

PRODUCT INFORMATION



Cefadroxil (hydrate)

Item No. 20365

CAS Registry No.: 66592-87-8
Formal Name: (6R,7R)-7-[[[(2R)-2-amino-2-(4-hydroxyphenyl)acetyl]amino]-3-methyl-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, monohydrate

MF: C₁₆H₁₇N₃O₅S • H₂O

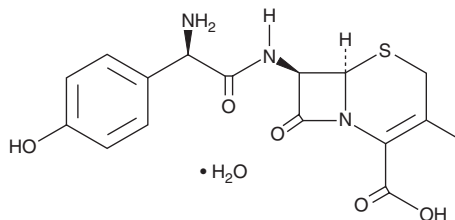
FW: 381.4

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

Cefadroxil (hydrate) is supplied as a solid. A stock solution may be made by dissolving the cefadroxil (hydrate) in the solvent of choice, which should be purged with an inert gas. Cefadroxil (hydrate) is slightly soluble in methanol (sonicated).

Cefadroxil (hydrate) is slightly soluble in aqueous solutions (warmed). To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Cefadroxil is a semisynthetic cephalosporin antibiotic.¹ It is active against a variety of Gram-positive and Gram-negative bacteria, including *B. anthracis*, *V. cholerae*, and *P. multocida* strains (MICs = 1, 2, and 4 µg/ml, respectively), among others. Cefadroxil is protective against infection by strains of *S. pyogenes*, *S. pneumoniae*, or *S. aureus* in mice with 50% protective dose (PD₅₀) values of 0.6-2.4, 18-24, and 0.3-29 mg/kg, respectively. Formulations containing cefadroxil have been used in the treatment of a variety of bacterial infections.

Reference

1. Buck, R.E. and Price, K.E. Cefadroxil, a new broad-spectrum cephalosporin. *Antimicrob. Agents Chemother.* **11**(2), 324-330 (1977).

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, 12/01/2022

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