

PRODUCT INFORMATION



Dabcyl-YVADAPV-EDANS

Item No. 24993

CAS Registry No.: 161877-70-9
Formal Name: N-[4-[[4-(dimethylamino)phenyl]azo]benzoyl]-L-tyrosyl-L-valyl-L-alanyl-L- α -aspartyl-L-alanyl-L-prolyl-N-[2-[[5-sulfo-1-naphthalenyl]amino]ethyl]-L-valinamide

Synonym: Dabcyl-ICE-EDANS

MF: C₆₁H₇₆N₁₂O₁₄S

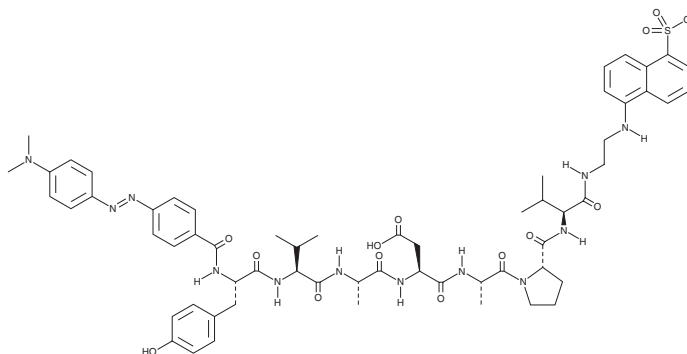
FW: 1,233.4

Purity: \geq 95%

Supplied as: A lyophilized powder

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

Dabcyl-YVADAPV-EDANS is supplied as a lyophilized powder. Aqueous solutions of Dabcyl-YVADAPV-EDANS can be prepared by directly dissolving the lyophilized powder in aqueous buffers. Dabcyl-YVADAPV-EDANS is soluble in a 5% aqueous ammonia solution at a concentration of approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dabcyl-YVADAPV-EDANS is a fluorogenic substrate for caspase-1.¹ Upon enzymatic cleavage by caspase-1, EDANS is separated from the Dabcyl quencher and can be used to quantify caspase-1 activity. EDANS displays excitation/emission maxima of 340/485 nm, respectively.

Reference

1. Pennington, M.W. and Thornberry, N.A. Synthesis of a fluorogenic interleukin-1 beta converting enzyme substrate based on resonance energy transfer. *Pept. Res.* **7(2)**, 72-76 (1994).

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