

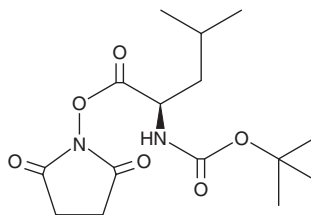
# PRODUCT INFORMATION



## Boc-D-Leu-OSu

Item No. 30543

**CAS Registry No.:** 60111-76-4  
**Formal Name:** N-[(1,1-dimethylethoxy)carbonyl]-D-leucine, 2,5-dioxo-1-pyrrolidiny ester  
**MF:** C<sub>15</sub>H<sub>24</sub>N<sub>2</sub>O<sub>6</sub>  
**FW:** 328.4  
**Purity:** ≥95%  
**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

Boc-D-Leu-OSu is supplied as a solid. A stock solution may be made by dissolving the Boc-D-Leu-OSu in the solvent of choice, which should be purged with an inert gas. Boc-D-Leu-OSu is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of Boc-D-Leu-OSu in these solvents is approximately 30 mg/ml.

### Description

Boc-D-Leu-OSu is an amino acid-containing building block.<sup>1,2</sup> It has been used in the synthesis of peptide gastrin antagonists, as well as analogs of the peptide antibiotic gramicidin S.

### References

1. Rodriguez, M., Fulcrand, P., Laur, J., et al. Synthesis of gastrin antagonists, analogues of the C-terminal tetrapeptide of gastrin, by introduction of a β-homo residue. *J. Med. Chem.* **32(3)**, 522-528 (1989).
2. Sato, K., Yamaguchi, Y., and Nagai, U. Design and synthesis of gramicidin S analogs with high antibiotic activity. *Bull. Chem. Soc. Jpn.* **86(1)**, 112-120 (2013).

#### 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