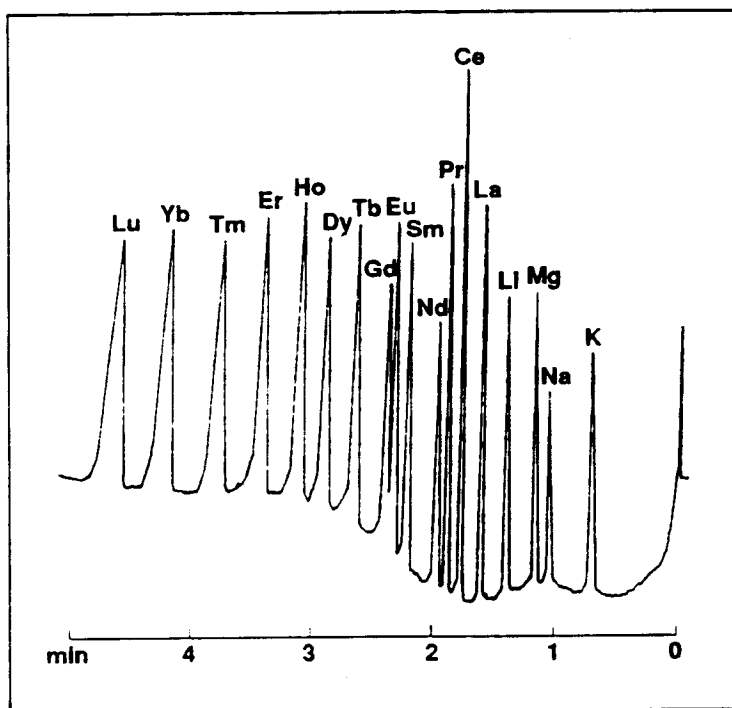


Rare Earth Metals

Application
Note: 29

Detection and quantitation of lanthanides is important in many environmental and industrial applications. The incorporation of complexing counterions with UV absorbing coions into the buffer allows indirect detection of cations by capillary electrophoresis. Eighteen different metals are resolved in less than five minutes in the electropherogram below.



Data provided by Frantisek Foret, Institute of Analytical Chemistry, Brno, Czechoslovakia

Capillary:	25 cm x 25 μ m, coated
Buffer:	0.03 M creatine-acetate, 0.004 M hydroxyisobutyric acid, pH 4.8
Load Conditions:	1.5 kV, 8 seconds
Run Conditions:	12 kV, constant voltage, \oplus - \ominus polarity
Detection:	UV, 220 nm, 0.1 AUFS

1575-29 1290 90-0291