

PRODUCT INFORMATION



Suc-AAP-Abu-pNA (trifluoroacetate salt)

Item No. 24704

Formal Name: N-(3-carboxy-1-oxopropyl)-L-alanyl-L-alanyl-L-prolyl-N-(4-nitrophenyl)-L-2-amino-butanamide, trifluoroacetate salt

Synonym: Elastase Colorimetric Substrate IV

MF: $C_{25}H_{34}N_6O_9 \cdot XCF_3COOH$

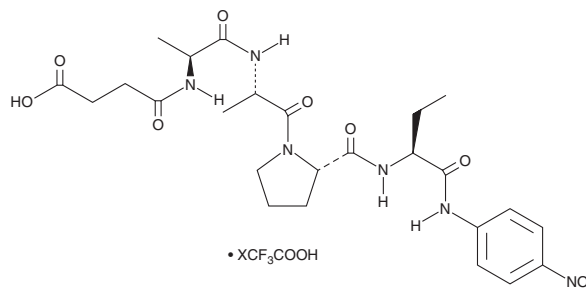
FW: 562.6

Purity: $\geq 95\%$

Supplied as: A lyophilized powder

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Suc-AAP-Abu-pNA (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Suc-AAP-Abu-pNA (trifluoroacetate salt) in water. The solubility of Suc-AAP-Abu-pNA (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Suc-AAP-Abu-pNA is a chromogenic peptide substrate for pancreatic elastase.^{1,2} It is cleaved by elastase to release *p*-nitroaldehyde (*p*-NA), which can be quantified by colorimetric detection at 405 nm as a measure of enzyme activity.

References

1. Del Mar, E.G., Largman, C., Brodrick, J.W., et al. Substrate specificity of human pancreatic elastase 2. *Biochemistry* **19**(3), 468-472 (1980).
2. Largman, C. Isolation and characterization of rat pancreatic elastase. *Biochemistry* **22**(16), 3763-3770 (1983).

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, 12/11/2023

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