

# PRODUCT INFORMATION



## Mca-PLGL-Dap(Dnp)-AR-NH<sub>2</sub> (trifluoroacetate salt)

Item No. 24149

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

**Synonyms:** Mca-Pro-Leu-Gly-Leu-Dap(Dnp)-Ala-Arg-NH<sub>2</sub>, 7-Methoxycoumarin-4-acetyl-P-L-G-L-β-(2,4-dinitrophenylamino)A-R amide

**MF:** C<sub>49</sub>H<sub>68</sub>N<sub>14</sub>O<sub>15</sub> • XCF<sub>3</sub>COOH

**FW:** 1,093.2

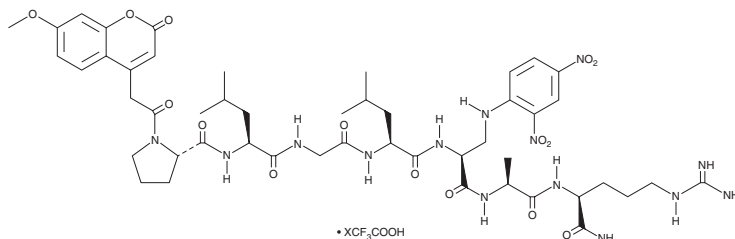
**Purity:** ≥95%

**Ex./Em. Max:** 328/420 nm

**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-PLGL-Dap(Dnp)-AR-NH<sub>2</sub> (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-PLGL-Dap(Dnp)-AR-NH<sub>2</sub> (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Mca-PLGL-Dap(Dnp)-AR-NH<sub>2</sub> (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

### Description

Mca-PLGL-Dap(Dnp)-AR-NH<sub>2</sub> is a fluorogenic peptide substrate for matrix metalloproteinases (MMPs).<sup>1</sup> Upon cleavage by an MMP, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify MMP activity. Mca has excitation and emission maxima of 328 and 420 nm, respectively.

### Reference

1. Edkins, T.J., Alhadeff, J.A., Kwok, V., *et al.* Development and validation of novel enzyme activity methods to assess inhibition of matrix metalloproteinases (MMPs) in human serum by antibodies against enzyme therapeutics. *J. Pharm. Biomed. Anal.* **70**, 408-414 (2012).

#### 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, 01/03/2023

#### 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