

PRODUCT INFORMATION



Sultamicillin

Item No. 46190

CAS Registry No.: 76497-13-7

Formal Name: (2S,5R,6R)-6-[[[(2R)-2-amino-2-phenylacetyl]amino]-3,3-dimethyl-7-oxo-4-thia-1-azabicyclo[3.2.0]heptane-2-carboxylic acid, [[[(2S,5R)-3,3-dimethyl-4,4-dioxido-7-oxo-4-thia-1-azabicyclo[3.2.0]hept-2-yl]carbonyl]oxy]methyl ester

Synonym:

CP 49,952

MF: $C_{25}H_{30}N_4O_9S_2$

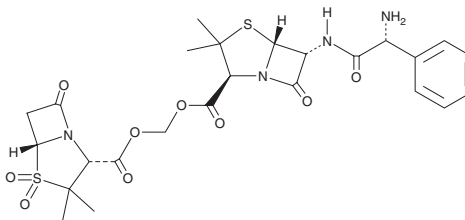
FW: 594.7

Purity: $\geq 98\%$

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

Sultamicillin is supplied as a solid. A stock solution may be made by dissolving the sultamicillin in the solvent of choice, which should be purged with an inert gas. Sultamicillin is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Sultamicillin is a prodrug form of ampicillin (Item Nos. 35740 | 14417) and sulbactam (Item No. 15024).¹ Formulations containing sultamicillin have been used in the treatment of bacterial infections.

Reference

1. English, A.R., Girard, D., and Haskell, S.L. Pharmacokinetics of dultamicillin in mice, rats, and dogs. *Antimicrob. Agents Chemother.* **25**(5), 599-602 (1984).

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, 08/19/2026

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