

PRODUCT INFORMATION



Desacetyl Cefapirin (sodium salt)

Item No. 27578

CAS Registry No.: 104557-24-6
Formal Name: (6R-*trans*)-3-(hydroxymethyl)-8-oxo-7-[[[4-pyridinylthio]acetyl]amino]-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, monosodium salt

MF: C₁₅H₁₄N₃O₅S₂ • Na
FW: 403.4

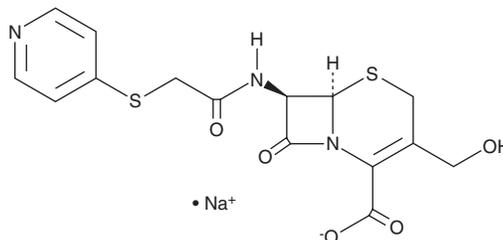
Purity: ≥95%

UV/Vis.: λ_{max}: 267 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

Desacetyl cefapirin (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the desacetyl cefapirin (sodium salt) in the solvent of choice, which should be purged with an inert gas. Desacetyl cefapirin (sodium salt) is soluble in DMSO and slightly soluble in methanol.

Desacetyl cefapirin (sodium salt) is slightly soluble in aqueous solutions. 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

Desacetyl cefapirin is an active metabolite of the cephalosporin antibiotic cefapirin (Item No. 17406).¹ Desacetyl cefapirin is active against *E. coli*, *K. pneumoniae*, *P. mirabilis*, and *S. aureus* with MIC values of 120, 24, 34, and 0.42 µg/ml, respectively.

Reference

1. Jones, R.N. and Packer, R.R. Cefotaxime, cephalothin, and cephalapirin: Antimicrobial activity and synergy studies of cephalosporins with significant in vivo desacetyl metabolite concentrations. *Diagn. Microbiol. Infect. Dis.* **2**(1), 65-68 (1984).

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, 08/10/2023

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