

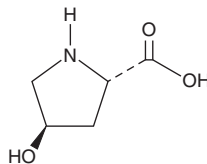
# PRODUCT INFORMATION



## *trans*-4-hydroxy L-Proline

Item No. 26061

**CAS Registry No.:** 51-35-4  
**Formal Name:** (4R)-hydroxy-L-proline  
**Synonyms:** (S)-(-)-*trans*-4-Hydroxyproline, 4(R)-Hydroxy-2(S)-pyrrolidinecarboxylic acid, NSC 46704  
**MF:** C<sub>5</sub>H<sub>9</sub>NO<sub>3</sub>  
**FW:** 131.1  
**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

*trans*-4-hydroxy L-Proline is supplied as a crystalline solid. Aqueous solutions of *trans*-4-hydroxy L-proline can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of *trans*-4-hydroxy L-proline in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### 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.<sup>1</sup> 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, 09/29/2022

#### 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