

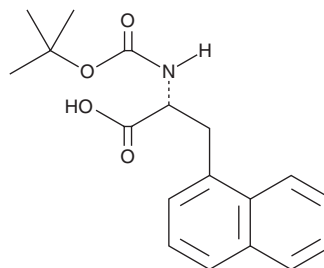
PRODUCT INFORMATION



Boc-D-1-Nal-OH

Item No. 36545

CAS Registry No.: 76932-48-4
Formal Name: α R-[[[(1,1-dimethylethoxy)carbonyl]amino]-1-naphthalenepropanoic acid
MF: $C_{18}H_{21}NO_4$
FW: 315.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 225 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Boc-D-1-Nal-OH is supplied as a solid. A stock solution may be made by dissolving the Boc-D-1-Nal-OH in the solvent of choice, which should be purged with an inert gas. Boc-D-1-Nal-OH is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Boc-D-1-Nal-OH in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

Description

Boc-D-1-Nal-OH is a building block.^{1,2} It has been used in the synthesis of melanocortin receptor 3 (MC3R) agonists and antimicrobial peptidomimetics.

References

1. Gozhina, O.V., Svendsen, J.B., and Lejon, T. Synthesis and antimicrobial activity of α -aminoboronic-containing peptidomimetics. *J. Pept. Sci.* **20(1)**, 20-24 (2014).
2. Doering, S.R., Freeman, K., Debevec, G., *et al.* Discovery of nanomolar melanocortin-3 receptor (MC3R)-selective small molecule pyrrolidine bis-cyclic guanidine agonist compounds via a high-throughput "unbiased" screening campaign. *J. Med. Chem.* **64(9)**, 5577-5592 (2021).

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, 05/31/2023

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