

PRODUCT INFORMATION



MMP Inhibitor I (trifluoroacetate salt)

Item No. 24708

Formal Name: (S)-1-((2-aminobenzoyl)glycyl)-N-((R)-1-(((R)-1-(hydroxyamino)-1-oxopropan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)pyrrolidine-2-carboxamide, trifluoroacetate salt

Synonyms: Abz-GP-DLeu-DAla-NH-OH, FN 439, Matrix Metalloproteinase Inhibitor I

MF: C₂₃H₃₅N₆O₆ • XCF₃COOH

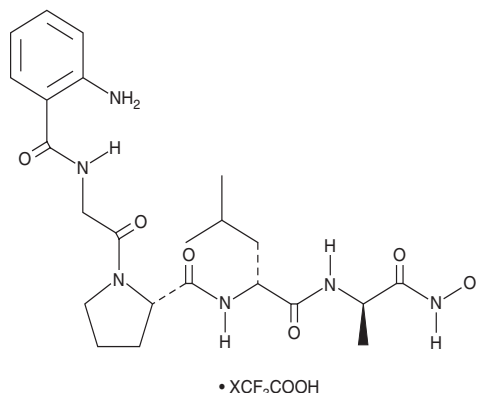
FW: 491.6

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

MMP inhibitor I (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the MMP inhibitor I (trifluoroacetate salt) in water. The solubility of MMP inhibitor I (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MMP inhibitor I is a peptide inhibitor of matrix metalloproteinase-1 (MMP-1), MMP-2, and MMP-3 (IC₅₀s = 1.3, 30, and 150 μM, respectively).¹ It is selective for MMP-1, MMP-2, and MMP-3 over the proteases thermolysin, urease, trypsin, α-chymotrypsin, plasmin, and elastase at concentrations up to 4 mM.

Reference

1. Otake, S., Morita, Y., Morikawa, T., *et al.* Inhibition of matrix metalloproteinases by peptidyl hydroxamic acids. *Biochem. Biophys. Res. Commun.* **199**(3), 1442-1446 (1994).

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