# **PRODUCT** INFORMATION



## Mca-PL (trifluoroacetate salt)

Item No. 24696

Formal Name:	N-[1-[(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]- L-prolyl]-L-leucine, trifluoroacetate salt	~0 ~ ~0 ~ 0
Synonyms:	Matrix Metalloproteinase-2/Matrix Metalloproteinase-7 Fluorogenic Substrate Control, Mca-Pro-Leu-OH, MMP-2/MMP-7 Fluorogenic Substrate Control	
MF:	$C_{23}H_{28}N_2O_7 \bullet XCF_3COOH$	
FW:	444.5	$\forall \uparrow \uparrow \uparrow \uparrow$
Purity:	≥95%	
Supplied as:	A lyophilized powder	
Storage:	-20°C	H O
Stability:	≥4 years	
Information represents the product experifications. Datch experific analytical results are provided on each cartificate of analysis		

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

### Laboratory Procedures

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

### Description

Mca-PL is a fluorogenic peptide that has been used as a building block in the synthesis of Mca-PLGL-Dpa-AR-NH<sub>2</sub> (Item No. 24695), a fluorogenic substrate for matrix metalloproteinase-2 (MMP-2) and MMP-7.1

#### Reference

1. Knight, C.G., Willenbrock, F., and Murphy, G. A novel coumarin-labelled peptide for sensitive continuous assays of the matrix metalloproteinases. FEBS Lett. 296(3), 263-266 (1992).

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.

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, 10/25/2022

## 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