W., Improved cyclodextrin chiral phases: a comparison and review, *J. Liq. Chromatogr.*, **9**, 407 (1986).
3. Blaschke, G., Chromatographic resolution of chiral drugs on polyamides and cellulose triacetate, *J. Liq. Chromatogr.*, **9**, 341 (1986).
4. Singh, N. H., F. N. Pasutto, F. N., Coutts, R. T. and Jamali, F., Gas chromatographic separation of optically active anti-inflammatory 2-arylpropionic acids using (+) or (-)-amphetamine as derivatizing reagent, *J. Chromatogr.*, **378**, 125 (1986).
5. Kobor, F. and Schomburg, G., 6-tert-butyl-dimethylsilyl-2,3-dimethyl- α -cyclodextrin, β -cyclodextrin, and γ -cyclodextrin, dissolved in polysiloxanes, as chiral selectors for gas chromatography – influence of selector concentration and polysiloxane matrix polarity on enantioselectivity, *HRC-J. High. Res. Chromatogr.*, **16**, 693 (1993).
6. Armstrong, D. W., Faulkner, Jr. and Han, S. M., Use of hydroxypropyl and hydroxyethyl-derivatized β -cyclodextrins for the thin-layer chromatographic separation of enantiomers and diastereomers, *J. Chromatogr.*, **452**, 323 (1988).
7. Terabe, S., Shibata, H. and Miyashita, Y., Chiral separation by electrokinetic chromatography with bile salt micelles, *J. Chromatogr.*, **480**, 403 (1989).

8. Guttman, A., Paulus, A., Cohen, A. S., Grinberg, N. and Karger, B. L., Use of complexing agents for selective separation in high performance capillary electrophoresis, chiral resolution via cyclodextrins incorporated within polyacrylamide gel column, *J. Chromatogr.*, **448**, 41 (1988).
9. Schmitt, T. and H. Engelhardt, H., Charged and uncharged cyclodextrins as chiral selectors in capillary electrophoresis, *Chromatographia*, **37**, 475 (1993).
10. Valtcheva, L., Mohammed, J., Pettersson, G. and Hjerten, S., Chiral separation of b-blockers by high performance capillary electrophoresis based on non-immobilized cellulase as enantioselective protein, *J. Chromatogr.*, **638**, 263 (1993).
11. Khun, R., Erni, F., Bereuter, T. and Hausler, J., Chiral recognition and enantiomeric resolution based on host-guest complexation with crown-ethers in capillary zone electrophoresis, *Anal. Chem.*, **64**, 2815 (1992).
12. Belder, D. and Schomburg, G., Chiral separations of basic and acidic compounds in modified capillaries using cyclodextrin-modified capillary zone electrophoresis, *J. Chromatogr. A.*, **666**, 351 (1994).
13. Fanali, S., Separation of optical isomers by capillary zone electrophoresis based on host-guest complexation, *J. Chromatogr.*, **474**, 441 (1989).
14. Fanali, S., and Bocek, P., Enantiomeric resolution by using capillary zone electrophoresis: resolution of racemic tryptophan and determination of the enantiomer composition of commercial pharmaceutical epinephrine, *Electrophoresis*, **11**, 757 (1990).
15. Fanali, S., Use of cyclodextrins in capillary zone electrophoresis. Resolution of terbutaline and propranolol enantiomers, *J. Chromatogr.*, **545**, 437 (1991).
16. Foret, F., Krivankova, L. and Bocek, P., Capillary Zone Electrophoresis, VCH Verlagsgesellschaft mbH, Weinheim-New York-Basel-Cambridge-Tokyo, 1993.
17. Snopek, J., Jelinek, I. and Smolkova-Keulemansova, E., Chiral separation by analytical electromigration methods, *J. Chromatogr.*, **609**, 1 (1992).
18. Snopek, J. and Smolkova-Keulemansova, E., in D. Duchene (Editor), New trends in cyclodextrins and derivatives, Edition de Santé, Paris, p. 483, 1991.
