

PRODUCT INFORMATION



Dabcyl-GABA-PQGL-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt)

Item No. 24842

Formal Name: 1-[4-[[4-[2-[4-(dimethylamino)phenyl]diazanyl]benzoyl]amino]-1-oxobutyl]-L-prolyl-L-glutaminyglycyl-L-leucyl-N-[2-[(5-sulfo-1-naphthalenyl)amino]ethyl]-L-glutaminy-L-alanyl-L-lysineamide, trifluoroacetate salt

Synonym: MMP Fluorogenic Substrate III

MF: C₆₃H₈₈N₁₆O₁₄S

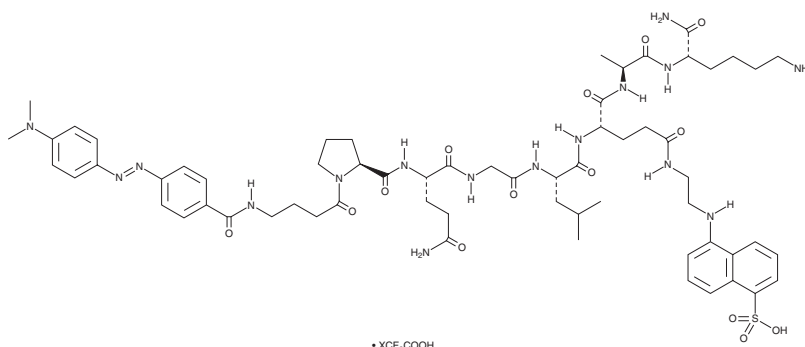
FW: 1,325.6

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-GABA-PQGL-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Dabcyl-GABA-PQGL-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dabcyl-GABA-PQGL-Glu(EDANS)-AK-NH₂ (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Dabcyl-GABA-PQGL-Glu(EDANS)-AK-NH₂ is a fluorogenic substrate for matrix metalloproteinase-1 (MMP-1), MMP-2, MMP-3, and MMP-9.¹ Upon enzymatic cleavage by MMP-1, MMP-2, MMP-3, or MMP-9 at the Gly-Leu site, EDANS is separated from the Dabcyl quencher and its fluorescence can be used to quantify MMP-1, MMP-2, MMP-3, and MMP-9 activity.^{1,2} EDANS displays excitation/emission maxima of 340/485 nm, respectively.

References

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).
2. Krismastuti, F.S.H., Pace, S., and Voelcker, N.H. Porous silicon resonant microcavity biosensor for matrix metalloproteinase detection. *Adv. Funct. Mater.* **24(23)**, 3639-3650 (2014).

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