

Product Data Sheet

Product Name: Spexin-2 (53-70), human/mouse/rat

Catalog Number: AS-65586 (1 mg) Lot Number: See label on vial

Sequence: Phe-Ile-Ser-Asp-Gln-Ser-Arg-Lys-Asp-Leu-Ser-Asp-Arg-Pro-Leu-Pro-

Glu-OH

FISDQSRRKDLSDRPLPE

Molecular Weight: 2159.5

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This peptide is shipped at ambient temperature. Upon receipt, store

lyophilized peptide at -20°C or lower. Reconstituted peptide can be

aliquoted and stored at -20°C or lower.

Description: This is Spexin-2 (53-70), non-amidated version of the novel peptide

hormone that is derived from prohormone proNPQ and is conserved

amongst mammalian species. It is a central modulator of cardiovascular and renal function. Spexin-2, when administered to rats, decreases heart rate

and increases urine flow rate.

Related Products: Spexin (36-49), amidated Cat# AS-65569

Spexin-2 (53-70), amidated Cat# AS-65570

References:

1. Gu L. et al. *Peptides* **71**, 232-239 (2015).

2. Toll L. et al. *FASEB* **26**, 947-954 (2012).

For Research Use Only