

# Glucagon

## Summary

Glucagon is a 29-amino-acid non-steroidal peptide hormone used in the treatment of acute hypoglycemia.

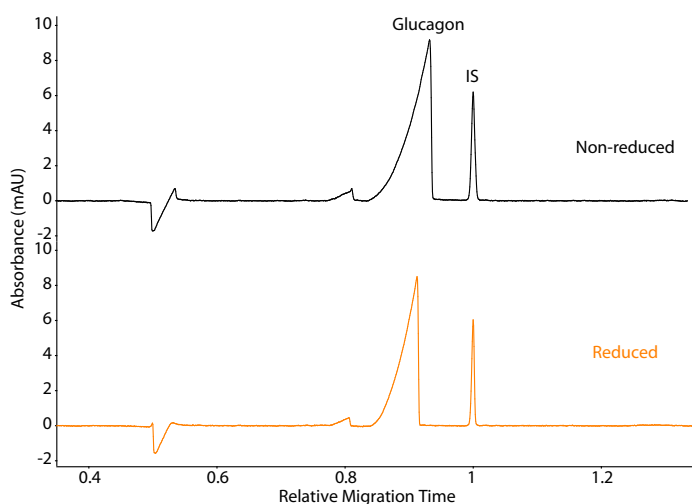
## EDQM catalogue number

Glucagon: Y0000191, Batch 2

## Maurice CE-SDS Method

**Sample preparation:** Glucagon was diluted to 1.0 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  $\beta$ ME (reduced).

**Running conditions:** Samples were injected for 8 seconds, followed by a 20-minute separation at 5750 V.



## Maurice icIEF method

**Sample preparation:** Glucagon was diluted to 0.06 mg/mL.

**Ampholyte solution:** Pharmalytes 3–10 (4%) containing 8 M urea

**Running conditions:** 1 minute at 1500 V, 10 minutes at 3000 V

**pI markers:** 5.85 and 7.05

**Imaging:** Absorbance and fluorescence (5 seconds)

