

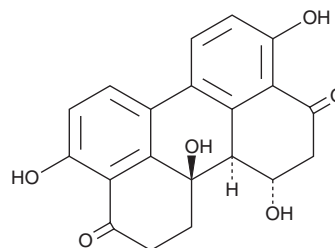
PRODUCT INFORMATION



Altertoxin I

Item No. 21129

CAS Registry No.: 56258-32-3
Formal Name: (1S,12aR,12bS)-1,2,11,12,12a,12b-hexahydro-1,4,9,12a-tetrahydroxy-3,10-perylene-1,2-dione
Synonym: Dihydroalterperyleneol
MF: C₂₀H₁₆O₆
FW: 352.3
Purity: ≥95%
Supplied as: A powder
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Altertoxin I is supplied as a powder. A stock solution may be made by dissolving the altertoxin I in the solvent of choice, which should be purged with an inert gas. Altertoxin I is soluble (≥10 mg/ml) in ethanol and dichloromethane.

Description

Altertoxin I is a natural mycotoxin first isolated from fungi of the genus *Alternaria*. It is a hydroxyperylenequinone that shows mutagenic activity against *S. typhimurium* and inhibits human topoisomerase II and bacterial gyrase.^{1,2} Altertoxin I completely inhibits replication of HIV-1 virus at 2.2 μM in a cell-based assay.³

References

1. Jarolim, K., Del Favero, G., Ellmer, D., *et al.* Dual effectiveness of *Alternaria* but not *Fusarium* mycotoxins against human topoisomerase II and bacterial gyrase. *Arch. Toxicol.* **91**(4), 2007-2016 (2016).
2. Stack, M.E. and Prival, M.J. Mutagenicity of the *Alternaria* metabolites altertoxins I, II, and III. *Appl. Environ. Microbiol.* **52**(4), 718-722 (1986).
3. Bashyal, B.P., Wellensiek, B.P., Ramakrishnan, R., *et al.* Altertoxins with potent anti-HIV activity from *Alternaria tenuissima* QUE1Se, a fungal endophyte of *Quercus emoryi*. *Bioorg. Med. Chem.* **22**(21), 6112-6116 (2014).

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, 03/20/2024

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