

PRODUCT INFORMATION



CH 275 (acetate)

Item No. 41050

Formal Name: L-cysteinyl-L-lysyl-L-phenylalanyl-L-phenylalanyl-D-tryptophyl-4-[[[(1-methylethyl)amino]methyl]-L-phenylalanyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L-cysteine, cyclic (1→11)-disulfide, acetate

Peptide Sequence: CLFFwXTFTSC-OH (X= 4-(N-isopropyl)-aminomethylphenylalanine)

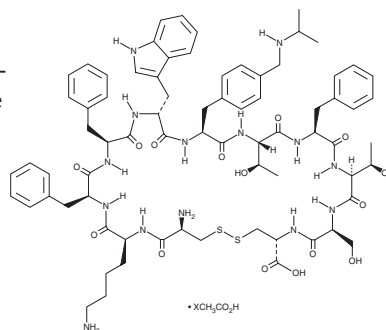
MF: C₇₄H₉₆N₁₄O₁₅S₂ • XC₂H₄O₂
FW: 1,485.8

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

CH 275 (acetate) is supplied as a solid. A stock solution may be made by dissolving the CH 275 (acetate) in the solvent of choice, which should be purged with an inert gas. CH 275 (acetate) is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

CH 275 (acetate) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

CH 275 is a cyclic peptide agonist of somatostatin receptor 1 (SST₁; IC₅₀ = 30.9 nM for the human receptor).¹ It is selective for SST₁ over SST₂, SST₃, SST₄, and SST₅ (IC₅₀s = >10,000, 345, >1,000, and >10,000 nM, respectively, for the human receptors). CH 275 (500 nM) inhibits basal growth hormone secretion from primary mouse pituitary cells.² It inhibits glucose-induced increases in electrical activity without inducing membrane hyperpolarization in a whole-cell patch-clamp assay using MIN6 pancreatic β-cells when used at a concentration of 100 nM.³

References

1. Rivier, J.E., Hoeger, C., Erchegyi, J., *et al.* Potent somatostatin undecapeptide agonists selective for somatostatin receptor 1 (sst1). *J. Med. Chem.* **44**(13), 2238-2246 (2001).
2. Kreienkamp, H.J., Akgün, E., Baumeister, H., *et al.* Somatostatin receptor subtype 1 modulates basal inhibition of growth hormone release in somatotrophs. *FEBS Lett.* **462**(3), 464-466 (1999).
3. Smith, P.A., Sellers, L.A., and Humphrey, P.P. Somatostatin activates two types of inwardly rectifying K⁺ channels in MIN-6 cells. *J. Physiol.* **532**(Pt 1), 127-142 (2001).

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, 07/15/2024

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