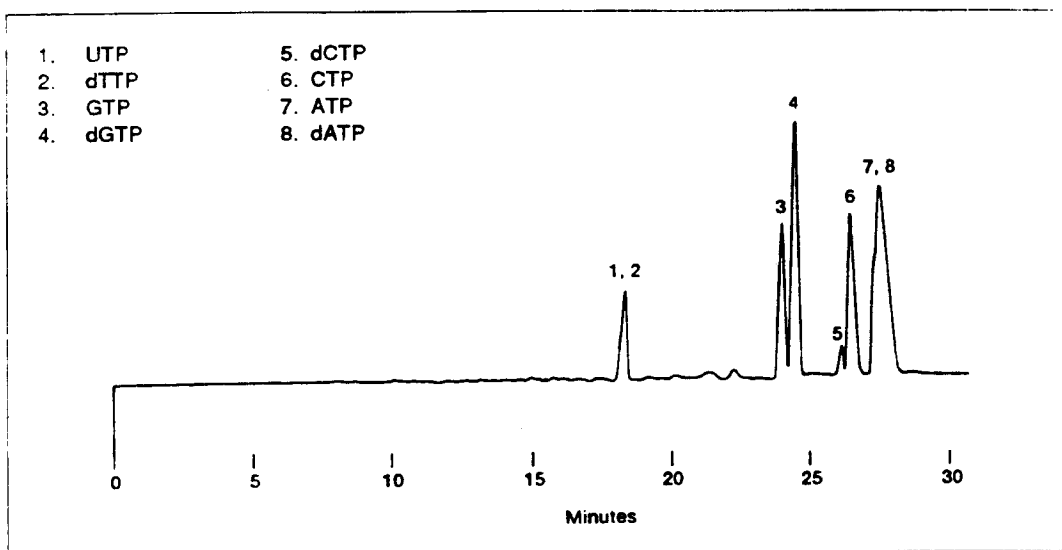


Single Nucleotides

Application
Note: 14

Separation of nucleotides in mixtures provides a means of looking at nucleotide pools as a way to characterize the physiological state of a cell. The separation shown here demonstrates nucleotides separated by capillary electrophoresis with visualization at 254 nm.



Capillary: 50 cm x 50 μ m, coated
Buffer: 0.1 M sodium phosphate, pH 2.5, 0.002 M EDTA
Load conditions: 8 kV, 8 seconds
Run conditions: 8 kV, constant voltage, $\ominus \rightarrow \oplus$ polarity
Detection: UV, 254 nm, 0.03 AUFS

1575-14 1290 90-0291