

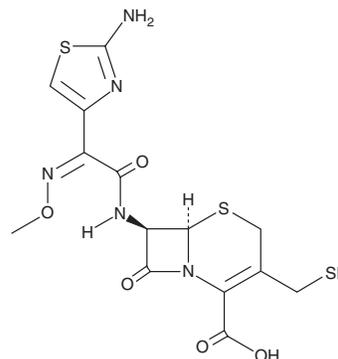
PRODUCT INFORMATION



Desfuoylceftiofur

Item No. 20360

CAS Registry No.: 120882-22-6
Formal Name: (6R,7R)-[[[(2Z)-2-(2-amino-4-thiazolyl)-2-(methoxyimino)acetyl]amino]-3-(mercaptomethyl)-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid
MF: C₁₄H₁₅N₅O₅S₃
FW: 429.5
Purity: ≥80%
UV/Vis.: λ_{max}: 236 nm
Supplied as: A solid
Storage: -80°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Desfuoylceftiofur is supplied as a solid. A stock solution may be made by dissolving the desfuoylceftiofur in the solvent of choice, which should be purged with an inert gas. Desfuoylceftiofur is slightly soluble in DMSO and methanol (sonicated).

Description

Desfuoylceftiofur is an active metabolite of the broad-spectrum cephalosporin antibiotic ceftiofur (Item No. 27301).¹ It is formed from ceftiofur by bovine kidney and liver esterases.² Desfuoylceftiofur is active against 539 veterinary isolates of Gram-negative bacteria and Gram-positive cocci bacteria (MIC_{50S} = 0.0078-1 and 0.015-4 µg/ml, respectively).

References

1. Salmon, S.A., Watts, J.L., and Yancey Jr., R.J. In vitro activity of ceftiofur and its primary metabolite, desfuoylceftiofur, against organisms of veterinary importance. *J. Vet. Diagn. Invest.* **8(3)**, 332-336 (1996).
2. Olson, S.C., Beconi-Barker, M.G., Smith, E.B., et al. In vitro metabolism of ceftiofur in bovine tissues. *J. Vet. Pharmacol. Ther.* **21(2)**, 112-120 (1998).

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, 08/05/2021

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