

PRODUCT INFORMATION



Chymotrypsin Substrate II, Fluorogenic

Item No. 41593

CAS Registry No.: 88467-45-2

Formal Name: N-(3-carboxy-1-oxopropyl)-L-alanyl-L-alanyl-L-prolyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-phenylalaninamide

Synonyms: Suc-AAPF-AMC,
Suc-Ala-Ala-Pro-Phe-AMC,
Suc-Ala₂-Pro-Phe-AMC,
Succinyl-Ala-Ala-Pro-Phe-AMC

MF: C₃₄H₃₉N₅O₉

FW: 661.7

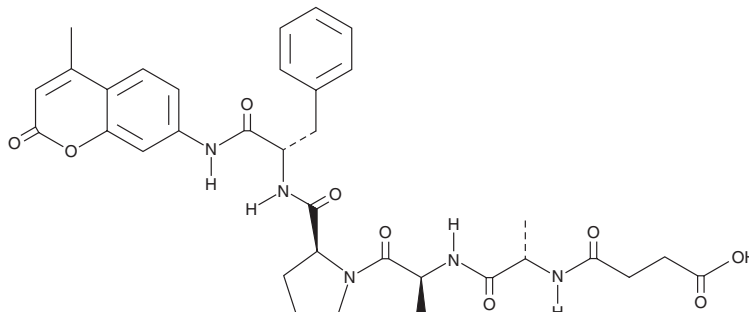
Purity: ≥98%

Ex./Em. Max: 340-360/440-460 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

Chymotrypsin substrate II, fluorogenic is supplied as a solid. A stock solution may be made by dissolving the chymotrypsin substrate II, fluorogenic in the solvent of choice, which should be purged with an inert gas. Chymotrypsin substrate II, fluorogenic is soluble (≥10 mg/ml) in DMSO and ethanol.

Description

Chymotrypsin substrate II, fluorogenic is a fluorogenic substrate for chymotrypsin and chymotrypsin-like proteases.^{1,2} Upon enzymatic cleavage by proteases, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify protease activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

References

1. Oshima, G. Interaction of α-chymotrypsin with dimyristoyl phosphatidylcholine vesicles. *J. Biochem.* **95**(4), 1131-1136. (1984).
2. Irvine, G.B., Ennis, M., and Williams, C.H. Visual detection of peptidase activity using fluorogenic substrates in a microtiter plate assay. *Anal. Biochem.* **185**(2), 304-307 (1990).

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, 09/09/2024

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