

PRODUCT INFORMATION



Mca-RPPGFSAFK(Dnp)-OH (trifluoroacetate salt)

Item No. 36887

Formal Name: N⁶-(2,4-dinitrophenyl)-N²-(2-(7-methoxy-2-oxo-2H-chromen-4-yl)acetyl)-L-arginyl-L-prolyl-L-prolylglycyl-L-phenylalanyl-L-seryl-L-alanyl-L-phenylalanyl-L-lysine, trifluoroacetate salt

Synonym: (7-methoxycoumarin-4-yl)acetyl-Arg-Pro-Pro-Gly-Phe-Ser-Ala-Phe-Lys(2,4-dinitrophenyl)-OH

MF: C₆₆H₈₁N₁₅O₁₉ • XCF₃COOH

FW: 1,388.4

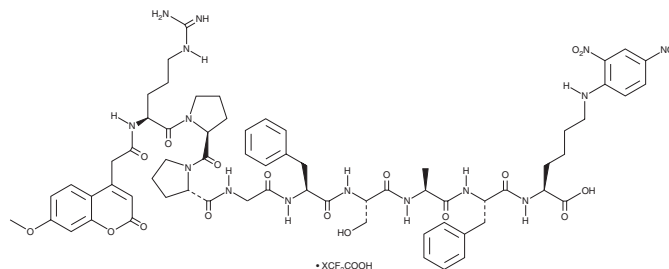
Ex./Em. Max: 328/420 nm

Purity: ≥95%

Supplied as: A solid

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-RPPGFSAFK(Dnp)-OH (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Mca-RPPGFSAFK(Dnp)-OH (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Mca-RPPGFSAFK(Dnp)-OH (trifluoroacetate salt) is soluble in 20% acetonitrile/water.

Description

Mca-RPPGFSAFK(Dnp)-OH is a fluorogenic substrate for endothelin-converting enzyme-1 (ECE-1).¹ Upon cleavage by ECE-1, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify ECE-1 activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

Reference

1. Johnson, G.D. and Ahn, K. Development of an internally quenched fluorescent substrate selective for endothelin-converting enzyme-1. *Anal. Biochem.* **286**(1), 112-118 (2000).

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, 07/05/2023

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