

PRODUCT INFORMATION



Neuropeptide Y (1-24) (human, rat) (trifluoroacetate salt)

Item No. 24715

Synonym:	NPY (1-24) (human, rat)	
MF:	$C_{116}H_{170}N_{30}O_{40}S \cdot XCF_3COOH$	H-Tyr-Pro-Ser-Lys-Pro-Asp-Asn-Pro-Gly-Glu-
FW:	2,656.8	Asp-Ala-Pro-Ala-Glu-Asp-Met-Ala-Arg-Tyr-
Purity:	≥95%	Tyr-Ser-Ala-Leu-NH ₂
Supplied as:	A lyophilized powder	• XCF ₃ COOH
Storage:	-20°C	
Stability:	≥4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Neuropeptide Y (NPY) (1-24) (human, rat) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the NPY (1-24) (human, rat) (trifluoroacetate salt) in water. The solubility of NPY (1-24) (human, rat) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

NPY (1-24) is an N-terminal fragment of NPY, a neuropeptide involved in controlling appetite, blood pressure, cardiac contractility, and intestinal secretion.¹ NPY (1-24) does not inhibit [³H]-NPY binding in rat CNS membranes at concentrations up to 10 μM and has no effect on the twitch response of isolated rat vas deferens either alone (EC₅₀ = >1 μM) or when used in combination with the C-terminal fragment NPY (25-36) at a concentration of 1 μM.^{2,3} NPY (1-24) stimulates histamine release from isolated rat peritoneal mast cells with a lower efficacy than full length NPY.⁴

References

1. Balasubramaniam, A. Neuropeptide Y family of hormones: Receptor subtypes and antagonists. *Peptides* **18(3)**, 445-457 (1997).
2. Martel, J.-C., Fournier, A., St-Pierre, S., *et al.* Comparative structural requirements of brain neuropeptide Y binding sites and vas deferens neuropeptide Y receptors. *Mol. Pharmacol.* **38(4)**, 494-502 (1990).
3. Kahl, U., Langel, Ü., Bartfai, T., *et al.* Functional effects and ligand binding of chimeric galanin-neuropeptide Y (NPY) peptides on NPY and galanin receptor types. *Br. J. Pharmacol.* **111(4)**, 1129-1134 (1994).
4. Grundemar, L., Krstenansky, J.L., and Håkanson, R. Neuropeptide Y and truncated neuropeptide Y analogs evoke histamine release from rat peritoneal mast cells. A direct effect on G proteins? *Eur. J. Pharmacol.* **258(1-2)**, 163-166 (1994).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/17/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM