

PRODUCT INFORMATION



Mca-PLA-Nva-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt)

Item No. 24694

Formal Name: 1-[(7-methoxy-2-oxo-2H-1-benzopyran-4-yl) acetyl]-L-prolyl-L-leucyl-L-alanyl-L-norvalyl-3-[[2,4-dinitrophenyl]amino]-L-alanyl-L-alanyl-L-argininamide, trifluoroacetate salt

Synonyms: Matrix Metalloproteinase-2 Fluorogenic Substrate, Mca-Pro-Leu-Ala-Nva-Dap(Dnp)-Ala-Arg-NH₂, MMP-2 Fluorogenic Substrate

MF: C₄₉H₆₈N₁₄O₁₅ • XCF₃COOH

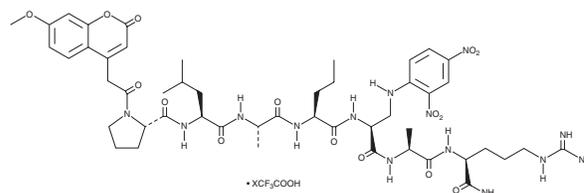
FW: 1,093.1

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

Mca-PLA-Nva-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-PLA-Nva-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Mca-PLA-Nva-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Mca-PLA-Nva-Dap(Dnp)-AR-NH₂ is a fluorogenic substrate for matrix metalloproteinase-2 (MMP-2).¹ Upon cleavage by MMP-2, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify MMP-2 activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

Reference

1. Hu, Y., Lin, X., Wang, P., et al. CRM197 in combination with shRNA interference of VCAM-1 displays enhanced inhibitory effects on human glioblastoma cells. *J. Cell. Physiol.* **230**(8), 1713-1728 (2015).

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