

# PRODUCT INFORMATION

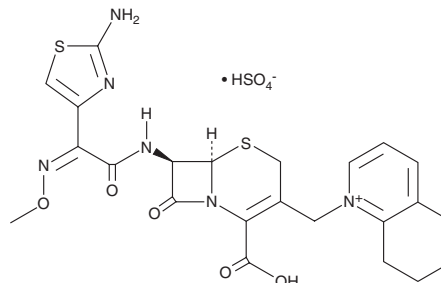


## Cefquinome (sulfate)

Item No. 34266

**CAS Registry No.:** 118443-89-3  
**Formal Name:** 1-[[[(6R,7R)-7-[[[(2Z)-2-(2-amino-4-thiazolyl)-2-(methoxyimino)acetyl]amino]-2-carboxy-8-oxo-5-thia-1-azabicyclo[4.2.0]oct-2-en-3-yl]methyl]-5,6,7,8-tetrahydro-quinolinium, monosulfate

**Synonym:** HR 111V sulfate  
**MF:** C<sub>23</sub>H<sub>25</sub>N<sub>6</sub>O<sub>5</sub>S<sub>2</sub> • HSO<sub>4</sub><sup>-</sup>  
**FW:** 626.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 267 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

Cefquinome (sulfate) is supplied as a solid. A stock solution may be made by dissolving the cefquinome (sulfate) in the solvent of choice, which should be purged with an inert gas. Cefquinome (sulfate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of cefquinome (sulfate) in these solvents is approximately 30 and 10 mg/ml, respectively.

Cefquinome (sulfate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cefquinome (sulfate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cefquinome (sulfate) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Cefquinome is a broad-spectrum cephalosporin antibiotic.<sup>1</sup> It is active against clinical isolates of various Gram-positive and Gram-negative bacteria, including *Salmonella*, *E. coli*, and methicillin-resistant *S. aureus* (MRSA; MICs = 0.06-0.5, 0.015-0.5, and 1-16 µg/ml, respectively). Intramammary administration of cefquinome (75 mg/animal) is curative in a bovine model of *E. coli*-induced mastitis.<sup>2</sup> Formulations containing cefquinome have been used in the treatment of coliform mastitis in veterinary medicine.

### References

1. Chin, N.-X., Gu, J.-W., Fang, W., *et al.* In vitro activity of cefquinome, a new cephalosporin, compared with other cephalosporin antibiotics. *Diagn. Microbiol. Infect. Dis.* **15**(4), 331-337 (1992).
2. Shpigel, N.Y., Levin, D., Winkler, M., *et al.* Efficacy of cefquinome for treatment of cows with mastitis experimentally induced using *Escherichia coli*. *J. Dairy Sci.* **80**(2), 318-323 (1996).

#### 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, 11/10/2022

#### 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