

PRODUCT INFORMATION



Cefradine

Item No. 26060

CAS Registry No.: 38821-53-3
Formal Name: (6R,7R)-7-[[[(2R)-2-amino-2-(1,4-cyclohexadien-1-yl)acetyl]amino]-3-methyl-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid

Synonym: SQ 11,436

MF: C₁₆H₁₉N₃O₄S

FW: 349.4

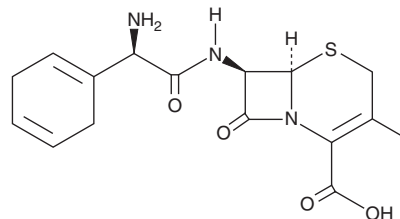
Purity: ≥90%

UV/Vis.: λ_{max}: 263 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.

Description

Cefradine is an orally bioavailable β-lactam cephalosporin antibiotic.¹ It is active against *S. pyogenes*, *E. coli*, *K. pneumoniae*, and *E. cloacae* (MICs = 0.04, 9.4, 9.4, and 6.3 μg/ml, respectively). Cefradine is also active against clinical isolates of *S. aureus* (MICs = 0.8-6 μg/ml) and *S. pyogenes* (MICs = 0.1-0.4 μg/ml), as well as *H. influenzae*, *E. coli*, and *K. pneumoniae* (MICs = 6.2-12.5 μg/ml).² It increases survival in mouse models of systemic lethal infection by *S. pyogenes*, *E. coli*, *K. pneumoniae*, or *E. cloacae* (ED₅₀s = 5, 37, 122, and 50 mg/kg, respectively), as well as by penicillin-susceptible or -resistant strains of *S. aureus* (ED₅₀s = 18 and 91 mg/kg, respectively).¹ Formulations containing cefradine have previously been used in the treatment of respiratory and urinary tract infections, skin infections, and otitis media.

References

1. Miraglia, G.J., Renz, K.J., and Gadebusch, H.H. Comparison of the chemotherapeutic and pharmacodynamic activities of cephradine, cephalothin, and cephaloridine in mice. *Antimicrob. Agents Chemother.* **3**(2), 270-273 (1973).
2. Shibl, A.M. and Durgham, S.M. *In-vitro* evaluation of a new oral cephalosporin, cefroxadine (CGP 9000). *J. Chemother.* **2**(1), 8-10 (1990).

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, 11/09/2022

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