

# PRODUCT INFORMATION



## Tebipenem (hydrate)

Item No. 33875

**Formal Name:** (4R,5S,6S)-3-[[1-(4,5-dihydro-2-thiazolyl)-3-azetidiny]thio]-6-[(1R)-1-hydroxyethyl]-4-methyl-7-oxo-1-azabicyclo[3.2.0]hept-2-ene-2-carboxylic acid, hydrate

**Synonyms:** LJC 11,036, SPR859, TBPM

**MF:**  $C_{16}H_{21}N_3O_4S_2 \cdot XH_2O$

**FW:** 383.5

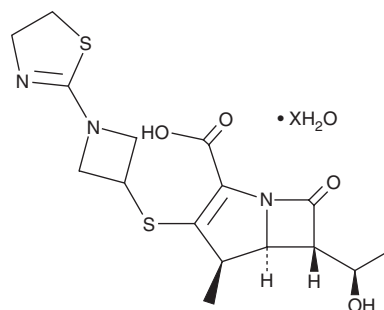
**Purity:**  $\geq 95\%$

**UV/Vis.:**  $\lambda_{max}$ : 210, 305 nm

**Supplied as:** A solid

**Storage:**  $-20^\circ C$

**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Tebipenem (hydrate) is supplied as a solid. A stock solution may be made by dissolving the tebipenem (hydrate) in the solvent of choice, which should be purged with an inert gas. Tebipenem (hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

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 tebipenem (hydrate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of tebipenem (hydrate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Tebipenem is a carbapenem antibiotic.<sup>1,2</sup> It is active against a panel of clinical isolates from a variety of bacterial species ( $MIC_{50s} = \leq 0.0039-8 \mu g/ml$ ), including methicillin-resistant strains of *S. aureus* and *S. epidermidis* and penicillin-resistant *S. pneumoniae*.<sup>1</sup> Tebipenem inhibits  $\beta$ -lactamase in a concentration-dependent manner.<sup>2</sup> It decreases the number of colony forming units (CFUs) in the lungs in a mouse model of penicillin-resistant *S. pneumoniae* infection when administered at doses ranging from 0.32 to 3.2 mg/kg.<sup>1</sup>

### References

1. Fujimoto, K., Takemoto, K., Hatano, K., *et al.* Novel carbapenem antibiotics for parenteral and oral applications: *In vitro* and *in vivo* activities of 2-aryl carbapenems and their pharmacokinetics in laboratory animals. *Antimicrob. Agents Chemother.* **57**(2), 697-707 (2013).
2. Hazra, S., Xu, H., and Blanchard, J.S. Tebipenem, a new carbapenem antibiotic, is a slow substrate that inhibits the  $\beta$ -lactamase from *Mycobacterium tuberculosis*. *Biochemistry* **53**(22), 3671-3678 (2014).

#### 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, 11/21/2022

#### 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](http://WWW.CAYMANCHEM.COM)