

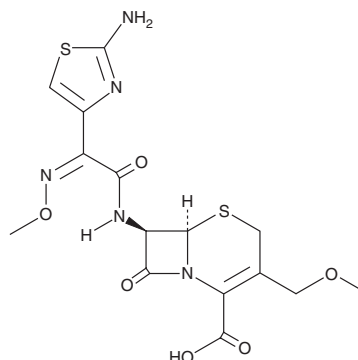
PRODUCT INFORMATION



Cefpodoxime

Item No. 16075

CAS Registry No.: 80210-62-4
Formal Name: (6R)-7R-[[[(2Z)-2-(2-amino-4-thiazolyl)-2-(methoxyimino)acetyl]amino]-3-(methoxymethyl)-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid
Synonym: R 3763
MF: C₁₅H₁₇N₅O₆S₂
FW: 427.5
Purity: ≥95%
UV/Vis.: λ_{max}: 235 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

Cefpodoxime is supplied as a crystalline solid. A stock solution may be made by dissolving the cefpodoxime in the solvent of choice, which should be purged with an inert gas. Cefpodoxime is soluble in organic solvents such as ethanol, methanol, DMSO, and acetonitrile. The solubility of cefpodoxime in ethanol, methanol, and DMSO is approximately 10 mg/ml and approximately 5 mg/ml in acetonitrile.

Description

Cefpodoxime is a metabolite of cefpodoxime proxetil, a third generation cephalosporin antibiotic, that is active against both Gram-positive and Gram-negative bacteria (MIC₉₀S = 0.015-8 µg/ml).¹

Reference

1. Angehrn, P., Hohl, P., Hubschwerlen, C., *et al.* Antibacterial properties of Ro 40-6890, a broad-spectrum cephalosporin, and its novel orally absorbable ester, Ro 41-3399. *Antimicrob. Agents Chemother.* **36(12)**, 2825-2834 (1992).

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, 12/09/2022

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