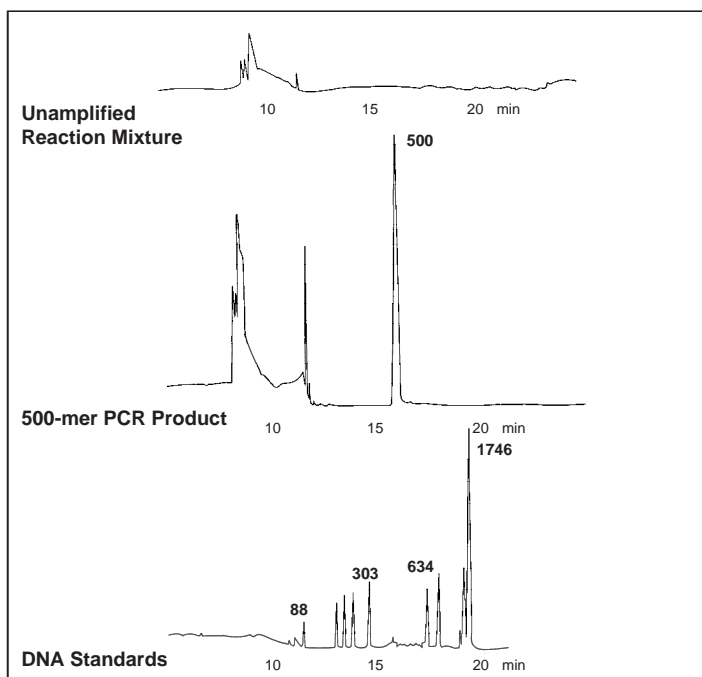


Non-Gel Sieving of PCR Products

The polymerase chain reaction (PCR), a rapid method for amplifying selected DNA sequences, has become a powerful tool in molecular biology. PCR products can range from several base pairs to kilobase lengths. Non-gel sieving by capillary electrophoresis provides a method for rapid analysis of PCR products. Analysis time for PCR products is typically less than 30 minutes, and on-tube UV detection permits quantitation. The figure below shows a 500 bp PCR product with a migration time corresponding to those of the DNA standards. Quantitative information on the yield of reaction products and side products should be useful in optimizing reaction conditions.



Capillary:	50 cm x 50 μ m, coated
Buffer:	0.1 M Tris borate, pH 8.3, 0.002 M EDTA, linear polymers
Load conditions:	8 kV, 8 seconds
Run conditions:	8 kV, constant voltage, negative polarity
Detection:	UV, 260 nm, 0.05 AUFS