

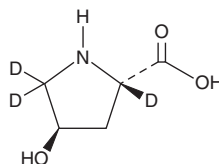
PRODUCT INFORMATION



trans-4-hydroxy L-Proline-d₃

Item No. 43922

CAS Registry No.: 1356016-86-8
Formal Name: (4R)-4-hydroxy-L-proline-2,5,5-d₃
Synonyms: (S)-(-)-*trans*-4-Hydroxyproline-d₃,
4(R)-Hydroxy-2(S)-pyrrolidinecarboxylic acid-d₃
MF: C₅H₆D₃NO₃
FW: 134.1
Purity: ≥98% (*trans*-4-hydroxy L-proline)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤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

trans-4-hydroxy L-Proline-d₃ is intended for use as an internal standard for the quantification of *trans*-4-hydroxy L-proline (Item No. 26061) 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

trans-4-hydroxy L-Proline is a non-essential amino acid that is commonly used as a chiral building block in the organic synthesis of various compounds.¹ It is a constituent of collagen and elastin and has been found in many bacterial secondary metabolites, including echinocandins.

Reference

1. Remuzon, P. *Trans*-4-hydroxy-L-proline, a useful and versatile chiral starting block. *Tetrahedron* **44**(52), 13803-13835 (1996).

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, 07/31/2025

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