

PRODUCT INFORMATION



Cefpodoxime-d₃

Item No. 25412

CAS Registry No.: 2477791-28-7
Formal Name: (6R,7R)-7-((Z)-2-(2-aminothiazol-4-yl)-2-((methoxy-d₃)imino)acetamido)-3-(methoxymethyl)-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid

MF: C₁₅H₁₄D₃N₅O₆S₂
FW: 430.5

Chemical Purity: ≥98% (Cefpodoxime)

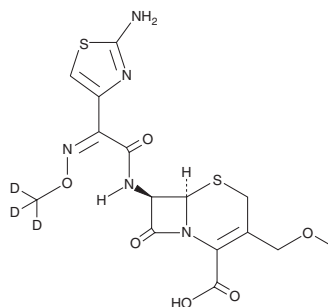
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

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

Cefpodoxime-d₃ is intended for use as an internal standard for the quantification of cefpodoxime (Item No. 16075) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Cefpodoxime-d₃ is supplied as a solid. A stock solution may be made by dissolving the cefpodoxime-d₃ in the solvent of choice, which should be purged with an inert gas. Cefpodoxime-d₃ is slightly soluble in DMSO and methanol.

Description

Cefpodoxime is a metabolite of cefpodoxime proxetil, a third generation cephalosporin antibiotic, that is active against both Gram-positive and Gram-negative bacteria (MIC_{90S} = 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, 01/13/2023

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