

PRODUCT INFORMATION



Ceftobiprole Medocaril (sodium salt)

Item No. 44248

CAS Registry No.: 252188-71-9

Formal Name: (6R,7R)-7-[[[(2Z)-2-(5-amino-1,2,4-thiadiazol-3-yl)-2-(hydroxyimino)acetyl]amino]-3-[(E)-[(3'R)-1'-[[[(5-methyl-2-oxo-1,3-dioxol-4-yl)methoxy]carbonyl]-2-oxo-[1,3'-bipyrrolidin]-3-ylidene)methyl]-8-oxo, 5-thia-1-azabicyclo[4.2.0]oct-2-ene-2-carboxylic acid, monosodium salt

Synonyms: BAL 5788, RO-5788

MF: $C_{26}H_{25}N_8O_{11}S_2 \cdot Na$

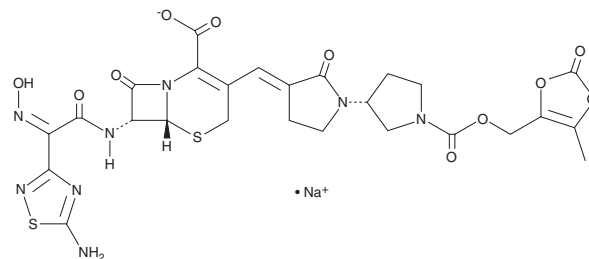
FW: 712.6

Purity: $\geq 90\%$

Supplied as: A 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

Ceftobiprole medocaril (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the ceftobiprole medocaril (sodium salt) in the solvent of choice, which should be purged with an inert gas.

Ceftobiprole medocaril (sodium salt) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Ceftobiprole medocaril is a water-soluble prodrug form of the cephalosporin antibiotic ceftobiprole.¹ *In vivo*, ceftobiprole medocaril (71 mg/kg) reduces the number of lung colony-forming units (CFUs) in mouse models of pneumonia induced by *H. influenzae*, *E. cloacae*, or *K. pneumoniae*. Formulations containing ceftobiprole medocaril have been used in the treatment of acute bacterial skin and skin structure infections and community-acquired bacterial pneumonia.

Reference

1. Rouse, M.S., Hein, M.M., Anguita-Alonso, P., *et al.* Ceftobiprole medocaril (BAL5788) treatment of experimental Haemophilus influenzae, Enterobacter cloacae, and Klebsiella pneumoniae murine pneumonia. *Diagn. Microbiol. Infect. Dis.* **55**(4), 333-336 (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, 10/17/2025

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