

PRODUCT INFORMATION



MMP-3 Fluorogenic Substrate (trifluoroacetate salt)

Item No. 37523

Formal Name: N²-[2-(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]-L-arginyl-L-prolyl-L-lysyl-L-prolyl-L-valyl-L-α-glutamyl-L-norvalyl-L-tryptophyl-L-arginyl-N⁶-(2,4-dinitrophenyl)-L-lysynamide, trifluoroacetate salt

Synonyms: Matrix Metalloproteinase-3 Fluorogenic Substrate, Mca-RPKPVE-Nval-WR-K(Dnp)-NH₂, Stromelysin-1 Fluorogenic Substrate

Peptide Sequence: McaRPKPVENvalWRK(Dnp)-NH₂

MF: C₇₈H₁₁₀N₂₂O₂₀ • XCF₃COOH

FW: 1675.9

Purity: ≥95%

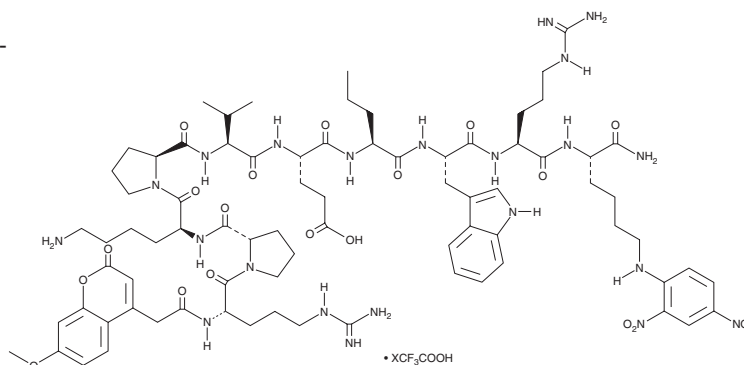
UV/Vis.: λ_{max}: 220 nm

Ex./EM. Max: 328/420 nm

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 fluorogenic substrate (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the MMP-3 fluorogenic substrate (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

MMP-3 fluorogenic substrate is a fluorogenic substrate for matrix metalloproteinase-3 (MMP-3).¹ Upon cleavage by MMP-3, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify MMP-3 activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

Reference

1. Nagase, H., Fields, C.G., and Fields, G.B. Design and characterization of a fluorogenic substrate selectively hydrolyzed by stromelysin 1 (matrix metalloproteinase-3). *J. Biol. Chem.* **269**(33), 20952-20957 (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, 06/15/2023

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