

PRODUCT INFORMATION



Chymotrypsin Substrate I, Colorimetric (trifluoroacetate salt)

Item No. 36194

Formal Name: N-(3-carboxy-1-oxopropyl) glycylglycyl-N-(4-nitrophenyl)-L-phenylalaninamide, trifluoroacetate salt

Synonyms: Suc-GGF-pNA, Suc-Gly-Gly-Phe-pNA, Suc-Gly-Gly-Phe-p-Nitroanilide

MF: C₂₃H₂₅N₅O₈ • XCF₃COOH

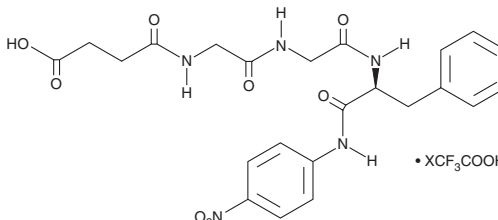
FW: 499.5

Purity: ≥95%

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 I, colorimetric (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the chymotrypsin substrate I, colorimetric (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Chymotrypsin substrate I, colorimetric (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of chymotrypsin substrate I, colorimetric (trifluoroacetate salt) in these solvents is approximately 10 and 5 mg/ml, respectively.

Description

Chymotrypsin substrate I, colorimetric is a colorimetric substrate for chymotrypsin.¹ Chymotrypsin binds and hydrolyzes chymotrypsin substrate I, colorimetric to release *p*-nitroanilide (*p*NA), which can be quantified by colorimetric detection at 405 nm as a measure of chymotrypsin activity.

Reference

1. Kasafírek, E. and Bartík, M. N-Acylated phenylalanine *p*-nitroanilides: New substrates for chymotrypsin. *Collect. Czech. Chem. Commun.* **45**, 442-451 (1980).

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