

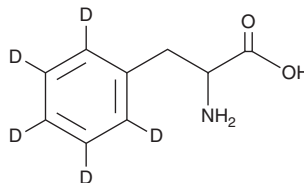
PRODUCT INFORMATION



DL-Phenyl-d₅-alanine

Item No. 34845

CAS Registry No.: 284664-89-7
Formal Name: phenylalanine-2,3,4,5,6-d₅
Synonym: (±)-Phenyl-d₅-alanine
MF: C₉H₆D₅NO₂
FW: 170.2
Chemical Purity: ≥98% (Phenylalanine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

DL-Phenyl-d₅-alanine is intended for use as an internal standard for the quantification of phenylalanine by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

DL-Phenylalanine is a racemic mixture of the essential amino acid L-phenylalanine (Item No. 31498) and its enantiomer D-phenylalanine (Item No. 31194). It has been used as a precursor in the synthesis of sulfonamide phenylalanine derivatives with antibacterial, antifungal, and anticancer activity.¹

Reference

1. Devi, K. and Awasthi, P. Sulfonamide phenylalanine (SPA) series of analogues as an antibacterial, antifungal, anticancer agents along with p53 tumor suppressor-DNA complex inhibitor - part 1. *J. Biomol. Struct. Dyn.* **38(14)**, 4018-4097 (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, 12/15/2022

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