

PRODUCT INFORMATION



FITC-C6-YVADAP-K(Dnp) (trifluoroacetate salt)

Item No. 27422

Formal Name: 5-(3-((2S,5S,8S,11S,14S)-1-((S)-2-(((S)-1-carboxy-5-((2,4-dinitrophenyl)amino)pentyl)carbamoyl)pyrrolidin-1-yl)-5-(carboxymethyl)-14-(4-hydroxybenzyl)-11-isopropyl-2,8-dimethyl-1,4,7,10,13,16-hexaoxo-3,6,9,12,15-pentaazahenicosan-21-yl)thioureido)-2-(6-hydroxy-3-oxo-3H-xanthen-9-yl)benzoic acid, trifluoroacetate salt

Synonym: FITC-C6-Tyr-Val-Ala-Asp-Ala-Pro-Lys(Dnp)-OH

MF: C₆₈H₇₈N₁₂O₂₁S • XCF₃COOH

FW: 1,431.5

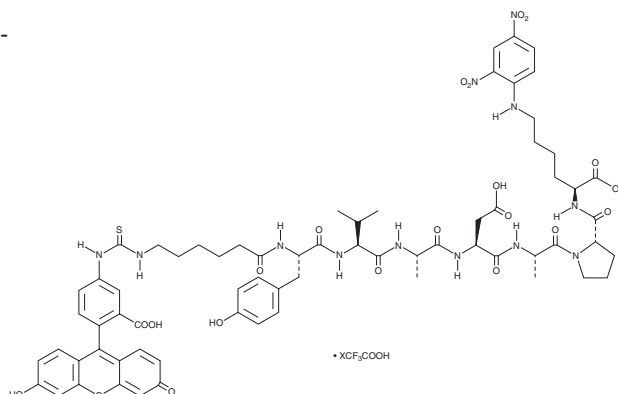
Purity: ≥95%

Ex./Em. Max: 485/535 nm

Supplied as: A solid

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-YVADAP-K(Dnp) (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of FITC-C6-YVADAP-K(Dnp) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. FITC-C6-YVADAP-K(Dnp) (trifluoroacetate salt) 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

FITC-C6-YVADAP-K(Dnp) is a caspase-1 fluorogenic substrate.¹ Upon enzymatic cleavage by caspase-1, FITC is released and its fluorescence can be used to quantify enzyme activity. FITC displays excitation/emission maxima of 485/535 nm, respectively.

References

1. Vickers, C., Hales, P., Kaushik, V., *et al.* Hydrolysis of biological peptides by human angiotensin-converting enzyme-related carboxypeptidase. *J. Biol. Chem.* **277**(17), 14838-14843 (2002).

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, 11/30/2022

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