

PRODUCT INFORMATION



Z-VEID-AFC (trifluoroacetate salt)

Item No. 27146

Formal Name: N-[(phenylmethoxy)carbonyl]-L-valyl-L- α -glutamyl-L-isoleucyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L- α -asparagine, trifluoroacetate salt

Synonyms: Z-Val-Glu-Ile-Asp-AFC, Z-Val-Glu-Ile-Asp-7-amino-4-trifluoromethylcoumarin

MF: C₃₈H₄₄F₃N₅O₁₂ • XCF₃COOH

FW: 819.8

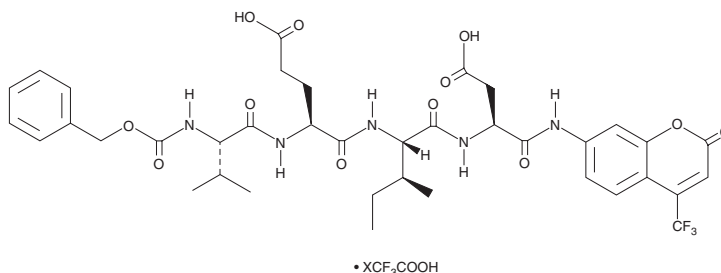
Purity: ≥95%

Ex./Em. Max: 400/505 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

Z-VEID-AFC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Z-VEID-AFC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Z-VEID-AFC (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Z-VEID-AFC is a fluorogenic substrate for caspase-6.¹ Upon enzymatic cleavage by caspase-6, 7-amino-4-trifluoromethylcoumarin (AFC) is released and its fluorescence can be used to quantify caspase-6 activity. AFC displays excitation/emission maxima of 400/505 nm, respectively.

Reference

1. Li, L., Prevette, D., Oppenheim, R.W., *et al.* Involvement of specific caspases in motoneuron cell death *in vivo* and *in vitro* following trophic factor deprivation. *Mol. Cell Neurosci.* **12(3)**, 157-167 (1998).

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, 11/18/2022

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