

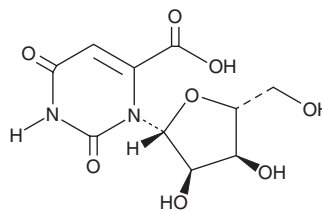
PRODUCT INFORMATION



Orotidine

Item No. 36401

CAS Registry No.: 314-50-1
Formal Name: 1,2,3,6-tetrahydro-2,6-dioxo-3-β-D-ribofuranosyl-4-pyrimidinecarboxylic acid
Synonyms: Orotic Acid riboside, 6-carboxy Uridine
MF: C₁₀H₁₂N₂O₈
FW: 288.2
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

Orotidine is supplied as a solid. A stock solution may be made by dissolving the orotidine in water. The solubility of orotidine in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Orotidine is a pyrimidine nucleoside composed of ribose and orotic acid.¹ Urine levels of orotidine are increased in rats following administration of the xanthine oxidase inhibitor allopurinol (Item No. 10012597).²

References

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. *Nucleotide biosynthesis*. Biochemistry, 5th ed., W.H. Freeman (2002).
2. Brown, G.K., Fox, R.M., and O'Sullivan, W.J. Alteration of quaternary structural behaviour of a hepatic orotate phosphoribosyltransferase-otridine-5'-phosphate decarboxylase complex in rats following allopurinol therapy. *Biochem. Pharmacol.* **21(18)**, 2469-2477 (1972).

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, 12/19/2022

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