19. Fanali, S., in N. A. Guzman (Editor), Capillary electrophoresis technology, Marcel Dekker, Inc, New York-Basel-Hong Kong, p. 731, 1993.
20. Ward, T. J., Chiral media for capillary electrophoresis, *Anal. Chem.*, **66**, A632 (1994).
21. Novotny, M., Soini, H. and Stefansson, M., Chiral separation through capillary electromigration methods, *Anal. Chem.*, **66**, 646A (1994).
22. Rogan, M. M., Altria, K. D. and Goodall, D. M., Enantioselective separations using capillary electrophoresis, *Chirality*, **6**, 25 (1994).
23. Terabe, S., Otsuka, K. and Nishi, H., Separation of enantiomers by capillary electrophoretic technique, *J. Chromatogr.*, **666**, 295 (1994).
24. Fanali, S., Cristalli, M., Vespalec, R. and Bocek, P., in A. Chrambach, M. J. Dunn and B. J. Radola (Editors), Advances in electrophoresis, VCH Verlagsgesellschaft mbH, Weinheim-New York-Basel-Cambridge-Tokyo, p. 3, 1994.
25. Fanali, S., Chiral separations by capillary electrophoresis, *Electrophoresis*, **15**, 753 (1994).
26. Cahn R. S., Ingold, C. K. and Prelog, V., Specification of molecular chirality, *Angew. Chem. Int. Ed. Engl.*, **5**, 385 (1966).

27. Schutzner, W., Fanali, S., Rizzi, A. and Kenndler, E., Separation of diastereomeric derivatives of enantiomers by capillary zone electrophoresis with a polymer network: use of polyvinylpyrrolidone as buffer additive, *J. Chromatogr.*, **639**, 375 (1993).
28. Schutzner, W., Caponecchi, G., Fanali, S., Rizzi, A. and Kenndler, E., Improved separation of diastereomeric derivatives of enantiomers by a physical network of linear polyvinylpyrrolidone applied as pseudophase in capillary zone electrophoresis, *Electrophoresis*, **15**, 769 (1994).
29. Terabe, S., Otsuka, K. and Nishi, H., Separation of enantiomers by capillary electrophoretic techniques, *J. Chromatogr. A.*, **666**, 295 (1994).
30. Nardi, A., Ossicini, L. and Fanali, S., Use of cyclodextrins in capillary zone electrophoresis for the separation of optical isomers. Resolution of racemic tryptophan derivatives, *Chirality*, **4**, 56 (1992).
31. Tanaka, M., Asano, S., Yoshinago, M., Kawaguchi, Y., Tetsumi, T. and Shono, T., Separation of racemates by capillary zone electrophoresis based on complexation with cyclodextrins, *Fresemitus J. Anal. Chem.*, **339**, 63 (1991).
32. Kuhn, R., Stoecklin, F. and Erni, F., Chiral separations by host-guest complexation with cyclodextrin and crown-ether in capillary zone electrophoresis, *Chromatographia*, **33**, 32 (1992).
33. Aumatell, A., Wells, R. J. and Wong, D. K. Y., Enantiomeric differentiation of a wide range of pharmacologically active substances by capillary electrophoresis using modified cyclodextrins, *J. Chromatogr. A*, **686**, 293 (1994).
34. Altria, K. D., Goodall, D. M. and Rogan, M. M., Chiral separation of β -amino alcohols by capillary electrophoresis using cyclodextrins as buffer additives. I. Effect of varying parameters, *Chromatographia*, **34**, 19 (1992).
35. Nielsen, M. W. F., Chiral separation of basic drugs using cyclodextrin-modified capillary zone electrophoresis, *Anal. Chem.*, **65**, 885 (1993).
36. Fanali, S., Fliieger, M., Steinerova, N. and Nardi, A., Use of cyclodextrins for the enantioselective separation of ergot alkaloids by capillary zone electrophoresis, *Electrophoresis*, **13**, 39 (1992).
