

PRODUCT INFORMATION



Satratoxin H

Item No. 30199

CAS Registry No.: 53126-64-0
Formal Name: (2'S,4E,9R,10E,12Z,16R,16aS,18R,19aR,23aR,25R)-6,7,16,16a,19a,22-hexahydro-25-hydroxy-9-[(1S)-1-hydroxyethyl]-16a,21-dimethyl-spiro[5,9:16,18-dimethano-1H,3H,23H-[1,6,12]trioxacyclooctadecino[3,4-d][1]benzopyran-17(18H),2'-oxirane]-3,14(9H)-dione

MF: C₂₉H₃₆O₉

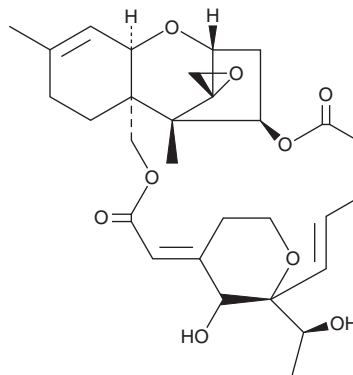
FW: 528.6

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Satratoxin H is supplied as a crystalline solid. A stock solution may be made by dissolving the satratoxin H in the solvent of choice, which should be purged with an inert gas. Satratoxin H is soluble in ethanol, DMSO, and dichloromethane.

Description

Satratoxin H is a trichothecene mycotoxin that has been found in *Stachybotrys*.¹ It induces production of reactive oxygen species (ROS), increases in malondialdehyde (MDA) levels, and induces apoptosis in PC12 cells when used at concentrations ranging from 5 to 100 nM. Satratoxin H is cytotoxic to human umbilical vein endothelial cells (HUVECs; IC₅₀ = 6.8 ng/ml), as well as HepG2, A549, A204, U937, and Jurkat cancer cells (IC₅₀s = 1.2-3.4 ng/ml).² *In vivo*, satratoxin H induces ulceration of the small intestine in, and is lethal to mice (LD₅₀ = 5.69 mg/kg).³

References

1. Nusuetrong, P., Pengsuparp, T., Meksuriyen, D., *et al.* Satratoxin H generates reactive oxygen species and lipid peroxides in PC12 cells. *Biol. Pharm. Bull.* **31(6)**, 1115-1120 (2008).
2. Nielsen, C., Casteel, M., Didier, A., *et al.* Trichothecene-induced cytotoxicity on human cell lines. *Mycotoxin Res.* **25(2)**, 77-84 (2009).
3. Yoshizawa, T., Ohtsubo, K., Sasaki, T., *et al.* Acute toxicities of satratoxins G and H in mice – a histopathological observation with special reference to the liver injury caused by satratoxin G. *Proc. Jpn. Assoc. Mycotoxicol.* **23**, 53-57 (1986).

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, 02/07/2020

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