

PRODUCT INFORMATION



Mca-KPLGL-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt)

Item No. 36965

CAS Registry No.: 1392300-20-7

Formal Name: N²-[2-(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]-L-lysyl-L-prolyl-L-leucylglycyl-L-leucyl-3-[(2,4-dinitrophenyl)amino]-L-alanyl-L-alanyl-L-argininamide, monotrifluoroacetate salt

Synonyms:

Mca-Lys-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH₂ (Dpa = N-3-(2,4-dinitrophenyl)-L-α,β-diaminopropionyl), FS-6

MF: C₅₅H₈₀N₁₆O₁₆ • CF₃COOH

FW: 1,335.4

Purity: ≥95%

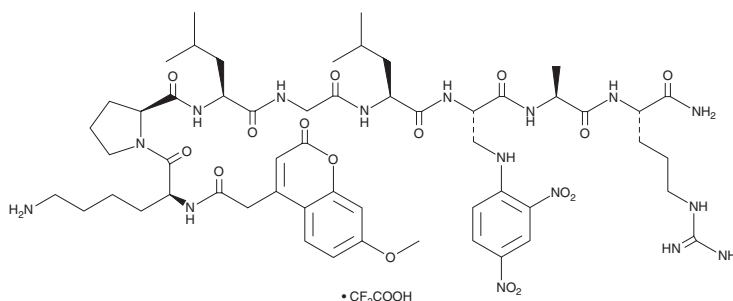
UV/Vis.: λ_{max}: 329 nm

Ex./Em. Max: 328/420 nm

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-KPLGL-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Mca-KPLGL-Dap(Dnp)-AR-NH₂ (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Mca-KPLGL-Dap(Dnp)-AR-NH₂ is a fluorogenic substrate for matrix metalloproteinases (MMPs).¹ Upon cleavage by MMPs, 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. Neumann, U., Kubota, H., Frei, K., *et al.* Characterization of Mca-Lys-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH₂, a fluorogenic substrate with increased specificity constants for collagenases and tumor necrosis factor converting enzyme. *Anal. Biochem.* **328**(2), 166-173 (2004).

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, 10/05/2023

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