

PRODUCT INFORMATION



3-hydroxy methyl Cefuroxime

Item No. 35079

CAS Registry No.: 56271-94-4

Formal Name: (6R,7R)-7-[[[(2Z)-2-(2-furanyl)-2-(methoxyimino)acetyl]amino]-3-(hydroxymethyl)-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid

Synonym: Descarbamoyl Cefuroxime

MF: C₁₅H₁₅N₃O₇S

FW: 381.4

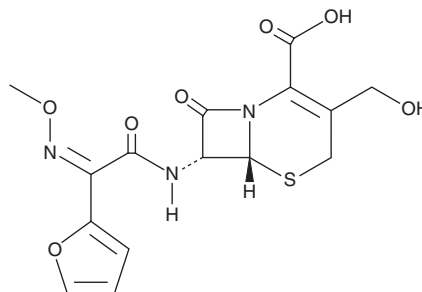
Purity: ≥95%

UV/Vis.: λ_{max}: 275 nm

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

3-hydroxy methyl Cefuroxime is supplied as a solid. A stock solution may be made by dissolving the 3-hydroxy methyl cefuroxime in the solvent of choice, which should be purged with an inert gas. 3-hydroxy methyl Cefuroxime is soluble in DMSO.

Description

3-hydroxy methyl Cefuroxime is an intermediate in the synthesis of cephalosporin antibiotic precursors.¹

Reference

1. Singh, S.K., Tiwari, R.P., and Jain, P. A novel and efficient synthesis of (6R,7R)-7-amino-3-hydroxymethyl cephalosporanic Acid: A versatile precursor of cefuroxime acid. *Synth. Comm.* **33**(14), 2475-2482 (2006).

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, 12/09/2022

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