37. Belder, D. and Schomburg, G., Modification of silica surfaces for CZE by adsorption of non-ionic hydrophilic polymers or use of radial electric fields, *J. High Resol. Chromatogr.*, **15**, 686 (1992).
38. Snopek, J., Soini, H., Novotny, M., Smolkova-Keulemansova, E. and Jelinek, I., Selected applications of cyclodextrin selectors in capillary electrophoresis, *J. Chromatogr.*, **559**, 215 (1991).
39. Sepaniak, M. J., Cole, R. D., and Clark, B. K., Use of native and chemically modified cyclodextrin for the capillary electrophoretic separation of enantiomers, *J. Liq. Chromatogr.*, **15**, 1023 (1992).
40. Soini, H., Riekkola, M. L. and Novotny, M. V., Chiral separation of basic drugs and quantitation of bupivacaine enantiomers in serum by capillary electrophoresis with modified cyclodextrin buffers, *J. Chromatogr.*, **608**, 265 (1992).
41. Peterson, T. E., Separation of drug stereoisomers by capillary electrophoresis with cyclodextrins, *J. Chromatogr.*, **630**, 353 (1993).
42. Wren, S. A. C. and Rowe, Theoretical aspects of chiral separation in capillary electrophoresis. III. Application to β -blockers, *J. Chromatogr.*, **635**, 113 (1993).
43. Mayer, S. and Schurig, V., Enantiomer separation by electrochromatography on capillaries coated with chirasil-dex, *J. High Resol. Chromatogr.*, **15**, 129 (1992).

44. Penn, S. G., Goodall, D. M. and Loran, J. S., Differential binding of tioconazole enantiomers to hydroxypropyl-beta-cyclodextrin studied by capillary electrophoresis, *J. Chromatogr.*, **636**, 149 (1993).
45. Cruzado, I. D. and Vigh, G., Chiral separations by capillary electrophoresis using cyclodextrin-containing gels, *J. Chromatogr.*, **608**, 421 (1992).
46. Ingelse, B. A., Everaerts, F. M., Desiderio, C. and Fanali, S., A study on the enantiomeric separation by capillary electrophoresis using a soluble neutral β -cyclodextrin polymer, *J. Chromatogr. A*, (1994).
47. Nardi, A., Eliseev, A., Bocek, P. and Fanali, S., Use of charged and neutral cyclodextrins in capillary zone electrophoresis: enantiomeric resolution of some 2-hydroxy acids, *J. Chromatogr.*, **638**, 247 (1993).
48. Dette, C., Ebel and, S. and Terabe, S., Neutral and anionic cyclodextrins in capillary electrophoresis: enantiomeric separation of ephedrine and related compounds, *Electrophoresis*, **15**, 799 (1994).
49. Chankvetadze, B., Endresz, G. and Blaschke, G., About some aspects of the use of charged cyclodextrins for capillary electrophoresis enantioseparation, *Electrophoresis*, **15**, 804 (1994).
50. Tait, R. J., Thompson, D. O., Stella, V. J. and Stobaugh, J. F., Sulfobutyl ether β -cyclodextrin as a chiral discriminator for use with capillary electrophoresis, *Anal. Chem.*, **66**, 4013 (1994).
51. Lurie, I. S., Klein, R. F. X., Dalcason, T. A., LeBelle, M. J., Brenneisen, R. and Weinberger, R. E., Chiral resolution of cationic drugs of forensic interest by capillary electrophoresis with mixtures of neutral and anionic cyclodextrins, *Anal. Chem.*, **66**, 4019 (1994).
52. Desiderio, C. and Fanali, S., Use of negatively charged sulfobutyl-ether- β -cyclodextrin for enantiomeric separation by capillary electrophoresis, *J. Chromatogr. A*, (submitted) (1995).
