

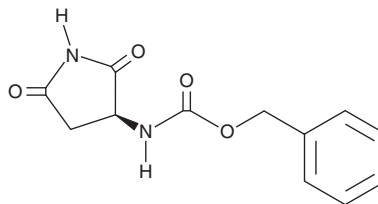
PRODUCT INFORMATION



(S)-3-N-Cbz-amino-succinimide

Item No. 28568

CAS Registry No.: 60846-91-5
Formal Name: [(3S)-2,5-dioxo-3-pyrrolidinyl]-
carbamic acid, phenylmethyl ester
MF: C₁₂H₁₂N₂O₄
FW: 248.2
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

(S)-3-N-Cbz-amino-succinimide is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-3-N-Cbz-amino-succinimide in the solvent of choice, which should be purged with an inert gas. (S)-3-N-Cbz-amino-succinimide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (S)-3-N-Cbz-amino-succinimide in these solvents is approximately 25 mg/ml.

(S)-3-N-Cbz-amino-succinimide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (S)-3-N-Cbz-amino-succinimide should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (S)-3-N-Cbz-amino-succinimide 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

(S)-3-N-Cbz-amino-succinimide is an anticonvulsant.¹ It inhibits tonic convulsions induced by pentylenetetrazole (Item No. 18682) and maximal electric shock (MES) in mice (ED₅₀s = 78.1 and 103 mg/kg, respectively).

Reference

1. Lee, J., Son, K., Jung, K., *et al.* The effect of N-substituted alkyl groups on the anticonvulsant activities of N-Cbz-α-amino-N-alkylsuccinimides. *Arch. Pharm. Res.* **20(1)**, 53-57 (1997).

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, 10/04/2022

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