

# PRODUCT INFORMATION



## Mca-EVKMDAEF-K(Dnp)-NH<sub>2</sub> (ammonium salt)

Item No. 24948

**Formal Name:** 4S-(((S)-1-(((S)-1-amino-6-((2,4-dinitrophenyl)amino)-1-oxohexan-2-yl)amino)-1-oxo-3-phenylpropan-2-yl)carbamoyl)-16S-(4-aminobutyl)-10S-(carboxymethyl)-19S-isopropyl-22S-(2-(7-methoxy-2-oxo-2H-chromen-4-yl)acetamido)-7S-methyl-13S-(2-(methylthio)ethyl)-6,9,12,15,18,21-hexaaxo-5,8,11,14,17,20-hexaazapentacosanedioic acid, ammonium salt

**Synonym:**  $\beta$ -Secretase Fluorogenic Substrate

**MF:** C<sub>66</sub>H<sub>88</sub>N<sub>14</sub>O<sub>23</sub>S • XNH<sub>3</sub>

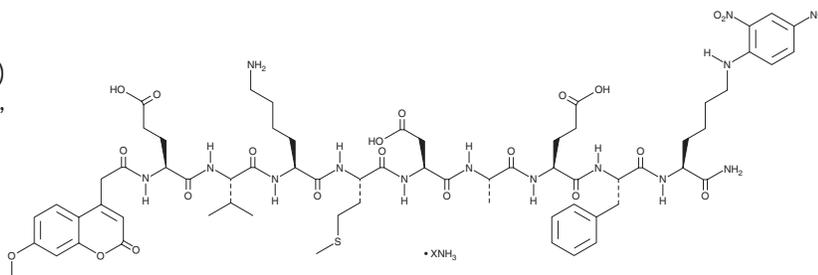
**FW:** 1,477.6

**Purity:**  $\geq$ 95%

**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Mca-EVKMDAEF-K(Dnp)-NH<sub>2</sub> (ammonium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-EVKMDAEF-K(Dnp)-NH<sub>2</sub> (ammonium salt) in a 5% solution of ammonium hydroxide. The solubility of Mca-EVKMDAEF-K(Dnp)-NH<sub>2</sub> (ammonium salt) in 5% ammonium hydroxide is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Mca-EVKMDAEF-K(Dnp)-NH<sub>2</sub> is a fluorogenic substrate for  $\beta$ -secretase.<sup>1</sup> Upon cleavage by  $\beta$ -secretase, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify  $\beta$ -secretase activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

### Reference

1. Gurney, M.E., Bienkowski, M.J., Henrikson, R.L., *et al.* Alzheimer's disease secretase, APP substrates therefor, and uses thereof. US6420534B1. (2002).

#### 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, 12/19/2022

#### 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