

PRODUCT INFORMATION



Indanomycin

Item No. 21761

CAS Registry No.: 66513-28-8

Formal Name: (α R,2R,5S,6R)-6-[(1E,3E)-1-ethyl-4-[[1S,3aR,4S,5S,7aS)-1-ethyl-2,3,3a,4,5,7a-hexahydro-4-(1H-pyrrol-2-ylcarbonyl)-1H-inden-5-yl]-1,3-butadien-1-yl]tetrahydro- α ,5-dimethyl-2H-pyran-2-acetic acid

Synonym: X 14547A

MF: $C_{31}H_{43}NO_4$

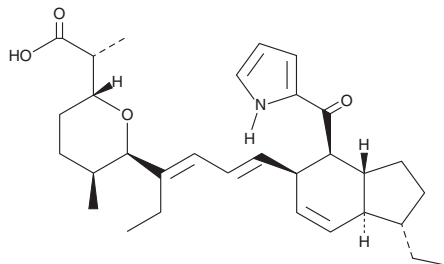
FW: 493.7

Purity: $\geq 95\%$

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

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

Description

Indanomycin is an antibiotic of the pyrrolaketoidane class, which includes an unusual indane ring structure.¹ It has bactericidal activity against Gram-positive bacteria, with minimum inhibitory concentration (MIC) values of $\leq 0.2 \mu\text{g/ml}$.²

References

1. Roege, K.E. and Kelly, W.L. Biosynthetic origins of the ionophore antibiotic indanomycin. *Org. Lett.* **11**(2), 397-300 (2009).
2. Liu, C.-M., Hermann, T.E., Liu, M., *et al.* X-14547A, a new ionophorous antibiotic produced by *Streptomyces antibioticus* NRRL 8167. Discovery, fermentation, biological properties and taxonomy of the producing culture. *J. Antibiot. (Tokyo)* **32**(2), 95-99 (1978).

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, 08/24/2023

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