



Analysis of Glutathione by Capillary Electrophoresis

Glutathione (L-glutamyl-L-cysteinyl-L-glycine) is a tripeptide found in animals, plants, and bacteria that acts as a biological redox agent, a coenzyme and cofactor, and serves as a substrate in certain coupling reactions. Oxidized and reduced forms can be separated by capillary electrophoresis with on-line detection at 200 nm.

Results

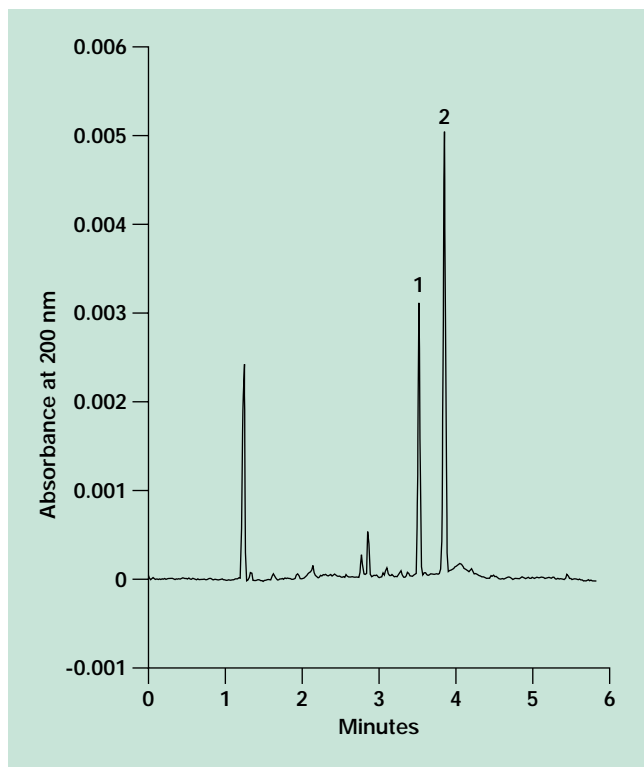


Fig. 1. Separation of oxidized (peak 1) and reduced (peak 2) glutathione by capillary zone electrophoresis.

Analysis Conditions

Instrument	BioFocus® 3000 system
Polarity	negative to positive
Capillary	24 cm x 50 µm, coated
Run buffer	0.05 M borate, pH 8.5
Capillary purge	60 seconds with run buffer
Injection	Electrokinetic at 4 kV for 4 seconds
Run voltage	10 kV
Detection	200 nm
Cartridge temperature	20 °C
Autosampler temperature	20 °C