

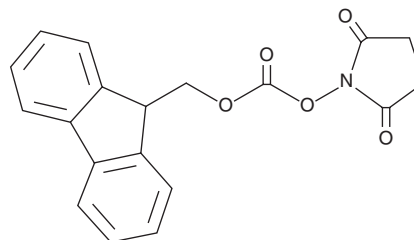
PRODUCT INFORMATION



Fmoc-Succinimide

Item No. 13861

CAS Registry No.: 82911-69-1
Formal Name: 2,5-dioxo-1-pyrrolidiny 9H-fluoren-9-ylmethyl ester-carbonic acid
Synonyms: N-(9-Fluorenylmethoxycarbonyloxy)succinimide, Fmoc-OSu
MF: C₁₉H₁₅NO₅
FW: 337.3
Purity: ≥98%
UV/Vis.: λ_{max}: 257 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-Succinimide (Fmoc-OSu) is supplied as a crystalline solid. A stock solution may be made by dissolving the Fmoc-OSu in the solvent of choice, which should be purged with an inert gas. Fmoc-OSu is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Fmoc-OSu in these solvents is approximately 20 and 25 mg/ml, respectively.

Description

Fmoc-OSu is a mixed carbamate reagent for the selective preparation of N-(9-fluorenylmethoxycarbonyl) derivatives of hydroxyamino acids and hydroxyamino acid esters in high yields.¹ Fmoc-OSu is a more preferable alternative to the use of the chloroformate, Fmoc-Cl, because it eliminates the formation of dipeptides.¹ Fmoc-OSu has also been employed in the synthesis of glycopeptides.

Reference

1. Paquet, A. Introduction of 9-fluorenylmethoxycarbonyl, trichloroethoxycarbonyl, and benzyloxycarbonyl amine protecting groups into O-unprotected hydroxyamino acids using succinimidyl carbonates. *Can. J. Chem.* **60**, 976-980 (1982).

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