

PRODUCT INFORMATION



MMP-3 Inhibitor

Item No. 24315

CAS Registry No.: 158841-76-0

Formal Name: N²-acetyl-L-arginyl-L-cysteinylglycyl-L-valyl-L-prolyl-L- α -asparagine

Synonyms: Matrix Metalloproteinase-3 Inhibitor, Stromelysin-1 Inhibitor

Peptide Sequence: H-RCGVDPD-OH

MF: C₂₇H₄₆N₁₀O₉S

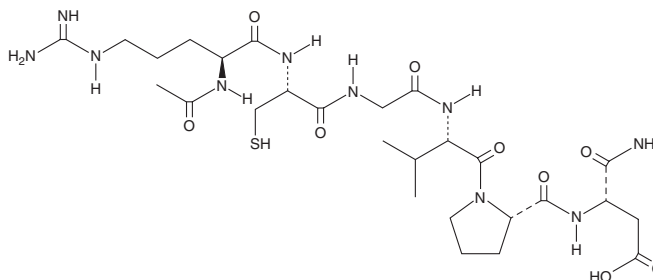
FW: 686.8

Purity: $\geq 95\%$

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

MMP-3 inhibitor is supplied as a solid. A stock solution may be made by dissolving the MMP-3 inhibitor in the solvent of choice, which should be purged with an inert gas. MMP-3 inhibitor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MMP-3 inhibitor in these solvents is approximately 30 mg/ml.

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 MMP-3 inhibitor can be prepared by directly dissolving the solid in aqueous buffers. The solubility of MMP-3 inhibitor in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MMP-3 inhibitor is a peptide inhibitor of matrix metalloproteinase-3 (MMP-3) with a K_i value of 95 nM.¹

Reference

1. Umezawa, N., Noro, Y., Ukai, K., *et al.* Photocontrol of peptide function: Backbone cyclization strategy with photocleavable amino acid. *ChemBioChem* **12**(11), 1694-1698 (2011).

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, 06/20/2023

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