53. Terabe, S., Electrokinetic chromatography: an interface between electrophoresis and chromatography, *Trends Anal. Chem.*, **8**, 129 (1989).
54. Schmitt, T. and Engelhardt, H., Derivatized cyclodextrins for the separation of enantiomers in capillary electrophoresis, *J. High Resol. Chromatogr.*, **16**, 525 (1993).
55. Aturki, Z. and Fanali, S., Use of beta-cyclodextrin polymer as a chiral selector in capillary electrophoresis, *J. Chromatogr. A*, **680**, 137 (1994).
56. Kuhn, R., Erni, F., Bereuter, T. and Hausler, J., Chiral recognition and enantiomeric resolution based on host-guest complexation with crown ethers in capillary zone electrophoresis, *Anal. Chem.*, **64**, 2815 (1992).
57. Gozel, P., Gassmann, E., Michelsen, H. and Zare, R. N., Electrokinetic resolution of amino acid enantiomers with copper(II)- aspartame support electrolyte, *Anal. Chem.*, **59**, 44 (1987).
58. Gassmann, E., Kuo, J. E. and Zare, R. N., Electrokinetic separation of chiral compounds, *Science*, **230**, 813 (1985).
59. Desiderio, C., Aturki, Z. and Fanali, S., Separation of α -hydroxyacid enantiomers by high performance capillary electrophoresis using copper(II)-L-amino acid and copper(II)-aspartame complexes as chiral selectors in the background electrolyte, *Electrophoresis*, **15**, 864 (1994).
60. D'Hulst, A. and Verbeke, N., Chiral separations by capillary electrophoresis with oligosaccharides, *J. Chromatogr.*, **608**, 275 (1992).
61. Nishi, H., Fukuyama, T., Matsuo, M. and Terabe, S., Chiral separation of optical isomeric drugs using micellar electrokinetic chromatography and bile salts, *J. Microcol. Sep.*, **1**, 234 (1989).

62. Cole, R. O., Sepaniak, M. J. and Hinze, W. L., Optimization of binaphthyl enantiomer separations by capillary zone electrophoresis using mobile phases containing bile salts and organic solvents, *J. High Resol. Chromatogr.*, **13**, 579 (1990).
63. Busch, S., Kraak, J.C. and Poppe, H., Chiral separations by complexation with proteins in capillary zone electrophoresis, *J. Chromatogr.*, **635**, 119 (1993).
64. Tanaka, Y., Matsubara, N. and Terabe, S., Separation of enantiomers by affinity electrokinetic chromatography using avidin, *Electrophoresis*, **15**, 848 (1994).
65. Yang, J. and Hage, D.S., Chiral separations in capillary electrophoresis using human serum albumin as a buffer additive, *Anal. Chem.*, **66**, 2719 (1994).
66. Armstrong, D.W., Rundlett, K. L. and Chen, J. R., Evaluation of the macrocyclic antibiotic vancomycin as a chiral selector for capillary electrophoresis, *Chirality*, **6**, 496 (1994).
67. Armstrong, D.W., Tang, Y. B., Chen, S. S., Zhou, Y. W., Bagwill, C. and Chen, J. R., Macrocyclic antibiotics as a new class of chiral selectors for liquid chromatography, *Anal. Chem.*, **66**, 1473 (1994).
68. Fanali, S., Ossicini, L., Foret, F. and Bocek, P., Resolution of optical isomers by capillary zone electrophoresis. Study of enantiomeric and diastereoisomeric Co(III)-complexes with ethylenediamine and amino acid ligands, *J. Microcol. Sep.*, **1**, 190 (1989).
69. Szejtli, J., Cyclodextrins and their inclusion complexes, Akademiai Kiado, Budapest, 1982.
70. Ward, T. J. and Armstrong, D. W., in M. Zief and L. J. Crane (Editors), Chromatographic chiral separation, Marcell Dekker, New York, p. 131, 1988.
