



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 electroosmosis (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

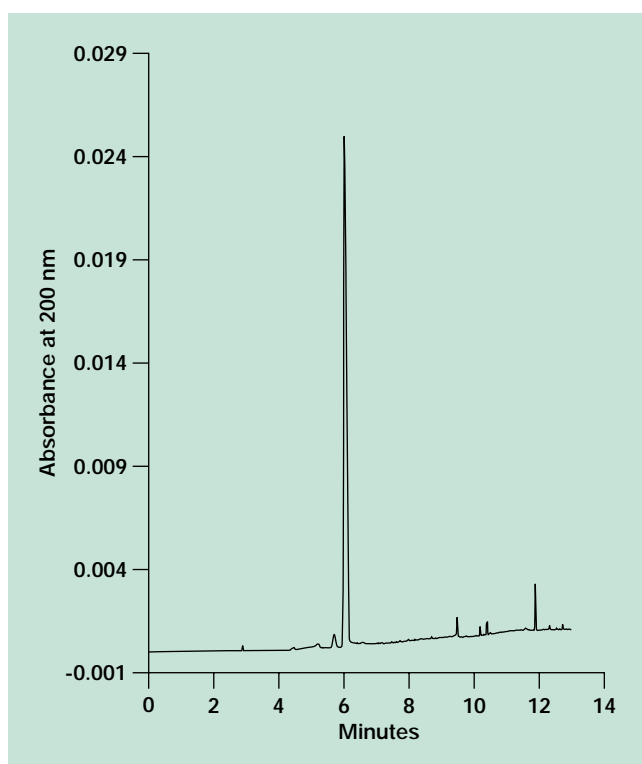


Fig. 1. Analysis of a standard solution of caffeine (1 mg/ml in water) by MEKC.

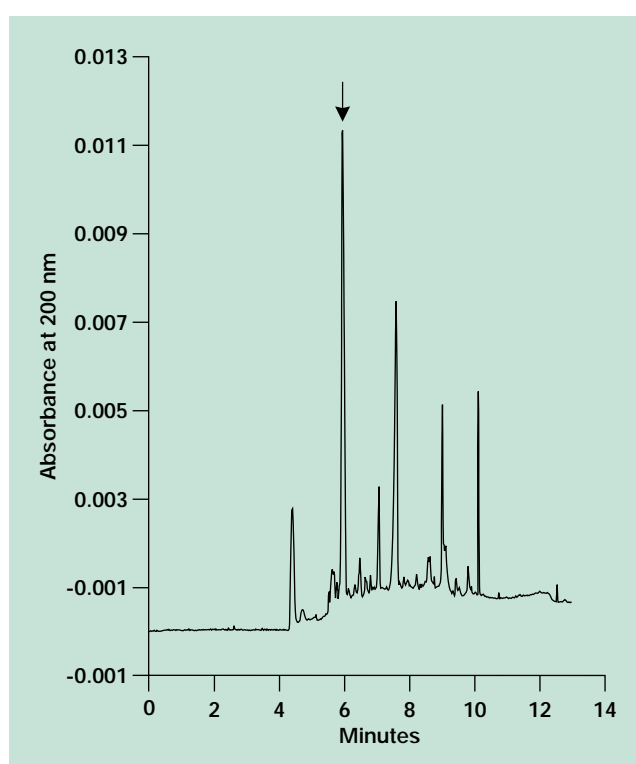


Fig. 2. Analysis of caffeine (indicated by arrow) in coffee by MEKC.

