

PRODUCT INFORMATION



Fmoc-Arg(Pbf)-OH

Item No. 36458

CAS Registry No.: 154445-77-9

Formal Name: N⁵-[[[(2,3-dihydro-2,2,4,6,7-pentamethyl-5-benzofuranyl)sulfonyl]amino]iminomethyl]-N²-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-ornithine

MF: C₃₄H₄₀N₄O₇S

FW: 648.8

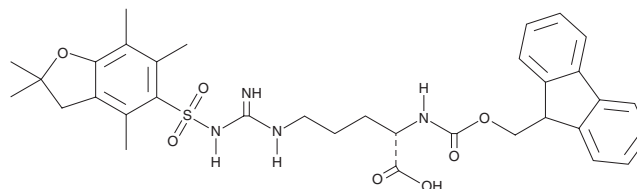
Purity: ≥98%

UV/Vis.: λ_{max}: 257 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-Arg(Pbf)-OH is supplied as a solid. A stock solution may be made by dissolving the Fmoc-Arg(Pbf)-OH in the solvent of choice, which should be purged with an inert gas. Fmoc-Arg(Pbf)-OH is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Fmoc-Arg(Pbf)-OH in ethanol is approximately 0.5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

Fmoc-Arg(Pbf)-OH is an amino acid-containing building block.¹ It has been used in the solid-phase synthesis of peptides containing L-arginine (Item No. 23703).

Reference

1. de la Torre, B.G., Kumar, A., Alhassan, M., *et al.* Successful development of a method for the incorporation of Fmoc-Arg(Pbf)-OH in solid-phase peptide synthesis using N-butylpyrrolidinone (NBP) as solvent. *Green Chem.* **10**(22), 3162 (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, 05/24/2023

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