71. Snopek, J., Jelinek, I. and Smolkova-Keulemansova, E., Micellar, Inclusion and metal-complex enantioselective pseudophases in high performance electromigration methods, *J. Chromatogr.*, **452**, 571 (1988).
72. Pharr, D. Y., Fu, Z. F., Smith, T. K. and Hinze, W. L., Solubilization of cyclodextrins for analytical applications, *Anal. Chem.*, **61**, 275 (1989).
73. Czugler, M., Eckle, E. and Stezowski, J., Crystal and molecular structure of a 2,6-tetradeca-O-methyl- β -cyclodextrin-adamantanol 1:1 inclusion complex, *J. Chem. Soc. Chem. Comm.*, **1291** (1981).
74. Croft, A. P. and Bartsch, R. A., Synthesis of chemically modified cyclodextrins, *Tetrahedron*, **39**, 1417 (1983).
75. Valko, I. E., Billiet, H. A. H., Frank, J. and Luyben, K. C. A. M., Effect of the degree of substitution of (2-hydroxy)propyl- β -cyclodextrin on the enantioseparation of organic acids by capillary electrophoresis, *J. Chromatogr. A.*, **678**, 139 (1994).
76. Yoshinaga, M., Asano, S., Tanaka, M. and Shono, T., Separation of racemates by capillary zone electrophoresis based on complexation with cyclodextrin-derivatives, *Analytical Science*, **7**, 257 (1991).
77. Wren, S. A. C. and Rowe, R. C., Theoretical aspects of chiral separation in capillary electrophoresis. I. Initial evaluation of a model, *J. Chromatogr.*, **603**, 235 (1992).
78. Terabe, S., Miyashita, Y., Ishihama, Y. and Shibata, O., Cyclodextrin-modified micellar electrokinetic chromatography – separation of hydrophobic and enantiomeric compounds, *J. Chromatogr.*, **636**, 47 (1993).
79. Ueda, T., Mitchell, R., Kitamura, F., Metcalf, T., Kuwana, T. and Nakamoto, A., Separation of naphthalene-2,3-dicarboxaldehyde-labeled amino acids by high-performance capillary electrophoresis with laser-induced fluorescence detection, *J. Chromatogr.*, **593**, 265 (1992).

80. Yamashoji, Y., Ariga, T., Asano, S. and Tanaka, M., Chiral Recognition and Enantiomeric Separation of Alanine *b*-naphthylamide by Cyclodextrins, *Anal. Chim. Acta*, **268**, 39 (1992).
81. Stefansson, M. and Novotny, M., Electrophoretic resolution of monosaccharide *e*-Enantiomers in borate-oligosaccharide complexation media, *J. Am. Chem. Soc.*, **115**, 11573 (1993).
82. Schmitt, T. and Engelhardt, H., Charged and uncharged cyclodextrins as chiral selectors in capillary electrophoresis, *Chromatographia*, **37**, 475 (1993).
83. Tait, R. J., Skanchy, D. J., Thompson, D. P., Chetwyn, N. C., Dunshee, D. A., Rajewsky, R. A., Stella, V. J. and Stobaugh, J. F., Characterization of sulfoalkyl ether derivatives of *b*-cyclodextrin by capillary electrophoresis with indirect UV detection, *J. Pharm. Biomed. Anal.*, **10**, 615 (1992).
84. Wren, S. A. C., Theory of chiral separation in capillary electrophoresis, *J. Chromatogr.*, **636**, 57 (1993).
85. Rawjee, Y. Y., Williams, R. L. and Vigh, G., Capillary electrophoretic chiral separations using cyclodextrin additives. III. Peak resolution surfaces for ibuprofen and homatropine as a function of the pH and the concentration of beta-cyclodextrin, *J. Chromatogr. A.*, **680**, 599 (1994).
86. Hinze, W. L., Applications of cyclodextrins in chromatographic separations and purification methods, *Sep. Purif. Meth.*, **10**, 159 (1981).
