

PRODUCT INFORMATION



MMP-2/MMP-9 Inhibitor III (trifluoroacetate salt)

Item No. 24697

Formal Name: L-cysteinyl-L-threonyl-L-threonyl-L-histidyl-L-tryptophylglycyl-L-phenylalanyl-L-threonyl-L-leucyl-L-cysteine, cyclic (1→10)-disulfide, 2,2,2-trifluoroacetate

Synonym: Matrix Metalloproteinase-2/Matrix Metalloproteinase-9 Inhibitor III

MF: C₅₂H₇₁N₁₃O₁₄S₂ • XCF₃COOH

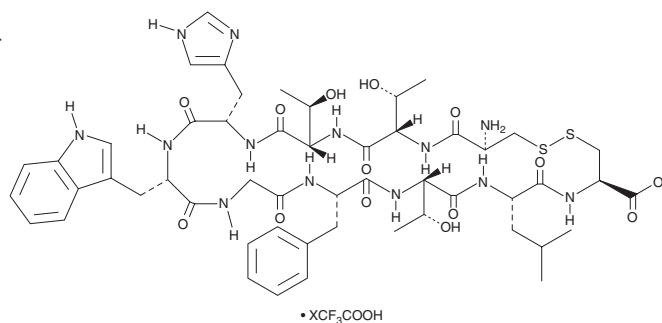
FW: 1,166.3

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-2/MMP-9 inhibitor III (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the MMP-2/MMP-9 inhibitor III (trifluoroacetate salt) in water. The solubility of MMP-2/MMP-9 inhibitor III (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MMP-2/MMP-9 inhibitor III is a cyclic peptide inhibitor of matrix metalloproteinase-2 (MMP-2) and MMP-9 (IC₅₀s = 10-20 μM for both).¹ It is selective for MMP-2 and MMP-9 over MT1-MMP, MMP-8, and MMP-13 as well as the serine proteases trypsin-2, neutrophil elastase, and cathepsin G when used at a concentration of 500 μM. MMP-2/MMP-9 inhibitor III reduces migration of HT1080 fibrosarcoma, C8161 melanoma, SKOV3 ovarian carcinoma, and KS1767 Kaposi's sarcoma cells *in vitro*. Pretreatment with MMP-2/MMP-9 inhibitor III delays tumor formation in an MDA-MB-435 mouse xenograft model. MMP-2/MMP-9 inhibitor III (100-200 μg per animal) also prevents growth of established tumors in MDA-MB-435 and KS1767 mouse xenograft models.

Reference

1. Koivunen, E., Arap, W., Valtanen, H., *et al.* Tumor targeting with a selective gelatinase inhibitor. *Nat. Biotechnol.* **17**(8), 768-774 (1999).

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, 03/09/2023

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