

# PRODUCT INFORMATION



## Mca-VDQVDGW-K(Dnp)-NH<sub>2</sub> (acetate)

Item No. 24567

**Formal Name:** (3S,6S,9S,12S,18S,21S)-18-((1H-indol-3-yl)methyl)-6-(3-amino-3-oxopropyl)-21-carbamoyl-12-(carboxymethyl)-25-((2,4-dinitrophenyl)amino)-9-isopropyl-3-((S)-2-(2-(7-methoxy-2-oxo-2H-chromen-4-yl)acetamido)-3-methylbutanamido)-4,7,10,13,16,19-hexaaxo-5,8,11,14,17,20-hexaazapentacosanoic acid, acetate

**Synonyms:** Caspase-7 Fluorogenic Substrate I, Mca-VDQVDGW-Lys(Dnp)-NH<sub>2</sub>

**MF:** C<sub>60</sub>H<sub>74</sub>N<sub>14</sub>O<sub>21</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

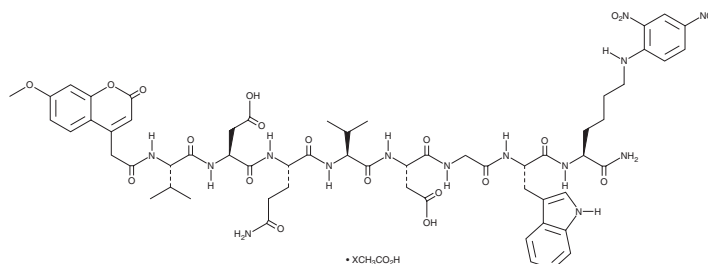
**FW:** 1,327.3

**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-VDQVDGW-K(Dnp)-NH<sub>2</sub> (acetate) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Mca-VDQVDGW-K(Dnp)-NH<sub>2</sub> (acetate) in water. The solubility of Mca-VDQVDGW-K(Dnp)-NH<sub>2</sub> (acetate) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Mca-VDQVDGW-K(Dnp)-NH<sub>2</sub> is a substrate for caspase-7.<sup>1</sup> Upon cleavage by caspase-7, 7-methoxycoumarin-4-acetyl (Mca) is released and its fluorescence can be used to quantify caspase-7 activity. Mca displays excitation/emission maxima of 328/420 nm, respectively.

### Reference

1. Yu, M.-S., Suen, K.-C., Kwok, N.-S., *et al.* Beta-amyloid peptides induces neuronal apoptosis via a mechanism independent of unfolded protein responses. *Apoptosis* **11**(5), 687-700 (2006).

#### 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, 05/30/2024

#### 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