

PRODUCT INFORMATION



Dabcyl-RGVVNASSRLA-EDANS (trifluoroacetate salt)

Item No. 24994

Formal Name: N²-[4-[2-[4-(dimethylamino)phenyl] diazenyl]benzoyl]-L-arginylglycyl-L-valyl-L-valyl-L-asparaginyll-L-alanyl-L-seryl-L-seryl-L-arginyl-L-leucyl-N-[2-[(5-sulfo-1-naphthalenyl)amino]ethyl]-L-alaninamide, trifluoroacetate salt
Synonym: Fluorogenic Human CMV Protease Substrate

MF: C₇₃H₁₀₉N₂₃O₁₈S • XCF₃COOH

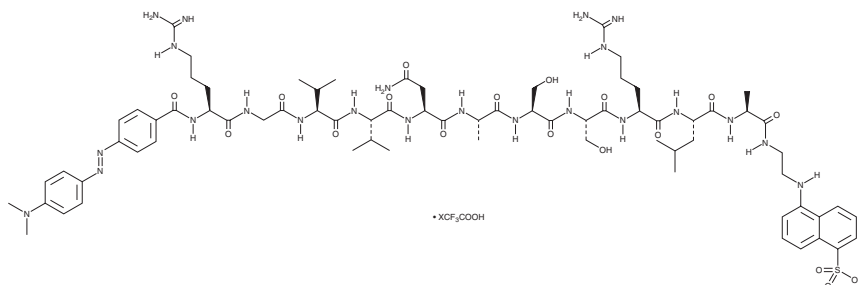
FW: 1,628.9

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

Dabcyl-RGVVNASSRLA-EDANS (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Dabcyl-RGVVNASSRLA-EDANS (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dabcyl-RGVVNASSRLA-EDANS (trifluoroacetate salt) is soluble in the organic solvent formic acid. The solubility of Dabcyl-RGVVNASSRLA-EDANS (trifluoroacetate salt) in formic acid is approximately 1 mg/ml.

Description

Dabcyl-RGVVNASSRLA-EDANS is a fluorogenic substrate for human cytomegalovirus (HCMV) protease.¹ Upon enzymatic cleavage by HCMV protease at the Ala-Ser site, EDANS is separated from the Dabcyl quencher and its fluorescence can be used to quantify HCMV protease activity. EDANS displays excitation/emission maxima of 340/485 nm, respectively.

Reference

1. Holskin, B.P., Bukhtiyarova, M., Dunn, B.M., *et al.* A continuous fluorescence-based assay of human cytomegalovirus protease using a peptide substrate. *Anal. Biochem.* **227**(1), 148-155 (1995).

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