

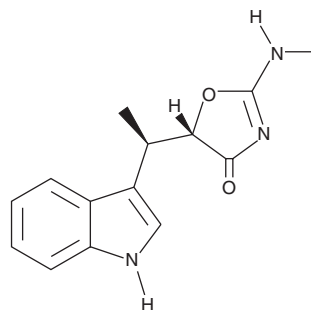
PRODUCT INFORMATION



Indolmycin

Item No. 15385

CAS Registry No.: 21200-24-8
Formal Name: 5S-[(1R)-1-(1H-indol-3-yl)ethyl]-2-(methylamino)-4(5H)-oxazolone
Synonyms: PA 155A, TAK-083
MF: C₁₄H₁₅N₃O₂
FW: 257.3
Purity: ≥95%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Indolmycin is supplied as a powder. A stock solution may be made by dissolving the indolmycin in the solvent of choice, which should be purged with an inert gas. Indolmycin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Indolmycin is an antibiotic isolated from various *Streptomyces* strains that functions by competitively inhibiting the prokaryotic tryptophanyl-tRNA synthetase (IC₅₀ = 30 μM) without significantly affecting the corresponding eukaryotic synthetase.^{1,2} It exhibits antibacterial activity against mupirocin- and fusidic acid-resistant strains of *Staphylococci*.³

References

1. Werner, R.G., Thorpe, L.F., Reuter, W., *et al.* Indolmycin inhibits prokaryotic tryptophanyl-tRNA ligase. *Eur. J. Biochem.* **68**(1), 1-3 (1976).
2. Bogosian, G., Haydock, P.V., and Somerville, R.L. Indolmycin-mediated inhibition and stimulation of transcription at the trp promoter of *Escherichia coli*. *J. Bacteriol.* **153**(2), 1120-1123 (1983).
3. Hurdle, J.G., O'Neill, A.J., and Chopra, I. Anti-staphylococcal activity of indolmycin, a potential topical agent for control of staphylococcal infections. *J. Antimicrob. Chemother.* **52**(4), 549-552 (2004).

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, 09/28/2022

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