

PRODUCT INFORMATION



Mca-YVADAP-K(Dnp)-OH (trifluoroacetate salt)

Item No. 24562

CAS Registry No.: 189696-01-3

Formal Name: N-[2-(7-methoxy-2-oxo-2H-1-benzopyran-4-yl)acetyl]-L-tyrosyl-L-valyl-L-alanyl-L- α -aspartyl-L-alanyl-L-prolyl-N⁶-(2,4-dinitrophenyl)-L-lysine, trifluoroacetate salt

Synonyms: Caspase-1 Fluorogenic Substrate V, Mca-Tyr-Val-Ala-Asp-Ala-Pro-Lys(DNP)-OH, Mca-YVADAP-Lys(Dnp)-OH, (7-Methoxycoumarin-4-yl)acetyl-YVADAPK(2,4-dinitrophenyl)-OH

MF: C₅₃H₆₄N₁₀O₁₉ • XCF₃COOH

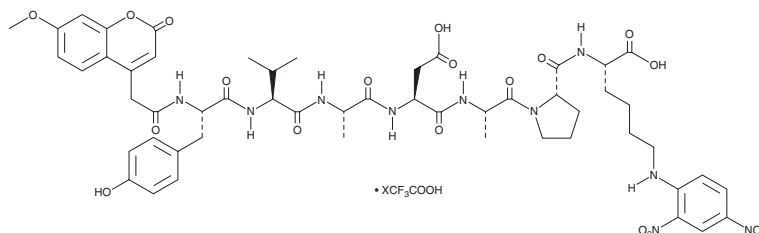
FW: 1,145.1

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-YVADAP-K(Dnp)-OH (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-YVADAP-K(Dnp)-OH (trifluoroacetate salt) in water. The solubility of Mca-YVADAP-K(Dnp)-OH (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Mca-YVADAP-K(Dnp)-OH is a substrate for caspase-1 and angiotensin-converting enzyme 2 (ACE2).^{1,2} Upon cleavage by caspase-1 or ACE2, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify enzyme activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

References

1. Vickers, C., Hales, P., Kaushik, V., *et al.* Hydrolysis of biological peptides by human angiotensin-converting enzyme-related carboxypeptidase. *J. Biol. Chem.* **277**(17), 14838-14843 (2002).
2. Enari, M., Talianian, R.V., Wong, W.W., *et al.* Sequential activation of ICE-like and CPP32-like proteases during Fas-mediated apoptosis. *Nature* **380**(6576), 723-726 (1996).

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, 11/07/2022

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