

PRODUCT INFORMATION



Moxalactam (sodium salt)

Item No. 15869

CAS Registry No.: 64953-12-4

Formal Name: (6R,7R)-7-[[[(2R)-2-carboxy-2-(4-hydroxyphenyl)acetyl]amino]-7-methoxy-3-[[[(1-methyl-1H-tetrazol-5-yl)thio]methyl]-8-oxo-5-oxa-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, disodium salt

Synonyms:

MF: $C_{20}H_{18}N_6O_9S \cdot 2Na$

FW: 564.4

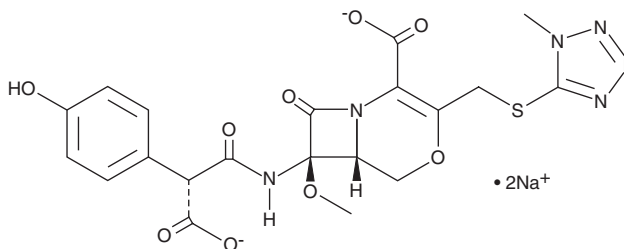
Purity: $\geq 85\%$ (mixture of isomers)

UV/Vis.: λ_{max} : 229, 271 nm

Supplied as: A crystalline solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Moxalactam (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the moxalactam (sodium salt) in the solvent of choice, which should be purged with an inert gas. Moxalactam (sodium salt) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of moxalactam (sodium salt) in DMSO and water is approximately 5 and 50 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Moxalactam is a β -lactam antibiotic in the cephalosporin class. It is classified as an oxacephem, as oxygen replaces sulfur in position one of the dihydrothiazolidine ring.¹ While moxalactam inhibits most commonly occurring gram positive, gram negative, and anaerobic bacteria, it is ineffective against methicillin-resistant *S. aureus* and *Enterococci*.¹ In certain situations, this antibiotic impairs normal hemostasis, most commonly manifested as a prolonged bleeding time.²

References

1. Neu, H.C. The in vitro activity, human pharmacology, and clinical effectiveness of new β -lactam antibiotics. *Annu. Rev. Pharmacol. Toxicol.* **22**, 599-642 (1982).
2. Sattler, F.R., Weitekamp, M.R., and Ballard, J.O. Potential for bleeding with the new beta-lactam antibiotics. *Ann. Intern. Med.* **105**(6), 924-931 (1986).

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