

PRODUCT INFORMATION



Fmoc-L-Cys(Mmt)-OH

Item No. 42177

CAS Registry No.: 177582-21-7

Formal Name: N-[(9H-fluoren-9-ylmethoxy) carbonyl]-S-[(4-methoxyphenyl) diphenylmethyl]-L-cysteine

MF: $C_{38}H_{33}NO_5S$

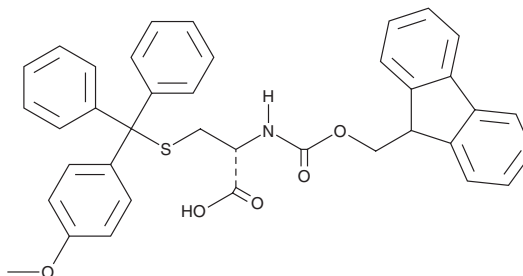
FW: 615.7

Purity: $\geq 98\%$

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

Fmoc-L-Cys(Mmt)-OH is supplied as a solid. A stock solution may be made by dissolving the Fmoc-L-Cys(Mmt)-OH in the solvent of choice, which should be purged with an inert gas. Fmoc-L-Cys(Mmt)-OH is sparingly soluble (1-10 mg/ml) in DMSO and ethanol.

Description

Fmoc-L-Cys(Mmt)-OH is an amino acid-containing building block.^{1,2} It has been used in the solid-phase synthesis of S-palmitoylated peptide antigens for use in macrophage uptake assays and stapled peptide dual agonists of the glucagon-like peptide 1 receptor (GLP-1R) and cholecystokinin 2 (CCK₂) receptor.

References

1. Nadège, P., Erwann, G., Judith, M.G., *et al.* Solid-phase synthesis of a biotin-labelled thiopalmitoylated myelin proteolipid protein epitope and application in the study of uptake of antigen by macrophages. *Int. J. Pept. Res. Ther.* **10**, 581-588 (2003).
2. Chen, X., Fu, J., Zhou, F., *et al.* Stapled and *Xenopus* glucagon-like peptide 1 (GLP-1)-based dual GLP-1/gastrin receptor agonists with improved metabolic benefits in rodent models of obesity and diabetes. *J. Med. Chem.* **63**(21), 12595-12613 (2020).

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, 11/14/2024

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