

Exenatide and Glu13-exenatide

Summary

Exenatide is a 39-amino-acid glucagon-like peptide-1 agonist used in the treatment of type II diabetes.

USP catalogue numbers

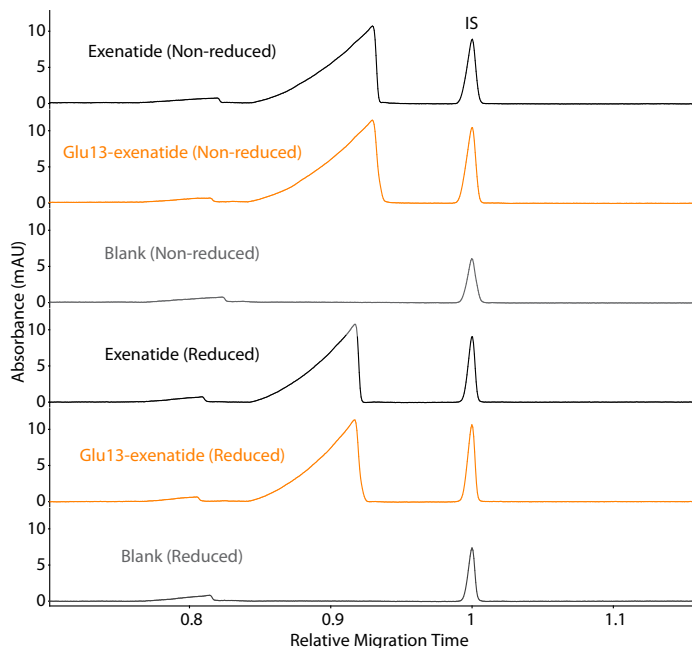
Exenatide: 1269105, Lot F052N0

Glu-13 exenatide: 1268998, Lot F04770

Maurice CE-SDS Method

Sample preparation: Exenatide and Glu13-exenatide were diluted to 0.5 mg/mL with 1X Sample Buffer prior to treatment for 5 minutes at 70 °C in the presence of either 11.5 mM IAM (non-reduced) or 650 mM β ME (reduced).

Running conditions: Samples were injected for 20 seconds, followed by a 15-minute separation at 4600 V.



Maurice icIEF method

Sample preparation: Exenatides were diluted to 0.05 mg/mL.

Ampholyte solution: Pharmalytes 3–10 (4%) containing 12.5 mM arginine, 12.5 mM IDA and 3.2 M urea

Running conditions: 1 minute at 1500 V, 6 minutes at 3000 V

pI markers: 4.05 and 9.99

Imaging: Absorbance and fluorescence (15 seconds)

