

Product Data Sheet

Product Name: Spexin (36-49), amidated, human/mouse/rat

Catalog Number: AS-65569 (1 mg) Lot Number: See label on vial

Sequence: Asn-Trp-Thr-Pro-Gln-Ala-Met-Leu-Tyr-Leu-Lys-Gly-Ala-Gln-NH2

NWTPQAMLYLKGAQ-NH2

Molecular Weight: 1620

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: Spexin is a novel peptide hormone that is derived from prohormone proNPQ

and is conserved amongst mammalian species. Spexin (36-49) is post-translationally amidated with a C-terminal amidation processing signal motif (GRR) following Q49 residue of the pro-hormone. Spexin reduces adipocyte uptake of long chain fatty acids and causes weight loss in rodents. Recently

it was suggested that this peptide may play a role in lipid and glucose

metabolism in patients with type-2 diabetes mellitus.

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

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).

3. Walewski JL., et.al., Obesity 22 (7), 1643-1652 (2014)

For Research Use Only