



Capillary Electrophoresis
Application Note 11

Analysis of Caffeine by Micellar Electrokinetic Capillary Chromatography

Micellar electrokinetic capillary chromatography (MEKC) is a chromatographic technique performed using capillary electrophoresis instrumentation. Analytes are separated by partitioning between an aqueous electrolyte and a pseudophase of surfactant molecules organized in micelles. Analytes, micelles, and the bulk electrolyte are transported toward the detector by electroendosmosis (EOF). The method described here employs uncoated capillaries which, when used with an alkaline electrolyte, exhibit a high EOF toward the cathode. Sodium dodecyl sulfate is used at a concentration well above the critical micelle concentration to form the pseudophase. The surfactant

monomers and micelles are anionic, so have an electrophoretic mobility counter to the direction of EOF. Therefore, the movement of neutral hydrophobic analytes will be retarded by partitioning into the micellar pseudophase.

In this note, MEKC was used to analyze caffeine in samples of coffee, tea, and a diet cola soda. Samples were analyzed directly with no sample preparation. Prior to use, the capillary was conditioned by washing at 100 psi second for 15 minutes with 0.1 N HCl, 5 minutes with deionized water, 15 minutes with 0.1 N NaOH, 5 minutes with deionized water, and 15 minutes with run buffer.

Results

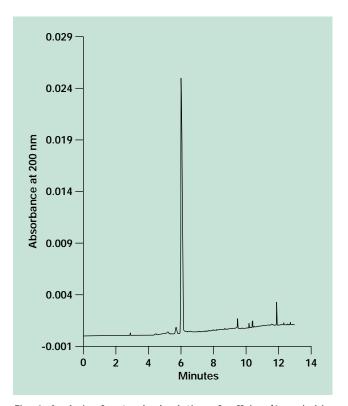


Fig. 1. Analysis of a standard solution of caffeine (1 mg/ml in water) by MEKC.

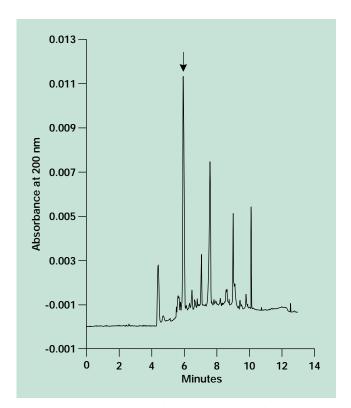


Fig. 2. Analysis of caffeine (indicated by arrow) in coffee by MEKC.

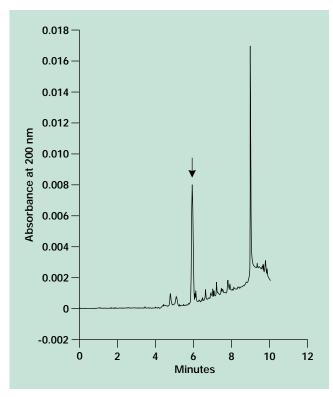


Fig. 3. Analysis of caffeine (indicated by arrow) in tea by MEKC.

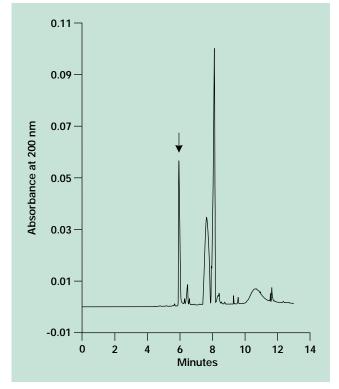


Fig. 4. Analysis of caffeine (indicated by arrow) in diet cola soda by MEKC.

Analysis Conditions

Instrument BioFocus® 3000 system
Polarity positive to negative
Capillary 50 cm x 50 µm, uncoated
Run buffer 0.05 M borate, pH 9.0

0.05 M SDS

Capillary purge 60 sec with run buffer Injection pressure at 10 psi*second

Run voltage 15 kV
Detection 200 nm
Cartridge temperature 20 °C
Autosampler temperature 20 °C



Bio-Rad Laboratories

150 9005 registered

erere.blo-radioom

Life Science Group Evo-Part Laboration via Main O Mon 2000 Amerika et Onine, Aeroeden, Oddronia 19647, ph. 1500 ph 5 5000, ph. 1500 ph 5 6000 Amerika et O 240 ph 24000, ph. 1524 ph 25000, ph. 1524 ph 2500 ph 2500

ıBulletin 1575-11 US/EG REV B 97-146 05/97 😋 🕬 🕮 7