87. Schutzner, W. and Fanali, S., Enantiomers resolution in capillary zone electrophoresis by using cyclodextrins, *Electrophoresis*, **13**, 687 (1992).
88. Li, S. and Lloyd, D. K., Packed-capillary electrochromatographic separation of the enantiomers of neutral and anionic compounds using beta-cyclodextrin as a chiral selector – effect of operating parameters and comparison with free-solution capillary electrophoresis, *J. Chromatogr. A.*, **666**, 321 (1994).
89. Snopek, J., Jelinek, I. and Smolkova-Keulemansova, E., VIII: Use of cyclodextrins in isotachopheresis. IV. The influence of cyclodextrins on the chiral resolution of ephedrine alkaloid enantiomers, *J. Chromatogr.*, **438**, 211 (1988).
90. Jelinek, I., Dohnal, J., Snopek, J. and Smolkova-Keulemansova, E., Use of cyclodextrins in isotachopheresis. VII. Resolution of structurally related and chiral phenothiazins, *J. Chromatogr.*, **464**, 139 (1989).
91. Jelinek, I., Snopek, J. and Smolkova-Keulemansova, E., Use of cyclodextrins in isotachopheresis. VIII. The separation of ketotifen and its polar intermediate enantiomers, *J. Chromatogr.*, **439**, 386 (1988).
92. Snopek, J., Jelinek, I. and Smolkova-Keulemansova, E., Use of cyclodextrin in isotachopheresis. VIII. Two dimensional chiral separation in isotachopheresis, *J. Chromatogr.*, **472**, 308 (1989).
93. Pedersen, C., Cyclic polyethers and their complexes with metal salts, *J. Am. Chem. Soc.*, **89**, 2495 (1967).
94. Kuhn, R., Wagner, J., Walbroehl, Y. and Bereuter, T., Potential and limitations of an optically active crown ether for chiral separation in capillary zone electrophoresis, *Electrophoresis*, **15**, 828 (1994).
95. Hohne, E., Kraus, G. J. and Gubitz, G., Capillary zone electrophoresis of the enantiomers of aminoalcohols based on host-guest complexation with a chiral crown ether, *J. High Resol. Chromatogr.*, **15**, 698 (1992).
96. Davankov, V. A. and Rogozhin, S. V., Ligand chromatography as a novel method for the investigation of mixed complexes: stereoselective effects in α -amino acid copper(II) complexes, *J. Chromatogr.*, **60**, 280 (1971).

97. Terabe, S., Otsuka, K., Ichikawa, K., Tsuchiya, A. and Ando, T., Electrokinetic separations with micellar solution and open-tubular capillaries, *Anal. Chem.*, **56**, 111 (1984).
98. Dobashi, A., Ono, T., Hara, S. and Yamaguchi, J., Enantioselective hydrophobic entanglement of enantiomeric solutes with chiral functionalized micelles by electrokinetic chromatography, *J. Chromatogr.*, **480**, 413 (1989).
99. Otsuka, K. and Terabe, S., Effects of methanol and urea on optical resolution of phenylthiohydantoin DL-amino acids by micellar electrokinetic chromatography with sodium-N-dodecanoyl-L-valinate, *Electrophoresis*, **11**, 982 (1990).
100. Otsuka, K. and Terabe, S., Enantiomeric resolution by micellar electrokinetic chromatography with chiral surfactants, *J. Chromatogr.*, **515**, 221 (1990).
101. Nishi, H., Fukuyama, T., Matsuo, M. and Terabe, S., Chiral separation of trimetoquinol hydrochloride and related compounds by micellar electrokinetic chromatography using sodium taurodeoxycholate solutions and application to optical isomers, *Anal. Chim. Acta*, **236**, 281 (1990).
102. Okafo, G. N., Bintz, C., Clarke, S. E. and Camilleri, P., Micellar electrokinetic capillary chromatography in a mixture of taurodeoxycholic acid and β -cyclodextrin, *J. Chem. Soc. Chem. Commun.*, **1189** (1992).
