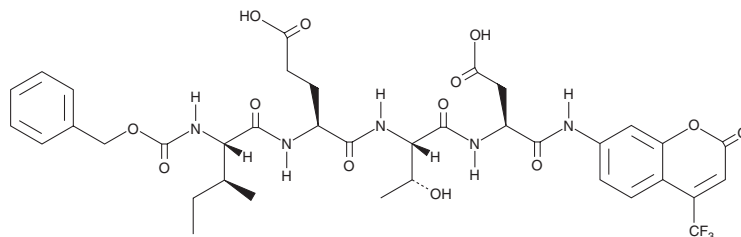


PRODUCT INFORMATION



Z-IETD-AFC Item No. 27101

CAS Registry No.: 219138-02-0
Formal Name: N-[(phenylmethoxy)carbonyl]-L-isoleucyl-L- α -glutamyl-L-threonyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L- α -asparagine
Synonym: Z-Ile-Glu-Thr-Asp-7-amino-4-Trifluoromethylcoumarin
MF: C₃₇H₄₂F₃N₅O₁₃
FW: 821.8
Purity: \geq 90%
Ex./Em. Max: 400/505 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Z-IETD-AFC is supplied as a solid. A stock solution may be made by dissolving the Z-IETD-AFC in the solvent of choice, which should be purged with an inert gas. Z-IETD-AFC is soluble in the organic solvent DMSO.

Description

Z-IETD-AFC is a fluorogenic substrate for caspase-8 and granzyme B.¹ Upon enzymatic cleavage by caspase-8 or granzyme B, 7-amino-4-trifluoromethylcoumarin (AFC) is released and its fluorescence can be used to quantify caspase-8 or granzyme B activity. AFC displays excitation/emission maxima of 400/505 nm, respectively.

Reference

1. Nie, Z., Phenix, B.N., Lum, J.J., *et al.* HIV-1 protease processes procaspase 8 to cause mitochondrial release of cytochrome c, caspase cleavage and nuclear fragmentation. *Cell Death Differ.* **9(11)**, 1172-1184 (2002).

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, 10/27/2022

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