



Product Data Sheet

Product Name:	[Leu31, Pro34]-Neuropeptide Y, human, rat	
Catalog Number:	AS-24366 (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-Leu-Thr-Arg-Pro-Arg-Tyr-NH ₂ (3-letter code) YPSKPDNPGEDAPAEDMARYYSALRHYINLLTRPRY-NH ₂ (1-letter code)	
Molecular Weight:	4240.8	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: [Leu31, Pro34]-Neuropeptide Y is freely soluble in water.

Storage: [Leu31, Pro34]-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.

Description: Like neuropeptide Y (NPY) and other peptides of the family, this peptide adopts a folded hairpin structure with the terminal segments in close proximity. An intracerebroventricular injection of NPY or [Leu31, Pro34]-NPY (non-Y₂ receptor agonist) given during middle cerebral artery occlusion increases the infarct volume and nitric oxide (NO) overproduction in the rat brain. Ref: Futilendorf, I. et al. *Proc. Natl. Acad. Sci. USA* **87**, 182 (1990). Khiat, AJ. *Pept. Res.* **51**, 317 (1998); Chen, S. and R.J. Cheung, *Biomed. Sci.* **12**, 267 (2005); Chen, S. and R.J. Cheung, *Biomed. Sci.* **11**, 781 (2004); Labelle, M. et al. *Peptides* **18**, 801 (1997)

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

On the day of the experiment, the animals were brought to the experiment room in which they acclimated to the surroundings. Two hours before the start of the experiment, 33-gauge internal injectors were inserted into the intracerebral cannula, which extended 1 mm beyond the tip of the guide cannula. A cannula connector was used to attach the injectors to a syringe pump for delivery of equal volumes of 0.5 μl vehicle (0.1% BSA in 0.9% saline) or 110 pmol (250 ng) of [Leu³¹pro³⁴]NPY (AnaSpec, San Jose, CA) into the PVN over a period of 5 min-[Dimitrov, EL. et al. *Endocrinology* **148**, 3666 \(2007\).](#)

Published Citations:

For Research Use Only

Kapoor, JR. and CD. Sladek, *Am J Physiol Regul Integr Comp Physiol* **280**, R69 (2001).
Dimitrov, EL. et al. *Endocrinology* **148**, 3666 (2007).

