

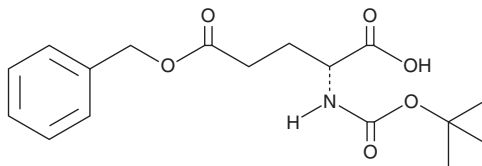
# PRODUCT INFORMATION



## Boc-Glu-OBzl

Item No. 30464

CAS Registry No.: 30924-93-7  
Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-L-glutamic acid, 1-(phenylmethyl) ester  
MF:  $C_{17}H_{23}NO_6$   
FW: 337.4  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Boc-Glu-OBzl is supplied as a solid. A stock solution may be made by dissolving the Boc-Glu-OBzl in the solvent of choice, which should be purged with an inert gas. Boc-Glu-OBzl is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Boc-Glu-OBzl in these solvents is approximately 30, 10, and 5 mg/ml, respectively.

Boc-Glu-OBzl is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Boc-Glu-OBzl should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Boc-Glu-OBzl has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Boc-Glu-OBzl is an amino acid building block.<sup>1,2</sup> It has been used in the synthesis of peptide-based inhibitors of human caspases and human rhinovirus (HRV) 3C protease that have enzyme inhibitory activity *in vitro*.

### References

1. Garcia-Calvo, M., Peterson, E.P., Leiting, B., *et al.* Inhibition of human caspases by peptide-based and macromolecular inhibitors. *J. Biol. Chem.* **273**(49), 32608-32613 (1998).
2. Dragovich, P.S., Webber, S.E., Babine, R.E., *et al.* Structure-based design, synthesis, and biological evaluation of irreversible human rhinovirus 3C protease inhibitors. 2. peptide structure-activity studies. *J. Med. Chem.* **41**(15), 2819-2934 (1998).

#### 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