

# PRODUCT INFORMATION



## Mca-PLAC(p-OMeBz)-WAR(Dpa)-NH<sub>2</sub> (trifluoroacetate salt)

Item No. 24703

**CAS Registry No.:** 2237216-67-8  
**Formal Name:** 1-[2-(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]-L-prolyl-L-leucyl-L-alanyl-S-[(4-methoxyphenyl)methyl]-L-cysteinyl-L-tryptophyl-L-alanyl-L-arginyl-3-[(2,4-dinitrophenyl)amino]-L-alaninamide, trifluoroacetate salt  
**Synonyms:** Matrix Metalloproteinase-14 Fluorogenic Substrate I, Mca-Pro-Leu-Ala-Cys(p-OMeBz)-Trp-Ala-Arg(Dpa)-NH<sub>2</sub>, MMP-14 Fluorogenic Substrate I

**MF:** C<sub>66</sub>H<sub>82</sub>N<sub>16</sub>O<sub>17</sub>S • CF<sub>3</sub>COOH

**FW:** 1,517.5

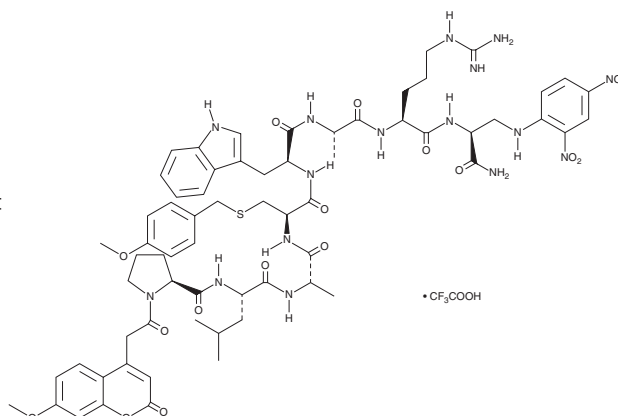
**Purity:** ≥95%

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

### Description

Mca-PLAC(p-OMeBz)-WAR(Dpa)-NH<sub>2</sub> is a fluorogenic substrate for matrix metalloproteinase-14 (MMP-14).<sup>1</sup> Upon cleavage by MMP-14, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify MMP activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

### Reference

1. Eckhouse, S.R., Akerman, A.W., Logdon, C.B., *et al.* Differential membrane type 1 matrix metalloproteinase substrate processing with ischemia-reperfusion: Relationship to interstitial microRNA dynamics and myocardial function. *J. Thorac. Cardiovasc. Surg.* **145**(1), 267-275 (2013).

#### 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, 11/18/2024

#### 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