

# PRODUCT INFORMATION



## Z-DEVD-AFC

Item No. 24564

**CAS Registry No.:** 1135416-11-3  
**Formal Name:** (5S,8S,11S,14S)-8-(2-carboxyethyl)-5-(carboxymethyl)-11-isopropyl-3,6,9,12-tetraoxo-14-((2-oxo-4-(trifluoromethyl)-2H-chromen-7-yl)carbamoyl)-1-phenyl-2-oxa-4,7,10,13-tetraazahexadecan-16-oic acid  
**Synonym:** Caspase-3 Fluorogenic Substrate IV

**MF:** C<sub>36</sub>H<sub>38</sub>F<sub>3</sub>N<sub>5</sub>O<sub>14</sub>

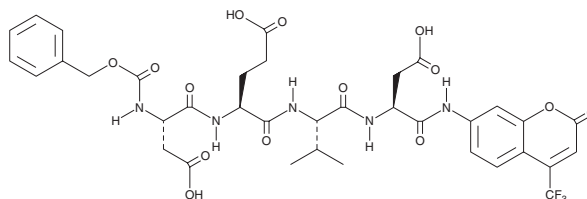
**FW:** 821.7

**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

Z-DEVD-AFC is supplied as a lyophilized powder. A stock solution may be made by dissolving the Z-DEVD-AFC in the solvent of choice. Z-DEVD-AFC is soluble in the organic solvent formic acid, which should be purged with an inert gas, at a concentration of approximately 1 mg/ml.

### Description

Z-DEVD-AFC is a fluorogenic substrate for caspase-3.<sup>1</sup> Upon cleavage by caspase-3, 7-amino-7-trifluoromethylcoumarin (AFC) is released and its fluorescence can be used to quantify caspase-3 activity. AFC displays excitation/emission maxima of 400/505 nm, respectively.

### Reference

1. Inayat-Hussain, S.H. Processing/activation of CPP32-like proteases is involved in transforming growth factor  $\beta_1$ -induced apoptosis in rat hepatocytes. *Hepatology* **25**(6), 1516-1526 (1997).

#### 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, 12/12/2022

#### 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](http://WWW.CAYMANCHEM.COM)