



Product Data Sheet

Product Name: Neuropeptide S, NPS, human

Catalog Number: 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

Peptide Purity: >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 (SFRNGVGS GAKKTSFRR AKQ)	61246	1 mg
Neuropeptide S, NPS, chimpanzee, dog (SFRNGVGTGMKKT SFRR AKS)	62617	1 mg
Neuropeptide S, NPS, rat (SFRNGVGS GVKKTSFRR AKQ)	61249	1 mg
Neuropeptide S, NPS, chicken (SFRNGVGS GIKKTSFRR AKS)	62618	1 mg
Neuropeptide S, hNPS (1-12), human (SFRNGVGTGMKK)	62608	1 mg
Neuropeptide S, NPS (1-10), rat (SFRNGVGS GV)	62599	1 mg

For Research Use Only