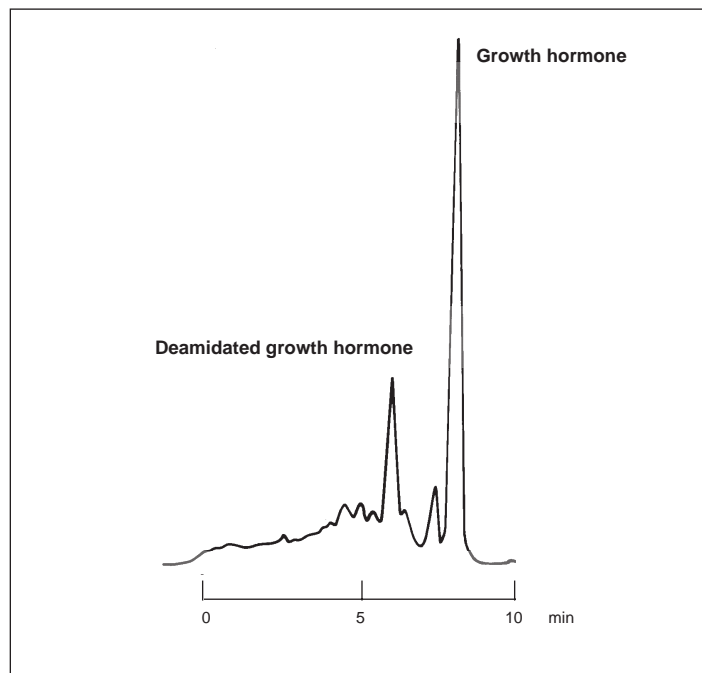


Measuring Deamidation of Human Growth Hormone

Capillary electrophoresis has captured the interest of the biopharmaceutical industry as an analytical tool for the quality control of polypeptide-based products. Here human growth hormone (HGH), extracted from pituitary tissue and subjected to elevated temperature and pH, is analyzed, revealing the presence of deamidated product. Natural degradation due to product aging can easily be monitored with capillary electrophoresis as shown by the separation below.



Capillary:	20 cm x 25 μ m, coated
Buffer:	0.05 M phosphate, pH 8.0, linear polymers, nonionic surfactant
Load conditions:	8 kV, 8 seconds
Run conditions:	8 kV, constant voltage, negative polarity
Detection:	UV, 200 nm, 0.05 AUFS