

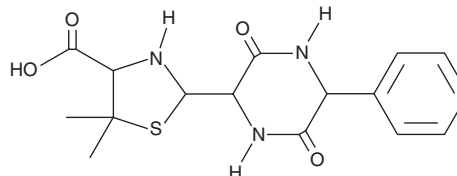
PRODUCT INFORMATION



Ampicillin Impurity C

Item No. 42900

CAS Registry No.: 49841-96-5
Formal Name: 2-(3,6-dioxo-5-phenyl-2-piperazinyl)-5,5-dimethyl-4-thiazolidinecarboxylic acid
Synonym: NSC 362657
MF: C₁₆H₁₉N₃O₄S
FW: 349.4
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

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

Description

Ampicillin impurity C is a metabolite of ampicillin (Item Nos. 14417 | 35740).^{1,2}

References

1. Suwanrumpha, S., Flory, D.A., Freas, R.B., *et al.* Tandem mass spectrometric studies of the fragmentation of penicillins and their metabolites. *Biomed. Environ. Mass Spectrom.* **16**(1-12), 381-386 (1988).
2. Suwanrumpha, S. and Freas, R.B. Identification of metabolites of ampicillin using liquid chromatography/thermospray mass spectrometry and fast atom bombardment tandem mass spectrometry. *Biomed. Environ. Mass Spectrom.* **18**(11), 983-994 (1989).

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, 06/17/2025

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