

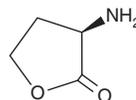
PRODUCT INFORMATION



D-Homoserine lactone

Item No. 31695

CAS Registry No.: 51744-82-2
Formal Name: 3R-aminodihydro-2(3H)-furanone
Synonym: D-HSL
MF: C₄H₇NO₂
FW: 101.1
Purity: ≥95%
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

D-Homoserine lactone is supplied as a solid. A stock solution may be made by dissolving the D-homoserine lactone in the solvent of choice, which should be purged with an inert gas. D-Homoserine lactone is soluble in the organic solvent DMSO. While D-homoserine lactone is also soluble in ethanol and other primary alcohols, their use is not recommended as they have been shown to open the lactone ring.

Description

D-homoserine lactone is an enantiomer of L-homoserine lactone (Item No. 31492) and an inhibitor of serine hydroxymethyltransferase (SHMT; K_i = 11 mM for the rabbit enzyme).¹

Reference

1. Wang, E.A., Kallen, R., and Walsh, C. Mechanism-based inactivation of serine transhydroxymethylases by D-fluoroalanine and related amino acids. *J. Biol. Chem.* **256(13)**, 6917-6926 (1981).

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, 12/02/2022

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