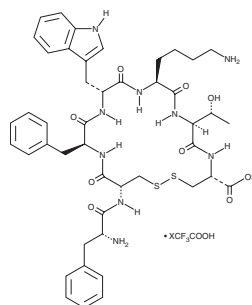


PRODUCT INFORMATION

Des-Threoninol-Octreotide (trifluoroacetate salt)

Item No. 42545

Formal Name: D-phenylalanyl-L-cysteinyl-L-phenylalanyl-D-tryptophyl-L-lysyl-L-threonyl-L-cysteine, cyclic (2→7)-disulfide, trifluoroacetate salt
Synonym: cyclo(D-Phe-Cys-Phe-D-Trp-Lys-Thr-Cys)
Peptide Sequence: cyclo(fCFwKTC)
MF: $C_{45}H_{57}N_9O_9S_2 \cdot XCF_3COOH$
FW: 932.1
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Des-threoninol-octreotide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the des-threoninol-octreotide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Des-threoninol-octreotide (trifluoroacetate salt) is soluble (≥10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of des-threoninol-octreotide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Des-threoninol-octreotide (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Des-threoninol-octreotide is a peptide ligand of somatostatin receptors (SSTs).¹ It binds Sst_{2B}, Sst₅, and SST₅ (IC₅₀s = 13.62, 10.63, and 10.95 nM for the mouse, rat, and human receptors, respectively).

Reference

- Ösapay, G., Prokai, L., Kim, H.S., *et al.* Lanthionine-somatostatin analogs: Synthesis, characterization, biological activity, and enzymatic stability studies. *J. Med. Chem.* **40**(14), 2241-2251 (1997).

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, 02/24/2025

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