

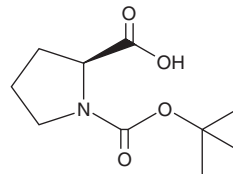
# PRODUCT INFORMATION



## N-Boc-L-Proline

Item No. 30386

**CAS Registry No.:** 15761-39-4  
**Formal Name:** (2S)-1,2-pyrrolidinedicarboxylic acid, 1-(1,1-dimethylethyl) ester  
**Synonyms:** BOC-L-Proline, 1-*tert*-butyloxycarbonyl-L-Proline, NSC 164660  
**MF:** C<sub>10</sub>H<sub>17</sub>NO<sub>4</sub>  
**FW:** 215.2  
**Purity:** ≥98%  
**Supplied as:** A 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-Boc-L-proline is supplied as a solid. A stock solution may be made by dissolving the N-Boc-L-proline in the solvent of choice, which should be purged with an inert gas. N-Boc-L-proline is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-Boc-L-proline in ethanol and DMSO is approximately 15 mg/ml and approximately 20 mg/ml in DMF.

N-Boc-L-proline is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-Boc-L-proline should first be dissolved in DMF and then diluted with the aqueous buffer of choice. N-Boc-L-proline has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

N-Boc-L-proline is a synthetic intermediate.<sup>1,2</sup> It has been used in the synthesis of enantioselective catalysts for aldol reactions and hepatitis C virus (HCV) NS5A inhibitors.

### References

1. Berkessel, A., Koch, B., and Lex, J. Proline-derived N-sulfonylcarboxamides: Readily available, highly enantioselective and versatile catalysts for direct Aldol reactions. *Adv. Synth. Catal.* **346(9-10)**, 1141-1146 (2004).
2. Shi, J., Zhou, L., Amblard, F., *et al.* Synthesis and biological evaluation of new potent and selective HCV NS5A inhibitors. *Bioorg. Med. Chem.* **22(10)**, 3488-3491 (2012).

#### 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