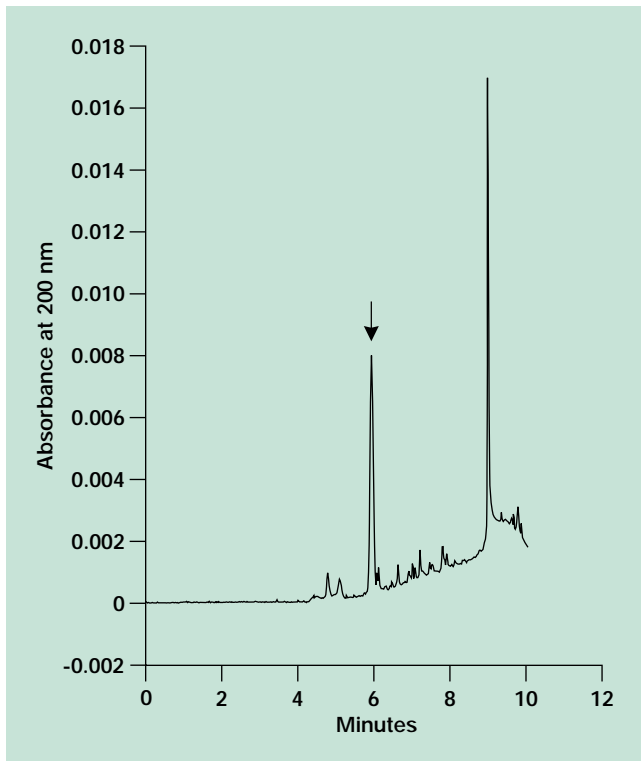


Fig. 3. Analysis of caffeine (indicated by arrow) in tea by MEKC.

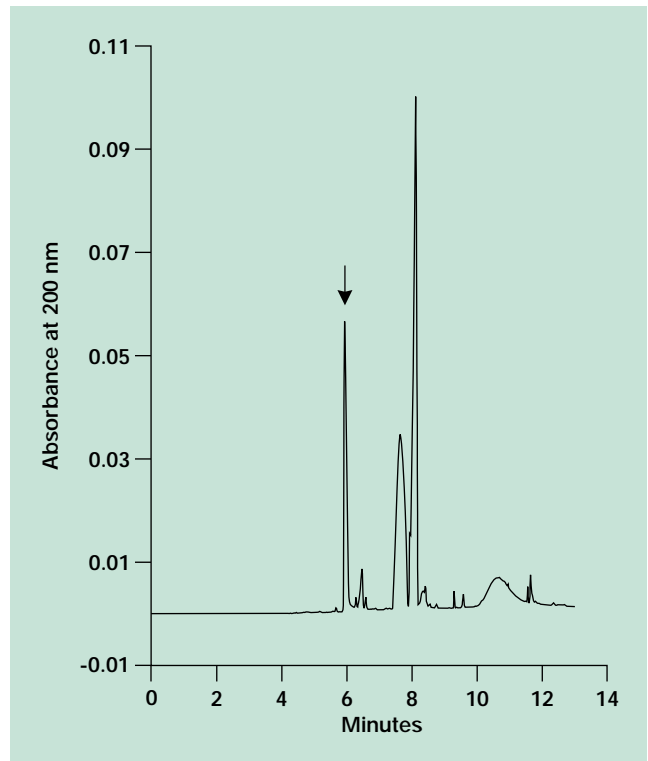


Fig. 4. Analysis of caffeine (indicated by arrow) in diet cola soda by MEKC.

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**

ISO 9001 registered www.bio-rad.com

Life Science
Group

Bio-Rad Laboratories North America 2000 Alfred Nobel Drive, Berkeley, California 94707, Ph. (510) 750-5000, Fx. (510) 750-5000
 Also in: Australia Ph. 02-4954-2300, Fx. 02-4954-2300 Austria Ph. (0)3 77 99 0 3, Fx. (0)3 76 98 29 Belgium Ph. 04-388 55 55, Fx. 04-388 65 54
 Canada Ph. (905) 752-2777, Fx. (905) 752-2900 China Ph. (86-30) 2096822, Fx. (86-30) 2055378 Denmark Ph. 39 579047, Fx. 39 27 5808
 Finland Ph. 00 304 2200, Fx. 00 304 3500 France Ph. (33) 49 00 46 00, Fx. (33) 49 75 24 67 Germany Ph. 089 38840, Fx. 089 38840 300
 Hong Kong Ph. 7398 800, Fx. 739 325 7 India Ph. 05-55465-0508, Fx. 05-55465-0765 Israel Ph. 08 953 4 327, Fx. 08 953 4 329
 Italy Ph. 02-2 900 3, Fx. 02-2 900 300 Japan Ph. 03-5355-8270, Fx. 03-5355-8272 The Netherlands Ph. 0618 38-240888, Fx. 0618 38-242296
 New Zealand Ph. 04-488 8000, Fx. 04-488 8007 Singapore Ph. (65) 2 22-9877, Fx. (65) 2 22-4883 Spain Ph. (34) 66 5 70 35, Fx. (34) 66 5 06 98
 Sweden Ph. 46 (0) 8 627 50 00, Fx. 46 (0) 8 627 54 00 Switzerland Ph. 05-800 55 55, Fx. 05-800 55 00
 United Kingdom Ph. 0800 38 5884, Fx. 05442 2295 38