

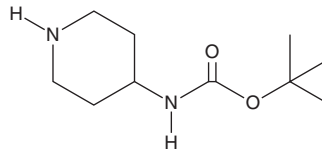
# PRODUCT INFORMATION



## 4-(N-Boc-amino)piperidine

Item No. 30398

**CAS Registry No.:** 73874-95-0  
**Formal Name:** N-4-piperidinyI-carbamic acid, 1,1-dimethylethyl ester  
**Synonyms:** 4-(t-Butoxycarbonylamino)piperidine, 4-(tert-Butoxycarbonylamino)piperidine  
**MF:** C<sub>10</sub>H<sub>20</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 200.3  
**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

4-(N-Boc-amino)piperidine is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-(N-Boc-amino)piperidine in the solvent of choice, which should be purged with an inert gas. 4-(N-Boc-amino)piperidine is soluble in chloroform at a concentration of approximately 30 mg/ml.

### Description

4-(N-Boc-amino)piperidine is an organic building block.<sup>1,2</sup> It has been used in the synthesis of aminopiperidine antiviral chemokine (C-C motif) receptor 5 (CCR5) antagonists and antibacterial agents.

### References

1. Burrows, J.N., Cumming, J.G., Fillery, S.M., *et al.* Modulators of the human CCR5 receptor. Part 1: Discovery and initial SAR of 1-(3,3-diphenylpropyl)-piperidinyI amides and ureas. *Bioorg. Med. Chem. Lett.* **15**(1), 25-28 (2005).
2. Reck, F., Alm, R., Brassil, P., *et al.* Novel N-linked aminopiperidine inhibitors of bacterial topoisomerase type II: Broad-spectrum antibacterial agents with reduced hERG activity. *J. Med. Chem.* **54**(22), 7834-7847 (2011).

#### 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, 10/27/2022

#### 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](http://WWW.CAYMANCHEM.COM)