

PRODUCT INFORMATION



CYN 154806 (trifluoroacetate salt)

Item No. 37343

Synonym: Ac-4NO₂-Phe-c(D-Cys-Tyr-D-Trp-Lys-Thr-Cys)-D-Tyr-NH₂

Peptide Sequence: Ac-4-NO₂-F-c(YwKTC)-y-NH₂

MF: C₅₆H₆₈N₁₂O₁₄S₂ • XCF₃COOH

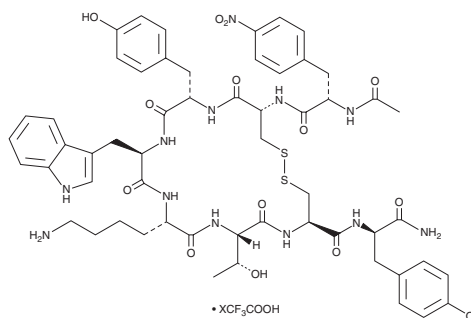
FW: 1,197.3

Purity: ≥90%

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

CYN 154806 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CYN 154806 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CYN 154806 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CYN 154806 (trifluoroacetate salt) in DMSO is approximately 10 mg/ml and approximately 25 mg/ml in ethanol and DMF.

Description

CYN 154806 is a cyclic peptide somatostatin receptor 2 (SST₂) antagonist (IC₅₀ = 2.62 nM).¹ It is selective for SST₂ over SST₁, SST₃, SST₄, and SST₅ (IC₅₀s = 3,890, 851, 1,737, and 331 nM, respectively). CYN 154806 inhibits SST-induced increases in extracellular acidification (EAR) in CHO-K1 cells expressing recombinant human SST₂ (pK_B = 7.92) and [³⁵S]GTPγS binding in CHO-K1 cell membranes expressing recombinant human SST₂ (pK_B = 7.81).

Reference

1. Feniuk, W., Jarvie, E., Luo, J., *et al.* Selective somatostatin sst₂ receptor blockade with the novel cyclic octapeptide, CYN-154806. *Neuropharmacology* **39**(8), 1443-1450 (2000).

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, 11/01/2023

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