

PRODUCT INFORMATION



[Orn⁸]-Urotensin II (human) (trifluoroacetate salt)

Item No. 24754

Formal Name: L- α -glutamyl-L-threonyl-L-prolyl-L- α -aspartyl-L-cysteinyl-L-phenylalanyl-L-tryptophyl-L-ornithyl-L-tyrosyl-L-cysteinyl-L-valine, cyclic (5 \rightarrow 10)-disulfide, trifluoroacetate salt

MF: C₆₃H₈₃N₁₃O₁₈S₂ • XCF₃COOH

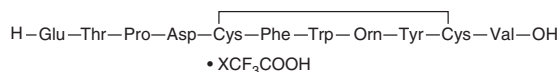
FW: 1,374.5

Purity: \geq 95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

[Orn⁸]-Urotensin II (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the [Orn⁸]-urotensin II (human) (trifluoroacetate salt) in water. The solubility of [Orn⁸]-urotensin II (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

[Orn⁸]-Urotensin II is a peptide analog of urotensin II (Item No. 24711) and agonist of the urotensin II receptor (UTR; EC₅₀s = 11.74 and 8.7 nM in HEK293 cells expressing the human and rat receptor, respectively).¹ It reduces urotensin II-induced contraction of isolated rat aorta (pA₂ = 6.56) but has no effect on aortic contraction when used alone at concentrations up to 10 μ M.

Reference

1. Camarda, V., Guerrini, R., Kostenis, E., *et al.* A new ligand for the urotensin II receptor. *Br. J. Pharmacol.* **137**(3), 311-314 (2002).

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, 11/17/2022

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