PRODUCT INFORMATION



Tyr —

Neuropeptide W-23 (human) (trifluoroacetate salt)

Item No. 24551

Formal Name:	L-tryptophyl-L-tyrosyl-L-lysyl-L-histidyl-L-valyl-L-alanyl- L-seryl-L-prolyl-L-arginyl-L-tyrosyl-L-histidyl-L-threonyl-	
	L-valylglycyl-L-arginyl-L-alanyl-L-alanylglycyl-L-leucyl-L- leucyl-L-methionylglycyl-L-leucine, trifluoroacetate salt	H-Trp-Tyr-Lys-His-Val-Ala-Ser-Pro-Arg-Tyr-
Synonym:	NPW23 (human)	His—Thr—Val—Gly—Arg—Ala—Ala—Gly—Leu—Leu—
MF:	C110H183N35O28S • XCF3COOH	Met_Glv_Leu_OH
FW:	2,584.0	Wet ony Lou off
Purity:	≥95%	• XCF ₃ COOH
Supplied as:	A lyophilized powder	
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 W-23 (NPW23) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the NPW23 (human) (trifluoroacetate salt) in water. The solubility of NPW23 (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

NPW23 is an endogenous neuropeptide ligand for the orphan G protein-coupled receptors GPR7 and GPR8 (K_i s = 31.8 and 20.7 pM, respectively).¹ It inhibits forskolin-induced cAMP accumulation in CHO cells expressing human GPR7 and GPR8 with IC₅₀ values of 0.025 and 0.178 nM, respectively. NPW23 (3 nmol) stimulates prolactin release in rats but has no effect on other pituitary hormones. NPW23 attenuates mechanical allodynia, but not thermal hyperalgesia, induced by carrageenan injection in rat paw in a dose-dependent manner.² It also decreases food intake and weight gain as well as increases body temperature and thermogenesis in rats.³

References

- 1. Shimomura, Y., Harada, M., Goto, M., et al. Identification of neuropeptide W as the endogenous ligand for orphan G-protein-coupled receptors GPR7 and GPR8. J. Biol. Chem. 277(39), 35826-35832 (2002).
- 2. Yamamoto, T., Saito, O., Shono, K., et al. Anti-hyperalgesic effects of intrathecally administered neuropeptide W-23, and neuropeptide B, in tests of inflammatory pain in rats. Brain Res. 1045(1-2), 97-106 (2005).
- 3. Naso, T., Shousha, S., and El-Kirdasy, A. Central neuropeptide W has anorexigenic effect in rats. J. Anim. Physiol. Anim. Nutr. (Berl.) 98(2), 228-234 (2014).

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

SAFFTY 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

uyer 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, 10/06/2022

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