

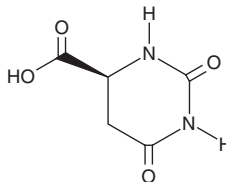
PRODUCT INFORMATION



L-Dihydroorotic Acid

Item No. 45688

CAS Registry No.: 5988-19-2
Formal Name: hexahydro-2,6-dioxo-4S-pyrimidinecarboxylic acid
Synonym: L-Dihydroorotate
MF: C₅H₆N₂O₄
FW: 158.1
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

L-Dihydroorotic acid is supplied as a solid. A stock solution may be made by dissolving the L-dihydroorotic acid in the solvent of choice, which should be purged with an inert gas. L-Dihydroorotic acid is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

L-Dihydroorotic acid is an intermediate in the *de novo* pyrimidine biosynthesis pathway.¹ It is formed from carbamoyl phosphate by dihydroorotase and converted into orotic acid by dihydroorotate dehydrogenase (DHODH).

Reference

1. Evans, D.R. and Guy, H.I. Mammalian pyrimidine biosynthesis: Fresh insights into an ancient pathway. *J. Biol. Chem.* **279**(32), 33035-33038 (2004).

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, 05/22/2026

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