



Product Data Sheet

Product Name:	Neuropeptide Y, human, rat	
Catalog Number:	AS-22464 (0.5 mg) AS-22465 (1 mg)	Lot Number: See label on vial
Sequence:	H-Tyr-Pro-Ser-Lys-Pro-Asp-Asn-Pro-Gly-Glu-Asp-Ala-Pro-Ala-Glu-Asp-Met-Ala-Arg-Tyr-Tyr-Ser-Ala-Leu-Arg-His-Tyr-Ile-Asn-Leu-Ile-Thr-Arg-Gln-Arg-Tyr-NH ₂ (3-letter code) YPSKPDNPGEDAPAEDMARYYSALRHYINLITRQRY-NH ₂ (1-letter code)	
Molecular Weight:	4271.8	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Neuropeptide Y is freely soluble in water.

Storage: Neuropeptide Y 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.

Reference: Minth, C. et al. *Proc. Natl. Acad. Sci. USA* **81**, 4577 (1984).

Description: Human NPY, rPP, D-Trp³⁴ NPY, D-Arg²⁵ NPY, C2-NPY, and cPP NPY were synthesized by AnaSpec (San Jose, CA). All peptides were dissolved in 0.9% sterile saline (Sigma, St. Louis, MO) and infused icv at 1 nmol, which was the minimal effective dose of D-Trp³⁴ NPY and D-Arg²⁵ NPY on feeding at the 1-h time point-Gao, J. et al. *Am J Physiol Endocrinol Metab* **287**, E706 (2004).

The doses of intra-hippocampal injection of NPY (10 $\mu\text{g}/\mu\text{l}$, Anaspec Company) and 5-HT (31.25 $\mu\text{g}/\mu\text{l}$, Sigma Company) were chosen according to previous studies. On days 1, 7, 14 and 21, rats received microinjection of 1 μl of 0.9% physiological saline for control and CUMS groups, 1 μl of NPY for NPY group, and 1 μl of 5-HT for 5-HT group. Each intra-hippocampal microinjection was made before stressing. The injection was manually delivered at the speed of 1 $\mu\text{l}/\text{min}$ -Luo, DD. et al. *Brain Res. Bulletin* **77**, 8 (2008).

Published Citations:

Kapoor, JR. and CD. Sladek. *Am J Physiol Regul Integr Comp Physiol* **280**, R69 (2001).
Taiwo, OB and BK, Taylor. *Pain* **96**, 353 (2002).
Gao, J. et al. *Am J Physiol Endocrinol Metab* **287**, E706 (2004).
Saurer, TB. et al. *J. of Neuroimmunol.* **177**, 18 (2006).
Pelz, KM. & J. Dark *Am J Physiol Regulatory Integrative Comp Physiol* **292**, R2299 (2007).

Taylor, BK. et al. *Pain* **131**, 83 (2007).
Intondi, AB. et al. *Pain* **137**, 352 (2008).
Luo, DD. et al. *Brain Res. Bullet.* **77**, 8 (2008).
Mullins, D. et al. *Euro. J. of Pharmacol.* **601**, 1 (2008)

For Research Use Only