



Product Data Sheet

Product Name: Neuropeptide S, NPS, human

Catalog Number: AS-61243 (1 mg) **Lot Number:** See label on vial

Sequence: H-Ser-Phe-Arg-Asn-Gly-Val-Gly-Thr-Gly-Met-Lys-Lys-Thr-Ser-Phe-Gln-Arg-Ala-Lys-Ser-OH (3-letter code)
SFRNGVGTGMKKT SFQRAKS (1-letter code)

Molecular Weight: 2188.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 µl to 1 mg Neuropeptide S.

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, mouse (SFRNGVGSGAKKTSFRRAKQ)	AS-61246	1 mg

For Research Use Only