

PRODUCT INFORMATION



Desacetylcefotaxime (potassium salt)

Item No. 33921

Formal Name: (6R,7R)-7-[[[(2Z)-2-(2-amino-4-thiazolyl)-2-(methoxyimino)acetyl]amino]-3-(hydroxymethyl)-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, monopotassium salt

Synonym: des-CTX

MF: C₁₄H₁₄N₅O₆S₂ • K

FW: 451.5

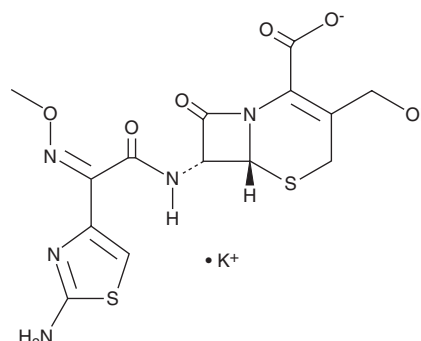
Purity: ≥95%

UV/Vis.: λ_{max}: 237 nm

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

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

Desacetylcefotaxime (potassium 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

Desacetylcefotaxime is an active metabolite of the cephalosporin antibiotic cefotaxime (Item No. 16040).¹ It is active against 60 clinical isolates derived from patients with meningitis, including *H. influenzae*, *S. pneumoniae*, *S. agalactiae*, and *N. meningitidis* (MIC₉₀^S = 0.008-0.12 µg/ml).

Reference

1. Jones, R.N. Cefotaxime and desacetylcefotaxime antimicrobial interactions. The clinical relevance of enhanced activity: A review. *Diagn. Microbiol. Infect. Dis.* **22**(1-2), 19-33 (1995).

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, 11/17/2022

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