

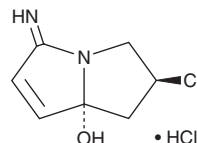
PRODUCT INFORMATION



Clazamycin A (hydrochloride)

Item No. 40224

CAS Registry No.: 71743-75-4
Formal Name: (2S-trans)-2-chloro-2,3-dihydro-5-imino-1H-pyrrolizin-7a(5H)-ol, monohydrochloride
Synonym: NSC 332188
MF: C₇H₉ClN₂O • HCl
FW: 209.1
Purity: ≥95%
Supplied as: A powder
Storage: -20°C (under argon atmosphere)
Stability: ≥1 year
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Clazamycin A (hydrochloride) is supplied as a powder. A stock solution may be made by dissolving the clazamycin A (hydrochloride) in the solvent of choice. Clazamycin A (hydrochloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas. Clazamycin A is slightly soluble in methanol. Clazamycin A (hydrochloride) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Clazamycin A is an antibiotic and a microbial metabolite originally isolated from *Streptomyces*.¹ It is active against a variety of bacteria, including *S. aureus*, *B. anthracis*, *B. subtilis*, *E. coli*, and *P. aeruginosa* (MICs = 100, 6.25, 100, 50, and 25 µg/ml, respectively).

Reference

1. Horiuchi, Y., Kondo, S., Ikeda, T., *et al.* New antibiotics clazamycins A and B. *J. Antibiot. (Tokyo)* **32(70)**, 762-764 (1979).

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, 11/08/2023

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