

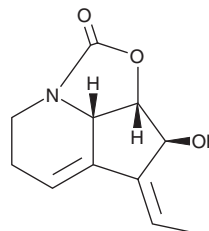
PRODUCT INFORMATION



Streptazolin

Item No. 27266

CAS Registry No.: 80152-07-4
Formal Name: 4-ethylidene-2aS,3S,4Z,6,7,7bS-hexahydro-3-hydroxy-1H-2-oxa-7a-azacyclopent[cd]inden-1-one
Synonym: (+)-Streptazolin
MF: C₁₁H₁₃NO₃
FW: 207.2
Purity: ≥70%
Supplied as: A solid
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

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

Description

Streptazolin is a fungal metabolite originally isolated from *S. viridochromogenes*.¹ It increases NF-κB activity when used at concentrations ranging from 60 to 130 μg/ml, at least in part, via PI3K signaling.² Streptazolin enhances TNF-α secretion induced by LPS and enhances IL-8 secretion when used alone or in combination with LPS in THP-1 Blue cells.

References

1. Drautz, H., Zaehner, H., Kupfer, E., *et al.* Metabolic products of microorganisms. 205. Isolation and structure of streptazolin. *Helv. Chim. Acta.* **64(6)**, 1752-1765 (1981).
2. Perry, J.A., Koteva, K., Verschoor, C.P., *et al.* A macrophage-stimulating compound from a screen of microbial natural products. *J. Antibiot. (Tokyo)* **68(1)**, 40-46 (2015).

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, 12/21/2022

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