



# Capillary Electrophoresis Application Notes Index

Molecule	Bulletin	Description
Glutathione	1575-1	Oxidized and reduced forms separated, coating essential
Hemoglobin	1575-2	4 variants analyzed, coating essential
500 mer PCR prod	1575-4	PCR check
Pentapeptides	1575-7	Resolution of reverse sequences
Substance P	1575-8	Peptide fragments separated
9 Small Peptides	1575-9	9 component mix (hormones and enkephalins)
Tears	1575-10	Size separation of 4 proteins
Caffeine	1575-11	Load by EEO sample of low inherent mobility
1hGH	1575-12	Deamidation
Angiotensin	1575-13	Deamidation
Nucleotides	1575-14	8 component mix (UDP,GTP,etc)
Vitamin C	1575-15	Isomers
BSA/Myog/Carb A	1575-16	Rapid,sharp resolution/no protein adsorbtion
BSA	1575-18	Coated vs uncoated
HSA	1575-19	CZE vs IEF Separation
B Vitamins	1575-22	Resolution of anionic, cationic and neutral species
Food Dyes	1575-23	Neutral molecules, purity check
Casein & Whey	1575-25	Dairy products application
Dipeptide Isomers	1575-26	Metal ions in buffer enhance separation
Porcine Insulin	1575-28	Purity check, coated capillary essential
Rare Earth Metals	1575-29	Rapid, indirect detection of 18 metal ions
IEF pI 4.2-9.6	1575-31	7 proteins, separation by isoelectric point
BSA-4 methods	1575-32	Different comparison of protein methods
Dipeptide	1575-38	2 peaks for same molecule, different configuration
Sandostatin	1575-39	Purity check
Human Insulin	1575-41	Purity check
TRH	1575-42	Hormone and analog
Organic Acids	1575-43	Indirect detection of organic acids in wine
Daunorubicin	1575-44	Analysis of daunorubicin by MECC with UV scanning detection
Antibiotics	1575-45	Analysis of sulfonamide antibiotics

(continued on back)

Molecule	Bulletin	Description
Amphotericin B	1575-46	Analysis of amphotericins by MECC with visible scanning detection
Spider Venoms	1575-47	CZE separation of tarantula venoms at low and high pH
Hop Acids	1575-49	Analysis of hop bitter acids
Enantiomers	1575-51	Separation of Sympathomimetic Drug Enantiomers by CE
Nucleic acids	1575-52	Analysis of Amplified Nucleic acids by CE using Polymer Solutions
Lead in Food	1575-53	Analysis of Lead in Foods using CE
Peptides	1575-54	Analysis of Peptides by CE/MS
Drugs	1575-55	Analysis of Chlorpheniramine, Codeine, and Methadone by CE/MS
Flavonoids	1575-57	Analysis of Flavonoids by MEKC with Scanning Detection

To request application notes, call 1(800)4BIORAD [1(800)424-6723] or your local Bio-Rad Representative.



**Bio-Rad  
Laboratories**

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) **Bio-Rad Laboratories Main Office** 2000 Alfred Nobel Drive, Hercules, CA 94547, Ph. (510) 747-7000, Fx. (510) 747-5800  
**Also in:** **Australia** Ph. (02) 9974 2800, Fx. (02) 9974 2889 **Austria** Ph. (03) 877 89 07, Fx. (07) 876 5629 **Belgium** Ph. 09-385 55 77, Fx. 09-385 65 54  
**Canada** Ph. (905) 772-2777, Fx. (905) 772-2390 **China** Ph. 86-70-62057850, Fx. 86-70-62057876 **Denmark** Ph. 45 33 7799 47, Fx. 45 33 27 76 98  
**Finland** Ph. 358 (09) 804 2200, Fx. 358 (09) 804 7700 **France** Ph. (01) 43 90 4690, Fx. (01) 46 77 24 67 **Germany** Ph. 089 378 84-0, Fx. 089 378 84-700  
**Hong Kong** Ph. 852-2700-3300, Fx. 852-2700-3257 **India** Ph. (011) 461-0703, Fx. (011) 461-0765 **Israel** Ph. 03 957 4727, Fx. 03 957 4729  
**Italy** Ph. 02 276007, Fx. 02 27600399 **Japan** Ph. 03-6207-6270, Fx. 03-6207-6272 **Korea** Ph. 82-2-3473-4460, Fx. 82-2-3472-7000  
**The Netherlands** Ph. 31 378-640886, Fx. 31 378-642276 **New Zealand** Ph. 61-9-4752280, Fx. 61-9-4752284  
**Singapore** Ph. 65-2729877, Fx. 65-2734835 **Spain** Ph. (91) 667 70 85, Fx. (91) 667 96 98 **Sweden** Ph. 46 (08) 62780 00, Fx. 46 (08) 62784 00  
**Switzerland** Ph. 01-800 55 55, Fx. 01-800 55 00 **United Kingdom** Ph. 0800-787754, Fx. 01442-259776