

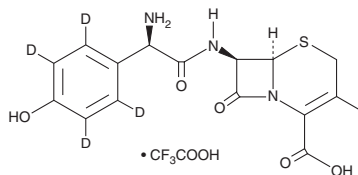
PRODUCT INFORMATION



Cefadroxil-d₄ (trifluoroacetate salt)

Item No. 33380

CAS Registry No.: 1426215-61-3
Formal Name: (6R,7R)-7-[[[(2R)-2-amino-2-(4-hydroxyphenyl-2,3,5,6-d₄)acetyl]amino]-3-methyl-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid
MF: C₁₆H₁₃D₄N₃O₅S • CF₃COOH
FW: 481.4
Chemical Purity: ≥95% (Cefadroxil)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤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

Cefadroxil-d₄ (trifluoroacetate salt) is intended for use as an internal standard for the quantification of cefadroxil (Item No. 20365) 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).

Cefadroxil-d₄ (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the cefadroxil-d₄ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Cefadroxil-d₄ (trifluoroacetate salt) is soluble in methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Cefadroxil is a semisynthetic cephalosporin antibiotic.¹ It is active against a variety of Gram-positive and Gram-negative bacteria, including *B. anthracis*, *V. cholerae*, and *P. multocida* strains (MICs = 1, 2, and 4 µg/ml, respectively), among others. Cefadroxil is protective against infection by strains of *S. pyogenes*, *S. pneumoniae*, or *S. aureus* in mice with 50% protective dose (PD₅₀) values of 0.6-2.4, 18-24, and 0.3-29 mg/kg, respectively. Formulations containing cefadroxil have been used in the treatment of a variety of bacterial infections.

Reference

1. Buck, R.E. and Price, K.E. Cefadroxil, a new broad-spectrum cephalosporin. *Antimicrob. Agents Chemother.* **11**(2), 324-330 (1977).

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, 10/05/2022

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