

# PRODUCT INFORMATION

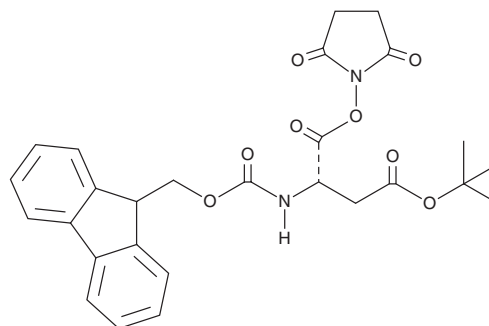


## Fmoc-Asp(OtBu)-OSu

Item No. 30537

**CAS Registry No.:** 78553-23-8  
**Formal Name:** 4-(1,1-dimethylethyl)N-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-aspartic acid, 1-(2,5-dioxo-1-pyrrolidinyl) ester

**MF:** C<sub>27</sub>H<sub>28</sub>N<sub>2</sub>O<sub>8</sub>  
**FW:** 508.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 265 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-Asp(OtBu)-OSu is supplied as a solid. A stock solution may be made by dissolving the Fmoc-Asp(OtBu)-OSu in the solvent of choice, which should be purged with an inert gas. Fmoc-Asp(OtBu)-OSu is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Fmoc-Asp(OtBu)-OSu in these solvents is approximately 20 and 10 mg/ml, respectively.

Fmoc-Asp(OtBu)-OSu is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Fmoc-Asp(OtBu)-OSu should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fmoc-Asp(OtBu)-OSu has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Fmoc-Asp(OtBu)-OSu is an amino acid-containing building block.<sup>1,2</sup> It has been used in the synthesis of peptaibols, as well as adenosine derivatives that inhibit the human ADP-ribosylhydrolase macrodomain-containing protein 1 (MacroD1).

### References

1. Jadhav, S.V., Bandyopadhyay, A., Benke, S.N., *et al.* A facile synthesis and crystallographic analysis of *N*-protected β-amino alcohols and short peptaibols. *Org. Biomol. Chem.* **99(11)**, 4182-4187 (2011).
2. Zhang, Y., Jumppanen, M., Maksimainen, M.M., *et al.* Adenosine analogs bearing phosphate isosteres as human MDO1 ligands. *Bioorg. Med. Chem.* **26(8)**, 1588-1597 (2018).

#### 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