103. Nishi, H., Fukuyama, T., Matsuo, M. and Terabe, S., Chiral separation of diltiazem, trimetoquinol and related compounds by micellar electrokinetic chromatography with bile salts, *J. Chromatogr.*, **515**, 233 (1990).
104. Ishihama, Y. and Terabe, S., Enantiomeric separation by micellar electrokinetic chromatography using saponins, *J. Liq. Chromatogr.*, **16**, 933 (1993).
105. Wang, J. A. and Warner, I. M., Chiral separations using micellar electrokinetic capillary chromatography and a polymerized chiral micelle, *Anal. Chem.*, **66**, 3773 (1994).
106. D'Hulst, A. and Verbeke, N., Quantitation in chiral capillary electrophoresis: theoretical and practical considerations, *Electrophoresis*, **15**, 854 (1994).
107. Stalcup, A. M. and Agyei, N. M., Heparin: A chiral mobile-phase additive for capillary zone electrophoresis, *Anal. Chem.*, **66**, 3054 (1994).
108. Carpenter, J. L., Camilleri, P., Dhanak, D. and Goodall, D. M., A study of the binding of vancomycin to dipeptides using capillary electrophoresis, *J. Chem. Soc. Chem. Commun.*, **804** (1992).
109. Chu, Y. H. and Whitesides, G. M., Affinity capillary electrophoresis can simultaneously measure binding constants of multiple peptides to vancomycin, *J. Org. Chem.*, **57**, 3524 (1992).
110. Armstrong, D. W., Rundlett, K. and Reid, G. L., Use of a macrocyclic antibiotic, rifamycin b, and indirect detection for the resolution of racemic amino alcohols by CE, *Anal. Chem.*, **66**, 1690 (1994).
111. Ueda, T., Kitamura, F., Mitchell, R., Metcalf, T., Kuwana, T. and Nakamoto, A., Chiral separation of naphthalene-2,3-dicarboxaldehyde labeled amino acid enantiomers by cyclodextrins modified micellar electrokinetic chromatography with laser induced fluorescence, *Anal. Chem.*, **63**, 2979 (1991).
112. Desiderio, C., Fanali, S., Kupfer, A. and Thormann, W., Analysis of mephentoin, 4-hydroxymephentoin and 4-hydroxyphenytoin enantiomers in human urine by cyclodextrin micellar electrokinetic capillary chromatography: Simple determination of a hydroxylation polymorphism in man, *Electrophoresis*, **15**, 87 (1994).
113. Altria, K. D., Essential peak area normalisation for quantitative impurity content determination by capillary electrophoresis, *Chromatographia*, **35**, 177 (1993).

114. Rogan, M. M., Drake, C. D., Goodall, D. M. and Altria, K. D., Enantioselective enzymatic biotransformation of 2'-deoxy-3'-thiacytidine (BCH 189) monitored by capillary electrophoresis, *Anal. Biochem.*, **208**, 343 (1993).
115. Altria, K. D., Walsh, A. R. and Smith, N. W., Validation of a capillary electrophoresis method for the enantiomeric purity testing of fluparoxan, *J. Chromatogr.*, **645**, 193 (1993).
116. Peterson, T. E. and Trowbridge, D., Quantitation of l-epinephrine and determination of the d/l- epinephrine enantiomers in a pharmaceutical formulation by capillary electrophoresis, *J. Chromatogr.*, **603**, 298 (1992).
117. Mazzeo, J. R., Grover, E. R., Swartz, M. E. and Petersen, J. S., Novel chiral surfactant for the separation of enantiomers by micellar electrokinetic capillary chromatography, *J. Chromatogr. A.*, **680**, 125 (1994).
118. Ong, C. P., Ng, C. L., Lee, H. K. and Li, S. F. Y., Determination of antihistamines in pharmaceutical by capillary electrophoresis, *J. Chromatogr.*, **588**, 335 (1991).
