

Product Data Sheet

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

Catalog Number: AS-65570 (1 mg) Lot Number: See label on vial

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

Glu-NH2

FISDQSRRKDLSDRPLPE-NH2

Molecular Weight: 2158.9

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 amidated Spexin-2 (53-70) peptide showing similar biological

function as that of its non- amidated version (Cat # 65586) as a central modulator of cardiovascular and renal function. Spexin-2 is a novel peptide hormone that is derived from prohormone proNPQ and is conserved

amongst mammalian species.

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

Spexin-2 (53-70) Cat# AS-65586

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