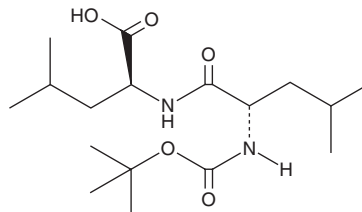


PRODUCT INFORMATION

Boc-Leu-Leu-OH

Item No. 30499

CAS Registry No.: 73401-65-7
Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-L-leucyl-L-leucine
Synonym: Leu-02
MF: C₁₇H₃₂N₂O₅
FW: 344.5
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

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

Boc-Leu-Leu-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Boc-Leu-Leu-OH should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Boc-Leu-Leu-OH has a solubility of approximately 0.25 mg/ml in a 1:3 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-Leu-Leu-OH is a peptide building block.¹ It has been used in the synthesis of a peptide-based curcumin derivative with anticancer activity *in vitro*. It has also been used in the synthesis of fluorescent peptides containing naphthalene and protoporphyrin IX.²

References

1. Ding, L., Ma, S., Lou, H., *et al.* Synthesis and biological evaluation of curcumin derivatives with water-soluble groups as potential antitumor agents: An *in vitro* investigation using tumor cell lines. *Molecules* **20**(12), 21501-21514 (2015).
2. Pispisa, B., Venanazi, M., Palleschi, A., *et al.* Intramolecular electronic energy transfer between naphthalene and protoporphyrin chromophores bound to sequential peptides. *Macromol.* **27**(26), 7800-7808 (1994).

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, 11/14/2022

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