

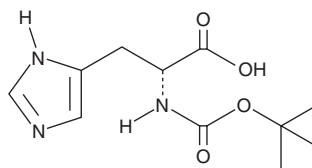
PRODUCT INFORMATION



N^α-Boc-D-Histidine

Item No. 30501

CAS Registry No.: 50654-94-9
Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-D-histidine
Synonyms: N-(*tert*-Butoxycarbonyl)-D-histidine, N-(*t*-Butoxycarbonyl)-D-histidine, N^α-*tert*-Butoxycarbonyl-D-histidine
MF: C₁₁H₁₇N₃O₄
FW: 255.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214 nm
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

N^α-Boc-D-Histidine is supplied as a crystalline solid. A stock solution may be made by dissolving the N^α-Boc-D-Histidine in the solvent of choice, which should be purged with an inert gas. N^α-Boc-D-Histidine is slightly soluble in ethanol and DMSO.

Description

N^α-Boc-D-Histidine is an amino acid building block.¹⁻³ It has been used in the synthesis of antimicrobial peptides, thyrotropin-releasing hormone enantiomers, and proteolysis-resistant peptide renin inhibitors.

References

1. Chen, Y., Mant, C.T., and Hodges, R.S. Preparative reversed-phase high-performance liquid chromatography collection efficiency for an antimicrobial peptide on columns of varying diameters (1 mm to 9.4 mm I.D.). *J. Chromatogr. A* **1140**(1-2), 112-120 (2007).
2. Paruszewski, R. and Strzalkowska, M. New enzymatically stable inhibitors of human renin. Part II. *Pol. J. Chem.* **64**(1-6), 149-156 (1990).
3. Flouret, G., Morgan, R.J., Gendrich, R., *et al.* Synthesis of the thyrotropin-releasing hormone enantiomer and some diastereoisomers and *in vitro* studies of their biological activity. *J. Med. Chem.* **16**(10), 1137-1140 (1973).

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