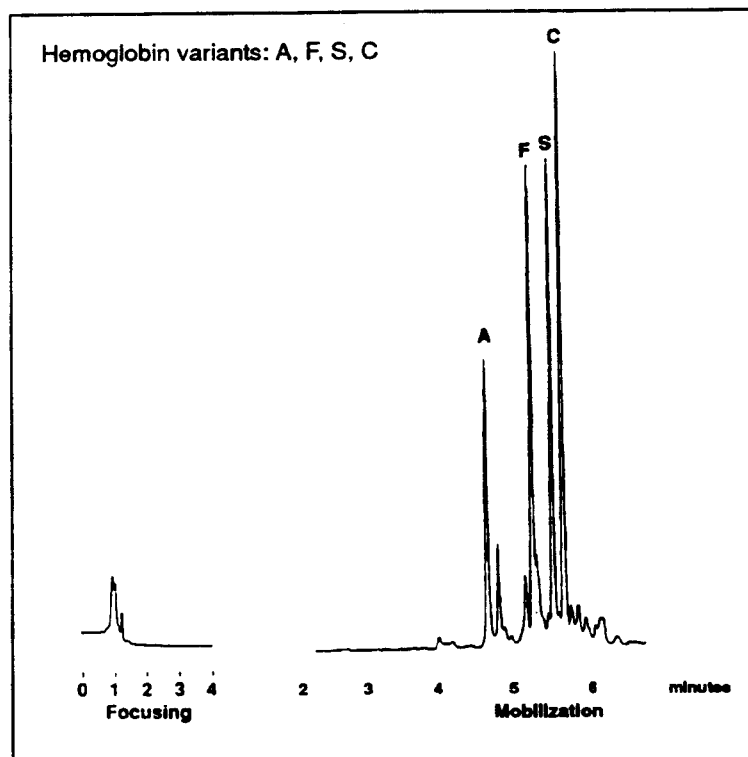


Isoelectric Focusing of Hemoglobin Variants

Application
Note: 2

High resolution separation of proteins based on their isoelectric points can be achieved using isoelectric focusing in coated capillaries. By changing the composition of the anolyte or catholyte solutions, the focused proteins can be mobilized through the detector to generate an electropherogram. The resolving power of high performance isoelectric focusing is demonstrated in this separation of hemoglobin variants A, F, S, and C.



Capillary:	12 cm x 25 μ m, coated
Buffer:	Bio-Lyte® ampholytes, pH 3-10
	Catholyte: 0.01 M phosphoric acid
	Anolyte: 0.02 M NaOH
	Mobilizer: 0.08 M NaCl and 0.02 M NaOH
Focusing conditions:	8 kV, constant voltage, $\oplus \rightarrow \ominus$ polarity
Mobilization:	8 kV, constant voltage, $\oplus \rightarrow \ominus$ polarity
Detection:	UV, 280 nm, 0.02 AUFS

1575-02 1290 90-0291