

PRODUCT INFORMATION



Mca-HQKLVFFA-K(Dnp)-NH₂ (trifluoroacetate salt)

Item No. 24946

Formal Name: (S)-2-((S)-3-(1H-imidazol-5-yl)-2-(2-(7-methoxy-2-oxo-2H-chromen-4-yl)acetamido)propanamido)-N¹-((5S,8S,11S,14S,17S,20S,23S)-27-amino-11,14-dibenzyl-5-carbamoyl-1-((2,4-dinitrophenyl)amino)-20-isobutyl-17-isopropyl-8-methyl-7,10,13,16,19,22-hexaoxo-6,9,12,15,18,21-hexaazaheptacosan-23-yl)pentanediamide, trifluoroacetate salt

Synonym: α -Secretase Fluorogenic Substrate

MF: C₇₃H₉₅N₁₇O₁₈ • XCF₃COOH

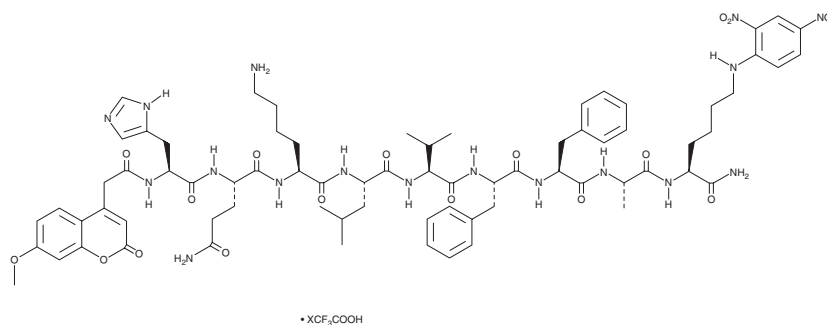
FW: 1,498.6

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-HQKLVFFA-K(Dnp)-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-HQKLVFFA-K(Dnp)-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Mca-HQKLVFFA-K(Dnp)-NH₂ (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

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

Reference

1. Fukumoto, H., Takahashi, H., Tarui, N., *et al.* A noncompetitive BACE1 inhibitor TAK-070 ameliorates A β pathology and behavioral deficits in a mouse model of Alzheimer's disease. *J. Neurosci.* **30(33)**, 11157-11166 (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, 11/18/2022

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