

PRODUCT INFORMATION



Mca-SEVNLDAEFR-K(Dnp)-RR-NH₂ (acetate)

Item No. 36879

Formal Name: N-[2-(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]-L-seryl-L- α -glutamyl-L-valyl-L-asparaginyl-L-leucyl-L- α -aspartyl-L-alanyl-L- α -glutamyl-L-phenylalanyl-L-arginyl-N⁶-(2,4-dinitrophenyl)-L-lysyl-L-arginyl-L-argininamide, acetate

MF: C₈₆H₁₂₅N₂₇O₂₉ • XC₂H₄O₂

FW: 2,001.1

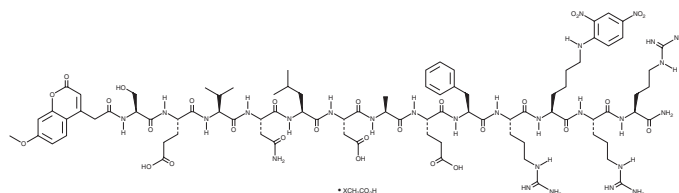
Purity: \geq 95%

Ex./Em. Max: 328/420 nm

Supplied as: A solid

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-SEVNLDAEFR-K(Dnp)-RR-NH₂ (acetate) is supplied as a solid. A stock solution may be made by dissolving the Mca-SEVNLDAEFR-K(Dnp)-RR-NH₂ (acetate) in the solvent of choice, which should be purged with an inert gas. Mca-SEVNLDAEFR-K(Dnp)-RR-NH₂ (acetate) is soluble in DMSO.

Description

Mca-SEVNLDAEFR-K(Dnp)-RR-NH₂ is a fluorogenic substrate for β -secretase (BACE).¹ Upon cleavage by BACE, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify BACE activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

Reference

1. Hong, H.-S., Hwang, J.Y., Son, S.-M., et al. FK506 reduces amyloid plaque burden and induces MMP-9 in A β PP/PS1 double transgenic mice. *J. Alzheimers Dis.* **22(1)**, 97-105 (2010).

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, 08/08/2023

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