

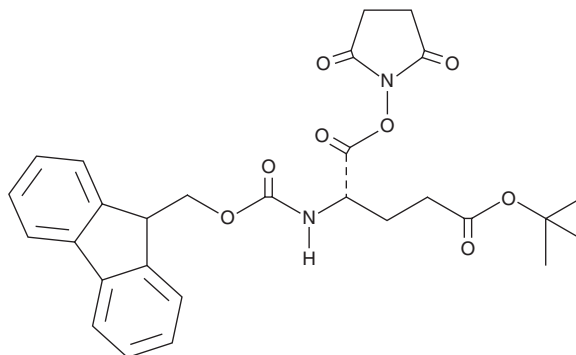
# PRODUCT INFORMATION



## Fmoc-Glu(OtBu)-OSu

Item No. 30545

**CAS Registry No.:** 101214-22-6  
**Formal Name:** N-[(9H-fluoren-9-ylmethoxy) carbonyl]-L-glutamic acid, 5-(1,1-dimethylethyl) 1-(2,5-dioxo-1-pyrrolidinyl) ester  
**MF:** C<sub>28</sub>H<sub>30</sub>N<sub>2</sub>O<sub>8</sub>  
**FW:** 522.6  
**Purity:** ≥95%  
**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-Glu(OtBu)-OSu is supplied as a crystalline solid. A stock solution may be made by dissolving the Fmoc-Glu(OtBu)-OSu in the solvent of choice, which should be purged with an inert gas. Fmoc-Glu(OtBu)-OSu is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Fmoc-Glu(OtBu)-OSu in these solvents is approximately 30 mg/ml.

Fmoc-Glu(OtBu)-OSu is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Fmoc-Glu(OtBu)-OSu should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fmoc-Glu(OtBu)-OSu has a solubility of approximately 0.33 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-Glu(OtBu)-OSu is an amino acid-containing building block.<sup>1</sup> It has been used in the synthesis of bis- and tris-branched carboxyl group-terminal PEGylating reagents. Fmoc-Glu(OtBu)-OSu has also been used in the synthesis of N<sub>3</sub>S chelators.<sup>2</sup>

### References

1. Felix, A.M. and Bandaranayake, R. Synthesis of bis- and tris-branched COOH-terminal pegylating reagents: Conjugation to NH<sub>2</sub>-terminal peptides. *J. Pept. Res.* **65**(1), 71-76 (2005).
2. Zhu, Z., Wang, Y., Zhang, Y., *et al.* A novel and simplified route to the synthesis of N3S chelators for <sup>99m</sup>Tc labeling. *Nucl. Med. Biol.* **28**(6), 703-708 (2001).

#### 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