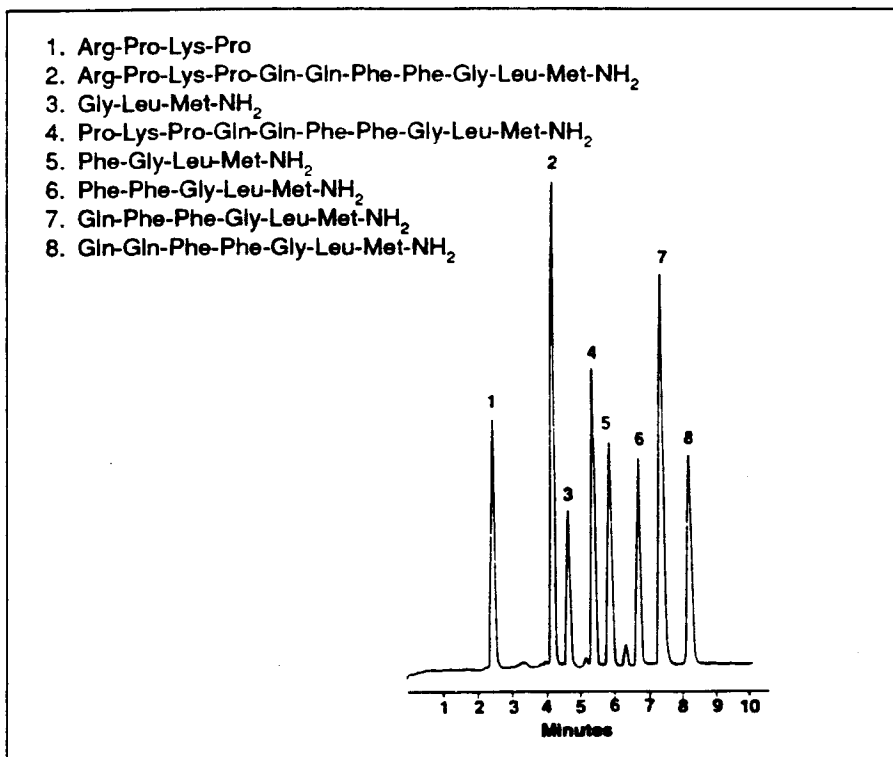


## Separation of Substance P and its Fragments

Application  
Note: 8

Substance P is an undecapeptide present in the brain and intestine. It acts as a neurotransmitter. The separation here was used to develop a standard curve to be used in the prediction of retention times for other proteins. For additional information, refer to Bulletin 1482.



<b>Capillary:</b>	20 cm x 25 $\mu$ m, coated
<b>Buffer:</b>	0.1 M sodium phosphate, pH 2.5
<b>Load conditions:</b>	8 kV, 8 seconds
<b>Run conditions:</b>	8 kV, constant voltage, $\oplus \rightarrow \ominus$ polarity
<b>Detection:</b>	UV, 200 nm, 0.02 AUFS