

Vasopressin and Lypressin

Summary

Vasopressin is known as the anti-diuretic hormone as it is responsible for regulating the retention of water in the body.

USP catalogue numbers

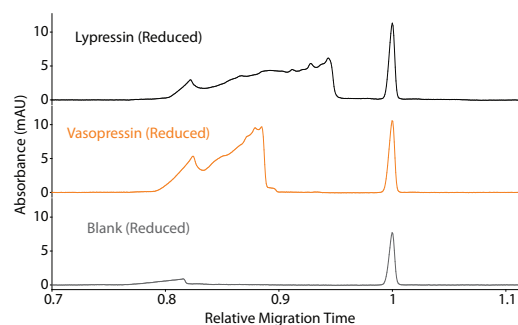
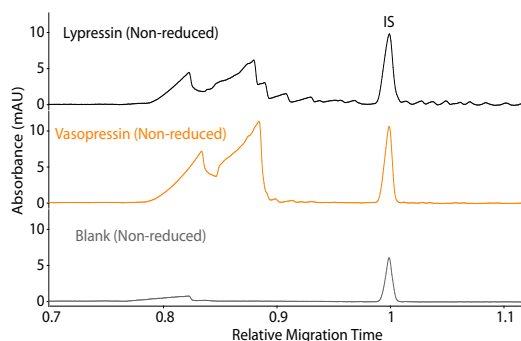
Lypressin: 1711100, Lot H0L444

Vasopressin: 1342321, Lot R016U0

Maurice CE-SDS Method

Sample preparation: Vasopressin and lypressin (purchased from USP) were diluted to 0.5 mg/mL in 1X Maurice 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).

Run conditions: Non-reduced vasopressin was injected for 20 seconds followed by a 15-minute separation at 4600 V. Reduced vasopressin was injected for 10 seconds followed by a 15-minute separation at 4600 V. Non-reduced lypressin was injected for 10 seconds followed by a 20-minute separation at 4600 V. Reduced lypressin was injected for 10 seconds followed by a 20-minute separation at 4600 V.



Maurice icIEF method

Sample preparation: Vasopressin and lypressin were diluted to 0.1 mg/mL.

Ampholyte solution: Pharmalytes 8–10.5 (2.25%) and Servalytes 9–11 (0.75%) containing 3.2 M urea and 1% TEMED

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

pI marker: 8.4

Imaging: Absorbance

