

PRODUCT INFORMATION

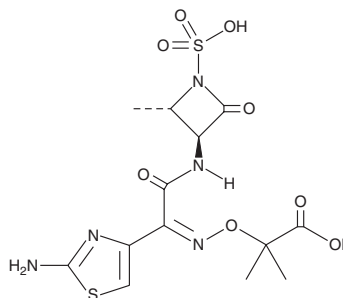


Aztreonam

Item No. 19784

CAS Registry No.: 78110-38-0
Formal Name: 2-[[[(Z)-[1-(2-amino-4-thiazolyl)-2-[(2S,3S)-2-methyl-4-oxo-1-sulfo-3-azetidiny] amino]-2-oxoethylidene]amino]oxy]-2-methyl-propanoic acid

Synonym: SQ 26,776
MF: C₁₃H₁₇N₅O₈S₂
FW: 435.4
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 260 nm
Supplied as: A crystalline 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

Aztreonam is supplied as a crystalline solid. A stock solution may be made by dissolving the aztreonam in the solvent of choice, which should be purged with an inert gas. Aztreonam is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of aztreonam in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of aztreonam can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of aztreonam in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Aztreonam is a synthetic β-lactam antibiotic of the monobactam class.^{31705} It is active against clinical isolates of *E. coli*, *S. marcescens*, *P. aeruginosa*, *Klebsiella* species, *Proteus* species, and *Enterobacter* species (MICs = <0.1, 1.2, 5.2, 3.1, <0.1, and 19.3 μg/ml, respectively) but is inactive against *S. pneumoniae*, *B. fragilis*, and *S. faecalis* (MICs = >100 μg/ml for all). It interferes with peptidoglycan synthesis in the bacterial cell wall of the Gram-negative bacteria *E. coli*, *P. vulgaris*, *E. cloacae*, *K. pneumoniae*, and *P. aeruginosa* by inhibiting penicillin-binding protein 3 (PBP3; IC₁₀₀ = 0.1 μg/ml for all). It increases survival in mice with systemic Gram-negative bacterial infections (ED₅₀s = 0.1-24.7 mg/kg). Formulations containing aztreonam have been used in the treatment of *P. aeruginosa* infections in individuals with cystic fibrosis.

References

1. Sykes, R. B., Bonner, D. P., Bush, K., et al. *Antimicrob. Agents Chemother.* **21(1)**, 85-92 (1982).
2. Brogden, R. N., and Heel, R. C., *Drugs* **31(2)**, 96-130 (1986).
3. Bush, K., *Clin. Microbiol. Rev.* **1(1)**, 109-123 (1988).
4. Drawz, S.M. and Bonomo, R.A., *Clin. Microbiol. Rev.* **23(1)**, 160-201 (2010).
5. Coudron, P.E., Moland, E.S., and Sanders, C.C., *J. Clin. Microbiol.* **35(10)**, 2593-2597 (1997).

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, 11/11/2022

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