

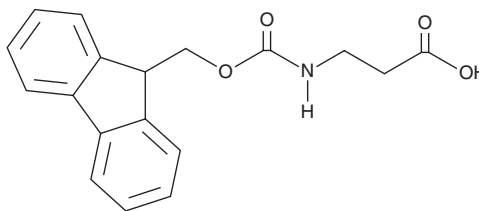
PRODUCT INFORMATION



Fmoc-β-Alanine

Item No. 41969

CAS Registry No.: 35737-10-1
Formal Name: N-[(9H-fluoren-9-ylmethoxy) carbonyl]-β-alanine
MF: C₁₈H₁₇NO₄
FW: 311.3
Purity: ≥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-β-alanine is supplied as a solid. A stock solution may be made by dissolving the Fmoc-β-alanine in the solvent of choice, which should be purged with an inert gas. Fmoc-β-alanine is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Fmoc-β-alanine is a building block.¹ It has been used in the solid-phase synthesis of glucagon receptor peptide antagonists and bradykinin B₁ and B₂ receptor pseudopeptide antagonists.^{1,2}

References

1. Lau, J., Behrens, C., Sidelmann, U.G., *et al.* New β-alanine derivatives are orally available glucagon receptor antagonists. *J. Med. Chem.* **50(1)**, 113-128 (2007).
2. Galoppini, C., Meini, S., Tancredi, M., *et al.* A new class of pseudopeptide antagonists of the kinin B₁ receptor containing alkyl spacers. *J. Med. Chem.* **42(3)**, 409-414 (1999).

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, 10/28/2024

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