

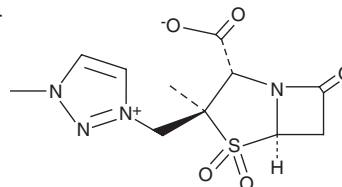
PRODUCT INFORMATION



Enmetazobactam

Item No. 30715

CAS Registry No.: 1001404-83-6
Formal Name: 3-[[[(2S,3S,5R)-2-carboxy-3-methyl-4,4-dioxido-7-oxo-4-thia-1-azabicyclo[3.2.0]hept-3-yl]methyl]-1-methyl-1H-1,2,3-triazolium, inner salt
Synonym: AAI-101
MF: C₁₁H₁₄N₄O₅S
FW: 314.3
Purity: ≥95%
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

Enmetazobactam is supplied as a solid. A stock solution may be made by dissolving the enmetazobactam in the solvent of choice, which should be purged with an inert gas. Enmetazobactam is soluble in DMSO.

Description

Enmetazobactam is a β -lactamase inhibitor.¹ It has greater activity against class A β -lactamases (IC_{50} s = 0.006-0.52 μ M) than class B or class D β -lactamases (IC_{50} s = 8.7->100 μ M). Enmetazobactam in combination with the cephalosporin antibiotic cefepime (Item No. 23633) is active against 223 *Enterobacteriaceae* clinical isolates, including multidrug-resistant *E. coli* and *K. pneumoniae* isolates, with an MIC_{50} value of 0.125 mg/L.² *In vivo*, enmetazobactam reduces mortality in mouse models of septicemia induced by *E. coli*, *K. pneumoniae*, or *E. cloacae* when used in combination with cefepime.¹

References

1. Papp-Wallace, K.M., Bethel, C.R., Caillon, J., *et al.* Beyond piperacillin-tazobactam: Cefepime and AAI101 as a potent β -lactam- β -lactamase inhibitor combination. *Antimicrob Agents Chemother.* **63(5)**, e00105-19 (2019).
2. Crandon, J.L. and Nicolau, D.P. *In vitro* activity of cefepime/AAI101 and comparators against cefepime non-susceptible *Enterobacteriaceae*. *Pathogens* **4(3)**, 620-625 (2015).

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, 11/03/2022

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