119. Otsuka, K. and Terabe, T., Optical resolution of chlorpheniramine by cyclodextrin added capillary zone electrophoresis and cyclodextrin modified micellar electrokinetic chromatography, *J. Liq. Chromatogr.*, **16**, 945 (1993).
120. Prunonosa, J., Obach, R., Diez-Coscon, A. and Gouesclou, L., Determination of cicletaline enantiomers in plasma by high performance capillary electrophoresis, *J. Chromatogr.*, **574**, 127 (1992).
121. Li, S. and Lloyd, D. K., Direct chiral separations by capillary electrophoresis using capillaries packed with an alpha(1)-acid glycoprotein chiral stationary phase, *Anal. Chem.*, **65**, 3684 (1993).
122. Nielen, M. W. F., (Enantio-)separation of phenoxy acid herbicides using capillary zone electrophoresis, *J. Chromatogr.*, **637**, 81 (1993).
123. Mayer, S. and Schurig, V., Enantiomer separation using mobile and immobile cyclodextrin derivatives with electromigration, *Electrophoresis*, **15**, 835 (1994).
124. Guttman, A. and Cooke, N., Practical aspects in chiral separation of pharmaceuticals by capillary electrophoresis. II. Quantitative separation of naproxen enantiomers, *J. Chromatogr. A*, **685**, 155 (1994).
125. Fanali, S. and Aturki, Z., Use of cyclodextrins in capillary electrophoresis for the chiral resolution of some 2-arylpropionic acid non-steroidal anti-inflammatory drugs, *J. Chromatogr. A*, (in press) (1995)
126. Aumatell, A. and Wells, R. J., Chiral differentiation of the optical isomers of racemethorphan and racemorphan in urine by capillary zone electrophoresis, *J. Chromatogr. Sci.*, **31**, 502 (1993).
127. Rogan, M. M., Altria, K. D. and Goodall, D. M., Enantiomeric separation of salbutamol and related impurities using capillary electrophoresis, *Electrophoresis*, **15**, 808 (1994).
128. Nishi, H., Nakamura, K., Nakai, H., Sato T. and Terabe, S., Enantiomeric separation of drugs by affinity electrokinetic chromatography using dextran sulphate, *Electrophoresis*, **15**, 1335 (1994).
129. Otsuka, K., Kawahara, J., Tatekawa, K. and Terabe, S., Chiral separations by micellar electrokinetic chromatography with sodium N-dodecanoyl-L-valinate, *J. Chromatogr.*, **559**, 209 (1991).
130. Gareil, P., Gramond, J. P. and Guyon, F., Separation and determination of warfarin enantiomers in human plasma samples by capillary zone electrophoresis using methylated b-cyclodextrin-containing electrolyte, *J. Chromatogr.*, **615**, 317 (1993).

Acknowledgments

We thank Drs. Zeineb Aturki, Emanuela Camera, Claudia Desiderio, Wolfgang Schutzner, Mr. Gabriele Caponecchi, and Michele Cristalli for their cooperation in the research work.



**Bio-Rad
Laboratories**

*Life Science
Group*

*U.S. (800) 4BIORAD California (510) 741-1000 New York (516) 756-2575 Australia
02-805-5000 Austria (1) 877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
China (01) 2046622 Denmark 39 17 9947 Finland 90 804 2200 France (1) 49 60 68
34 Germany 089 31884-0 India 91-11-461-0103 Italy 02-21609 1 Japan 03-5811-
6270 Hong Kong 7893300 The Netherlands 0318-540666 New Zealand 09-443
3099 Singapore (65) 4432529 Spain (91) 661 70 85 Sweden 46 (0) 735 83 00
Switzerland 01-809 55 55 United Kingdom 0800 181134*

Bulletin 1973 US/EG REVA

95-0284 0695 SIG 051995 Printed in USA