

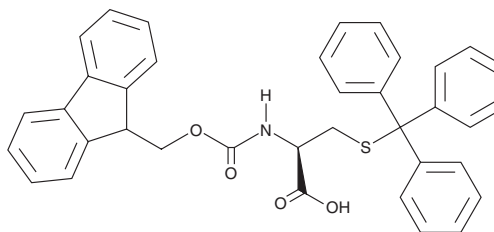
PRODUCT INFORMATION



Fmoc-Cys(Trt)-OH

Item No. 32828

CAS Registry No.: 103213-32-7
Formal Name: N-[(9H-fluoren-9-ylmethoxy)carbonyl]-S-(triphenylmethyl)-L-cysteine
MF: C₃₇H₃₁NO₄S
FW: 585.7
Purity: ≥98%
Supplied as: A crystalline 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

Fmoc-Cys(Trt)-OH is supplied as a crystalline solid. A stock solution may be made by dissolving the Fmoc-Cys(Trt)-OH in the solvent of choice, which should be purged with an inert gas. Fmoc-Cys(Trt)-OH is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Fmoc-Cys(Trt)-OH in these solvents is approximately 30 mg/ml.

Description

Fmoc-Cys(Trt)-OH is a building block.^{1,2} It has been used in the solid-phase synthesis of peptide thioesters.

References

1. Ingenito, R., Bianchi, E., Fattori, D., *et al.* Solid phase synthesis of peptide C-terminal thioesters by Fmoc/t-Bu chemistry. *J. Am. Chem. Soc.* **121(49)**, 11369-11374 (1999).
2. Blanco-Canosa, J.B., Nardone, B., Alberico, F., *et al.* Chemical protein synthesis using a second-generation N-acylurea linker for the preparation of peptide-thioester precursors. *J. Am. Chem. Soc.* **137(22)**, 7197-7209 (2015).

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, 11/10/2022

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