

PRODUCT INFORMATION



(Z-DEVD)₂-Rh 110 (trifluoroacetate salt)

Item No. 27432

Formal Name: 4,4'-[9-(2-carboxyphenyl)xanthylum-3,6-diyl]bis[N-[(phenylmethoxy)carbonyl]-L-α-aspartyl-L-α-glutamyl-L-valyl-L-α-asparagine, inner salt, trifluoroacetate salt

Synonym: (Z-Asp-Glu-Val-Asp)₂-Rhodamine 110

MF: C₇₂H₇₈N₁₀O₂₇ • XCF₃COOH

FW: 1,515.4

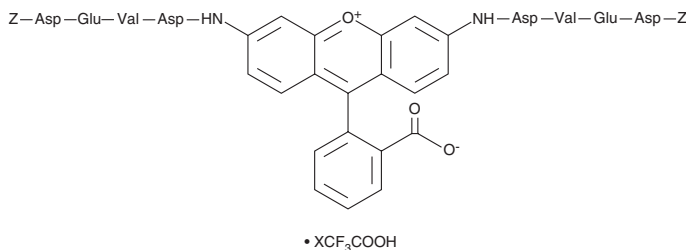
Purity: ≥95%

Ex./Em. Max: 496/520 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-DEVD)₂-Rh 110 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the (Z-DEVD)₂-Rh 110 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. (Z-DEVD)₂-Rh 110 (trifluoroacetate salt) is soluble in the formic acid at a concentration of approximately 1 mg/ml.

Description

(Z-DEVD)₂-Rh 110 is a fluorogenic substrate for caspase-3.¹ Upon enzymatic cleavage by caspase-3, rhodamine 110 (Item No. 19061) is released and its fluorescence can be used to quantify caspase-3 activity. Rhodamine 110 displays excitation/emission maxima of 496/520 nm, respectively.

Reference

1. Liu, F., Wang, J., Yang, X., *et al.* Discovery of a highly selective STK16 kinase inhibitor. *ACS Chem. Biol.* **11**(6), 1537-1543 (2016).

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, 10/06/2022

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