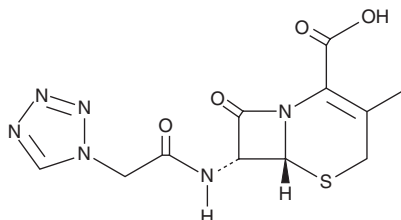


PRODUCT INFORMATION

7-[1-(1H)-Tetrazolylacetamido]desacetoxycephalosporanic Acid

Item No. 36228

CAS Registry No.: 56842-77-4
Formal Name: (6R-*trans*)-3-methyl-8-oxo-7-[(1H-tetrazol-1-ylacetyl)amino]-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid
Synonym: Cefazolin Impurity H
MF: C₁₁H₁₂N₆O₄S
FW: 324.3
Purity: ≥95%
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

7-[1-(1H)-Tetrazolylacetamido]desacetoxycephalosporanic acid is supplied as a solid. A stock solution may be made by dissolving the 7-[1-(1H)-tetrazolylacetamido]desacetoxycephalosporanic acid in the solvent of choice, which should be purged with an inert gas. 7-[1-(1H)-Tetrazolylacetamido]desacetoxycephalosporanic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 7-[1-(1H)-tetrazolylacetamido]desacetoxycephalosporanic acid in these solvents is approximately 10 mg/ml.

Description

7-[1-(1H)-Tetrazolylacetamido]desacetoxycephalosporanic acid is a potential impurity in commercial preparations of the cephalosporin antibiotic cefazolin.¹

Reference

1. Liu, S., Yao, S., Zhang, H., *et al.* Determination of relative response factors of cefazolin impurities by quantitative NMR. *AAPS PharmSciTech* **18**(6), 1895-1900 (2017).

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, 01/03/2023

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