

PRODUCT INFORMATION



DabcyI-GABA-RPKPVE-Nva-WR-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt)
Item No. 24843

Formal Name: N²-[4-[4-[2-[4-(dimethylamino)phenyl]diazanyl]benzoyl]amino]-1-oxobutyl]-L-arginyl-L-prolyl-L-lysyl-L-prolyl-L-valyl-L-α-glutamyl-L-norvalyl-L-tryptophyl-L-arginyl-N-[2-[(5-sulfo-1-naphthalenyl)amino]ethyl]-L-glutamyl-L-alanyl-L-lysineamide, trifluoroacetate salt

MF: C₉₉H₁₄₄N₂₈O₂₀S • XCF₃COOH

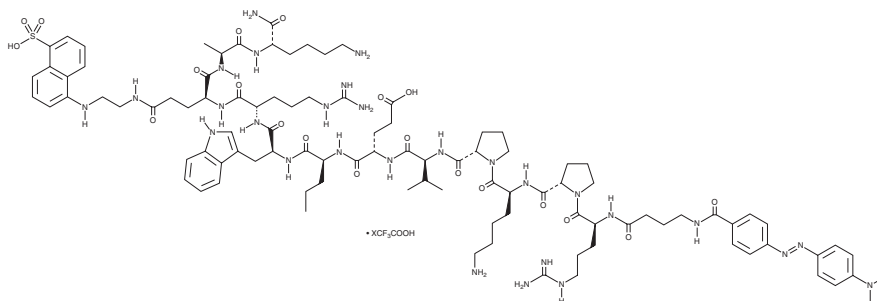
FW: 2,078.5

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

DabcyI-GABA-RPKPVE-Nva-WR-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the DabcyI-GABA-RPKPVE-Nva-WR-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) in water. The solubility of DabcyI-GABA-RPKPVE-Nva-WR-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DabcyI-GABA-RPKPVE-Nva-WR-Glu(EDANS)-AK-NH₂ is a fluorogenic substrate for matrix metalloproteinase-3 (MMP-3).¹ Upon enzymatic cleavage by MMP-3 at Glu-Nva, EDANS is separated from the DabcyI quencher and its fluorescence can be used to quantify MMP-3 activity. EDANS displays excitation/emission maxima of 340/485 nm, respectively.

Reference

1. Beekman, B., Drijfhout, J.W., Bloemhoff, W., *et al.* Convenient fluorometric assay for matrix metalloproteinase activity and its application in biological media. *FEBS Lett.* **390(2)**, 221-225 (1996).

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, 04/09/2018

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