

PRODUCT INFORMATION



Cefaclor

Item No. 23626

CAS Registry No.: 53994-73-3
Formal Name: (6R,7R)-7-[[[(2R)-2-amino-2-phenylacetyl]amino]-3-chloro-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid

MF: C₁₅H₁₄ClN₃O₄S

FW: 367.8

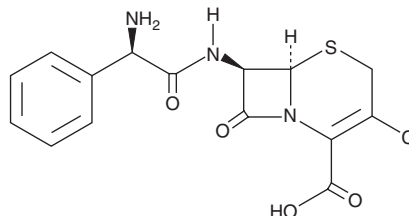
Purity: ≥95%

UV/Vis.: λ_{max}: 265 nm

Supplied as: A crystalline 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

Cefaclor is supplied as a crystalline solid. Aqueous solutions of cefaclor can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cefaclor in PBS, pH 7.2, is approximately 1.6 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cefaclor is a cephalosporin antibiotic that is active against *S. pyogenes*, *S. pneumoniae*, *S. aureus*, *P. mirabilis*, *S. typhi*, *E. coli*, and *H. influenzae* (MICs = 0.25, 0.25-2, 1-2, 1, 0.5, 1, and 2-4 µg/ml, respectively).¹ It is protective against *S. pyogenes*, *S. pneumoniae*, *S. aureus*, *P. mirabilis*, *S. typhi*, *E. coli*, and *H. influenzae* infections in mice (ED₅₀s = 0.08-30.2 mg/kg). Formulations containing cefaclor have been used in the treatment of various bacterial infections.

Reference

1. Preston, D.A. Summary of laboratory studies on the antibacterial activity of cefaclor. *Infection* 7(Suppl. 6), 557-561 (1979).

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, 10/11/2022

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