

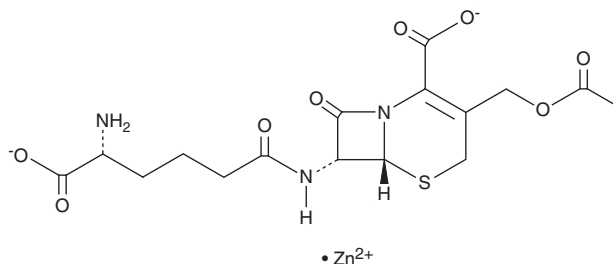
PRODUCT INFORMATION



Cephalosporin C (zinc)

Item No. 33262

CAS Registry No.: 59143-60-1
Formal Name: 3-[(acetyloxy)methyl]-7-[[[(5R)-5-amino-5-carboxy-1-oxopentyl]amino]-8-oxo-(6R,7R)-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, zinc salt
MF: C₁₆H₁₉N₃O₈S • Zn
FW: 478.8
Purity: ≥95%
UV/Vis.: λ_{max}: 258 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Cephalosporin C is a fungal metabolite and cephalosporin antibiotic that has been found in *Cephalosporium*.¹ It is active against a variety of bacteria, including *S. pyogenes*, *C. diphtheriae*, *E. coli*, and *H. influenzae* (MICs = 31, 15.6, 125, and 31 µg/ml, respectively), as well as clinical isolates of *N. gonorrhoeae* (MICs = 1.9-15.6 µg/ml).² Formulations containing cephalosporin C have been used in the treatment of various bacterial infections.

References

1. Newton, G.G. and Abaham, E.P. Cephalosporin C, a new antibiotic containing sulphur and D-alpha-aminoadipic acid. *Nature* **175(4456)**, 548 (1955).
2. Jago, M. and Heatley, N.G. Some biological properties of cephalosporin C and a derivative. *Br. J. Pharmacol. Chemother.* **16(2)**, 170-179 (1961).

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, 01/03/2023

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