

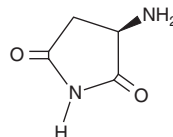
PRODUCT INFORMATION



(R)-Aspartimide

Item No. 31699

CAS Registry No.: 404887-22-5
Formal Name: 3R-amino-2,5-pyrrolidinedione
MF: $C_4H_6N_2O_2$
FW: 114.1
Purity: $\geq 95\%$
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

(R)-Aspartimide is supplied as a solid. A stock solution may be made by dissolving the (R)-aspartimide in the solvent of choice, which should be purged with an inert gas. (R)-Aspartimide is soluble in methanol and DMSO.

Description

(R)-Aspartimide is a byproduct formed during solid-phase peptide synthesis (SPPS) of aspartic acid-containing peptides.¹ Aspartimide is formed by cyclization of aspartic acid during Fmoc removal or peptide coupling.²

References

1. Ruczyński, J., Lewandowska, B., Mucha, P., *et al.* Problem of aspartimide formation in Fmoc-based solid-phase peptide synthesis using Dmab group to protect side chain of aspartic acid. *J. Pept. Sci.* **14(3)**, 335-341 (2008).
2. Neumann, K., Farnung, J., Baldauf, S., *et al.* Prevention of aspartimide formation during peptide synthesis using cyanosulfurylides as carboxylic acid-protecting groups. *Nat. Commun.* **11(1)**, 982 (2020).

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, 11/10/2022

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