

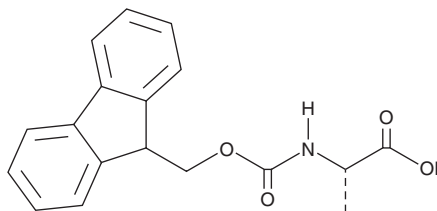
PRODUCT INFORMATION



Fmoc-L-Alanine

Item No. 36457

CAS Registry No.: 35661-39-3
Formal Name: N-[(9H-fluoren-9-ylmethoxy) carbonyl]-L-alanine
Synonyms: Fmoc-Alanine, NPC 14688, NSC 334296
MF: C₁₈H₁₇NO₄
FW: 311.3
Purity: ≥98%
UV/Vis.: λ_{max}: 266 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

Fmoc-L-alanine is supplied as a solid. A stock solution may be made by dissolving the Fmoc-L-alanine in the solvent of choice, which should be purged with an inert gas. Fmoc-L-alanine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Fmoc-L-alanine in these solvents is approximately 10 mg/ml.

Description

Fmoc-L-alanine is a building block.^{1,2} It has been used in the synthesis of antimicrobial and fluorescent peptides.

References

1. Wang, Y., Zhu, G., Wang, W., *et al.* Rational design of HJH antimicrobial peptides to improve antimicrobial activity. *Bioorg. Med. Chem. Lett.* **83**, 129176 (2023).
2. Qin, S., Liu, Q., Li, K., *et al.* Neuropilin 1-targeted near-infrared fluorescence probes for tumor diagnosis. *Bioorg. Med. Chem. Lett.* **84**, 129196 (2023).

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, 05/15/2023

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