

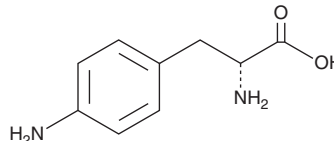
PRODUCT INFORMATION



4-Amino-D-phenylalanine

Item No. 41971

CAS Registry No.: 102281-45-8
Formal Name: (R)-2-amino-3-(4-aminophenyl)propanoic acid
Synonyms: *p*-Amino-D-phenylalanine, *para*-Amino-D-phenylalanine
MF: C₉H₁₂N₂O₂
FW: 180.2
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

4-Amino-D-phenylalanine is supplied as a solid. A stock solution may be made by dissolving the 4-amino-D-phenylalanine in the solvent of choice, which should be purged with an inert gas. 4-Amino-D-phenylalanine is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

4-Amino-D-phenylalanine is a synthetic amino acid and building block.¹ It has been used in the synthesis of chemokine (C-X-C motif) receptor 4 (CXCR4) antagonists.

Reference

1. Tamamura, H., Esaka, A., Ogawa, T., *et al.* Structure-activity relationship studies on CXCR4 antagonists having cyclic pentapeptide scaffolds. *Org. Biomol. Chem.* **3**(24), 4392-4394 (2005).

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, 10/17/2024

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