

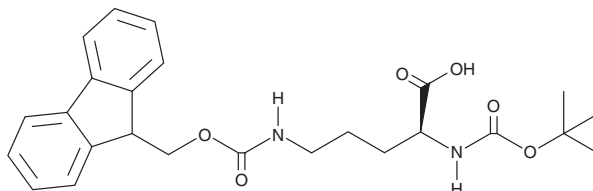
PRODUCT INFORMATION



Boc-Orn(Fmoc)-OH

Item No. 30495

CAS Registry No.: 150828-96-9
Formal Name: N²-[(1,1-dimethylethoxy)carbonyl]-N⁵-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-ornithine
MF: C₂₅H₃₀N₂O₆
FW: 454.5
Purity: ≥98%
UV/Vis.: λ_{max}: 265 nm
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

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

Boc-Orn(Fmoc)-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Boc-Orn(Fmoc)-OH should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Boc-Orn(Fmoc)-OH has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Boc-Orn(Fmoc)-OH is an amino acid-containing building block.¹ It has been used to introduce δ-linked ornithine turns into synthetic macrocyclic β-sheets.^{1,2}

References

1. Spencer, R., Chen, K.H., Manuel, G., *et al.* Recipe for β-sheets: Foldamers containing amyloidogenic peptide sequences. *Eur. J. Org. Chem.* **2013**(17), 3523–3528 (2013).
2. Nowick, J.S., Brower, J.O., Khakshoor, O., *et al.* Cyclic Modular β-Sheets. *Understanding Biology Using Peptides*. Sylvie E. Blondelle, editor, American Peptide Society, Irvine, CA (2005).

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, 11/23/2022

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