

Product Data Sheet

Product Name: Neuropeptide S, NPS, mouse

Catalog Number: AS-61246 (1 mg) Lot Number: See label on vial

Sequence: H-Ser-Phe-Arg-Asn-Gly-Val-Gly-Ser-Gly-Ala-Lys-Lys-Thr-Ser-Phe-Arg-

Arg-Ala-Lys-Gln-OH (3-letter code)

SFRNGVGSGAKKTSFRRAKQ (1-letter code)

Molecular Weight: 2183.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Neuropeptide S is freely soluble in H₂O.

Storage: Neuropeptide S 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: A neuropeptide that plays a role in arousal and fear responses. Ref: Xu, Y. et al. *Neuron* **43**, 487 (2004).

Related Products:

Name Cat # Size

Neuropeptide S, NPS, human
(SFRNGVGTGMKKTSFQRAKS)

AS-61243 1 mg

For Research Use Only