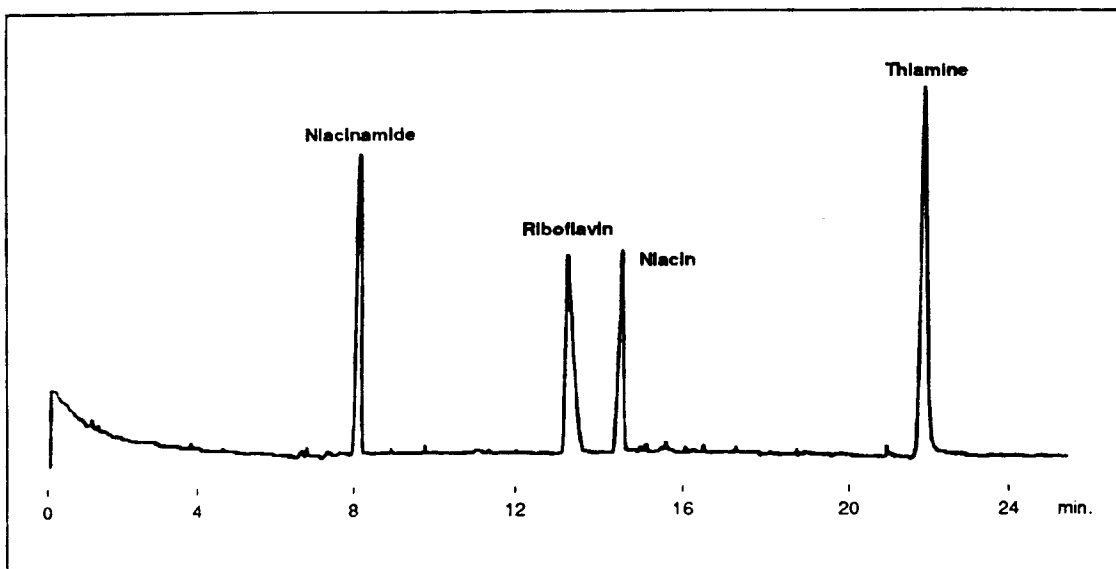


B Vitamins separated by MECC

Application
Note: 22

Micellar electrokinetic capillary chromatography (MECC) is a separation technique which permits resolution of anions, cations, and neutral species through the combined effects of electrophoretic migration and micellar partitioning. The electropherogram below shows a separation of 4 water soluble vitamins: niacinamide (neutral, 25 µg/ml), riboflavin (neutral, 100 µg/ml), niacin (anionic, 25 µg/ml), and thiamine (cationic, 100 µg/ml).



Capillary:	50 cm x 50 µm, uncoated
Buffer:	0.01 M sodium borate, 0.005 M sodium phosphate, 0.05 M SDS, pH 8.0
Load Conditions:	10 kV, 10 seconds
Run Conditions:	10 kV, constant voltage, ⊕-⊖ polarity
Detection:	UV, 200 nm, 0.032 AUFS