

PRODUCT INFORMATION



FITC-C6-PLALWAR-Lys(Biotin)-NH₂ (trifluoroacetate salt)

Item No. 24844

Formal Name: 5-(3-(6-((S)-2-(((4S,7S,10S,13S,16S,19S,22S)-13-((1H-indol-3-yl)methyl)-22-carbamoyl-19-(3-guanidinopropyl)-10-isobutyl-2,7,16-trimethyl-5,8,11,14,17,20,28-heptaoxo-32-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)-6,9,12,15,18,21,27-heptaazadotriacontan-4-yl)carbamoyl)pyrrolidin-1-yl)-6-oxohexyl)thioureido)-2-(6-hydroxy-3-oxo-3H-xanthen-9-yl)benzoic acid, trifluoroacetate salt

Synonyms: FITC-C6-Pro-Leu-Ala-Leu-Trp-Ala-Arg-Lys(Biotin)-NH₂, MMP-1/MMP-2/MMP-3 Fluorogenic Substrate

MF: C₈₃H₁₁₂N₁₈O₁₆S₂ • XCF₃COOH

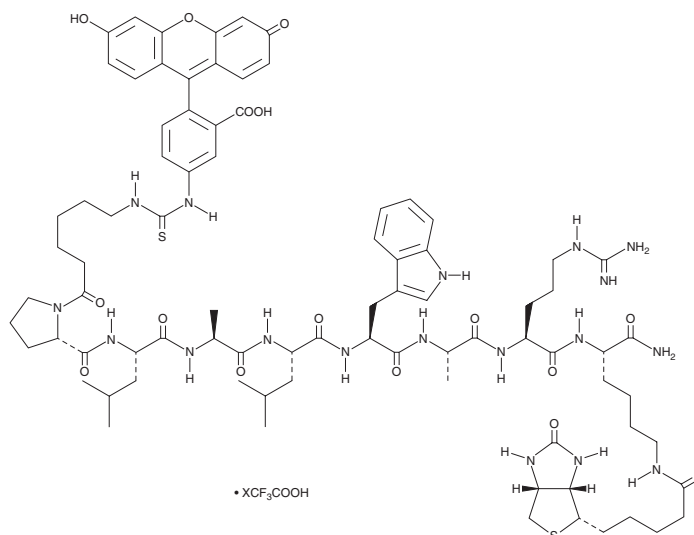
FW: 1,682.0

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

FITC-C6-PLALWAR-Lys(Biotin)-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the FITC-C6-PLALWAR-Lys(biotin)-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. FITC-C6-PLALWAR-Lys(Biotin)-NH₂ (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

FITC-C6-PLALWAR-Lys(Biotin)-NH₂ is a substrate for matrix metalloproteinase-1 (MMP-1), MMP-2, and MMP-3 that can be sequestered by streptavidin via its biotin group.¹ Upon cleavage by MMP-1, MMP-2, or MMP-3, FITC is released and its fluorescence can be used to quantify MMP-1, MMP-2, and MMP-3 activity. FITC displays excitation/emission maxima of 485/535 nm, respectively.

Reference

1. Jacobsen, E.J., Mitchell, M.A., Hendges, S.K., *et al.* Synthesis of a series of stromelysin-selective thiadiazole urea matrix metalloproteinase inhibitors. *J. Med. Chem.* **42(9)**, 1525-1536 (1999).

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, 12/19/2022

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