

PRODUCT INFORMATION

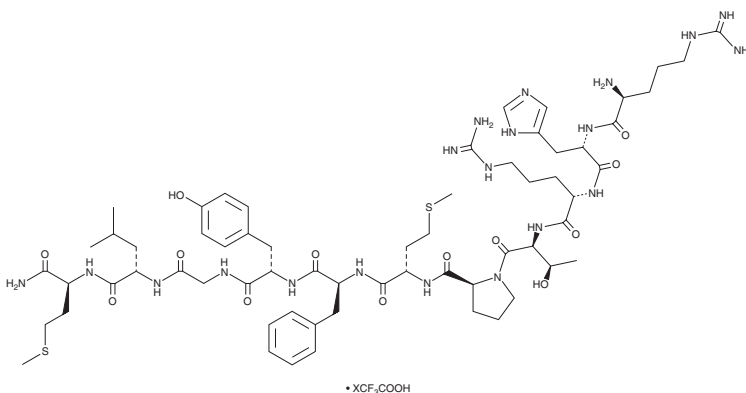


RHRTPMFYGLM-NH₂ (trifluoroacetate salt)

Item No. 43089

Formal Name: (S)-1-(L-arginyl-L-histidyl-L-arginyl-L-threonyl)-N-((S)-1-(((S)-1-(((S)-1-(((S)-1-(((S)-1-(((S)-1-amino-4-(methylthio)-1-oxobutan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)amino)-2-oxoethyl)amino)-3-(4-hydroxyphenyl)-1-oxopropan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)amino)-4-(methylthio)-1-oxobutan-2-yl)pyrrolidine-2-carboxamide, trifluoroacetate salt

Peptide Sequence: RHRTPMFYGLM-NH₂
MF: C₆₃H₉₈N₂₀O₁₃S₂ • XCF₃COOH
FW: 1,407.7
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

RHRTPMFYGLM-NH₂ (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the RHRTPMFYGLM-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. RHRTPMFYGLM-NH₂ (trifluoroacetate salt) is soluble (≥10 mg/ml) in 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 RHRTPMFYGLM-NH₂ (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. RHRTPMFYGLM-NH₂ (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

RHRTPMFYGLM-NH₂ is a derivative of the endogenous neurokinin-1 (NK₁) receptor agonist C14TKL-1 that contains leucine at position 10 instead of tyrosine.^{1,2} It contains the typical tachykinin motif FXGLM-NH₂.

References

1. Jiang, Y., Gao, G., Fang, G., *et al.* PepPat, a pattern-based oligopeptide homology search method and the identification of a novel tachykinin-like peptide. *Mamm. Genome* **14**(5), 341-349 (2003).
2. Page, N.M. New challenges in the study of the mammalian tachykinins. *Peptides* **26**(8), 1356-1368 (2005).

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, 05/16/2025

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