

Teriparatide

Summary

Teriparatide is a recombinant form of parathyroid hormone. Teriparatide is a 34-amino-acid bioactive protein. It is used in the treatment of some forms of osteoporosis.

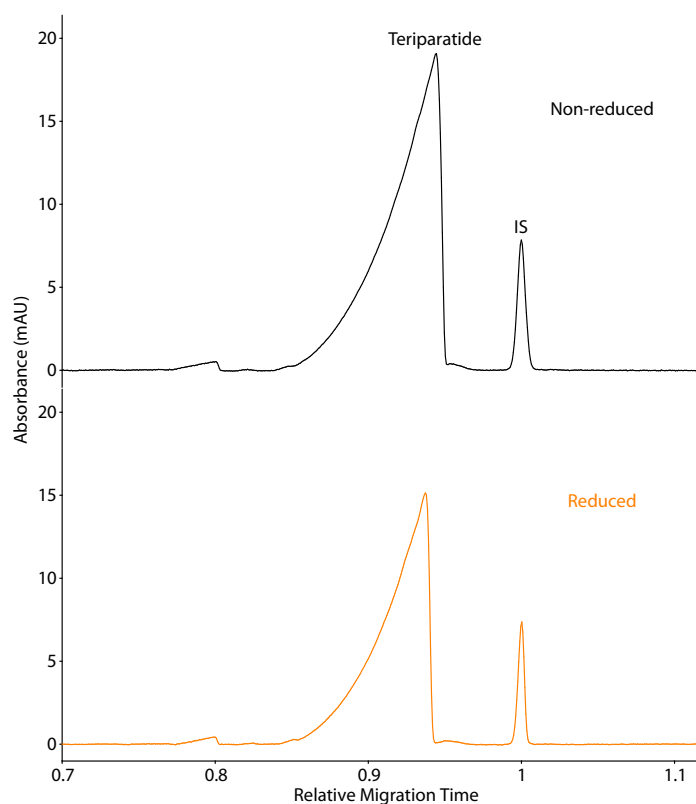
EDQM catalogue number

Teriparatide: Y0001916, Batch 1

Maurice CE-SDS Method

Sample preparation: Teriparatide was diluted to 1 mg/mL with 1X Sample Buffer prior to treatment for 10 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 8 seconds at 4600 V, followed by a 17-minute separation at 5750 V.



Maurice icIEF method

Sample preparation: Teriparatide was diluted to 0.065 mg/mL.

Ampholyte solution: Pharmalytes 3–10 (4%) containing 20 mM arginine and 8 M urea

Running conditions: 1 minute at 1500 V, 15 minutes at 3000 V

pI markers: 7.05 and 10.17

Imaging: Absorbance and fluorescence (10 seconds)

