

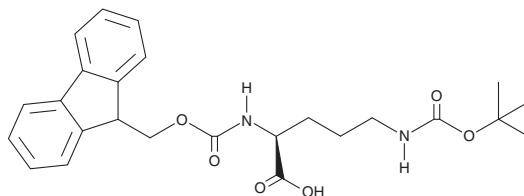
PRODUCT INFORMATION



Fmoc-Orn(Boc)-OH

Item No. 30471

CAS Registry No.: 109425-55-0
Formal Name: N⁵-[(1,1-dimethylethoxy) carbonyl]-N²-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-ornithine
MF: C₂₅H₃₀N₂O₆
FW: 454.5
Purity: ≥98%
UV/Vis.: λ_{max}: 266 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

Fmoc-Orn(Boc)-OH is supplied as a crystalline solid. A stock solution may be made by dissolving the Fmoc-Orn(Boc)-OH in the solvent of choice, which should be purged with an inert gas. Fmoc-Orn(Boc)-OH is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Fmoc-Orn(Boc)-OH in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

Fmoc-Orn(Boc)-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Fmoc-Orn(Boc)-OH should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Fmoc-Orn(Boc)-OH has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fmoc-Orn(Boc)-OH is an ornithine-containing amino acid building block¹ It has been used to conjugate ornithine to GFP-labeled peptides, enhancing cell permeability when compared to unconjugated GFP-labeled peptides.²

References

1. Wender, P.A., Rothbard, J.B., Jessop, T.C., *et al.* Oligocarbamate molecular transporters: Design, synthesis, and biological evaluation of a new class of transporters for drug delivery. *J. Am. Chem. Soc.* **124**(45), 13382-13383 (2002).
2. Kumar, J., Yellepeddi, V.K., Davies, G.E., *et al.* Enhanced gene transfection efficiency by polyamidoamine (PAMAM) dendrimers modified with ornithine residues. *Int. J. Pharm.* **392**(1-2), 294-303 (2010).

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, 11/03/2022

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