

PRODUCT INFORMATION

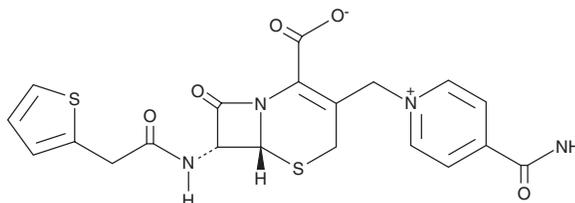


Cefalonium

Item No. 34468

CAS Registry No.: 5575-21-3
Formal Name: 4-(aminocarbonyl)-1-[[[(6R,7R)-2-carboxy-8-oxo-7-[[2-(2-thienyl)acetyl]amino]-5-thia-1-azabicyclo[4.2.0]oct-2-en-3-yl]methyl]-pyridinium, inner salt

Synonym: Cephalonium
MF: C₂₀H₁₈N₄O₅S₂
FW: 458.5
Purity: ≥98%
UV/Vis.: λ_{max}: 266 nm
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

Cefalonium is supplied as a solid. A stock solution may be made by dissolving the cefalonium in the solvent of choice, which should be purged with an inert gas. Cefalonium is soluble in the organic solvent DMSO. The solubility of cefalonium in DMSO approximately 25 mg/ml.

Cefalonium is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cefalonium should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cefalonium has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cefalonium is a cephalosporin antibiotic.¹ It is active against pooled isolates of *S. aureus* from the milk of lactating dairy cows with subclinical mastitis (MIC₉₀ = 2 µg/ml). Formulations containing cefalonium have previously been used in dry cow therapy in the treatment and prevention of mastitis.

Reference

1. Moroni, P., Pisoni, G., Antonini, M., *et al.* Short communication: Antimicrobial drug susceptibility of *Staphylococcus aureus* from subclinical bovine mastitis in Italy. *J. Dairy Sci.* **89**(8), 2973-2976 (2006).

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, 11/22/2022

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