

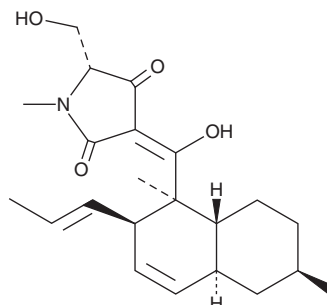
PRODUCT INFORMATION



Epiequisetin

Item No. 25454

CAS Registry No.: 255377-45-8
Formal Name: 5R-(hydroxymethyl)-3E-[hydroxy[(1S,2R,4aS,6R,8aR)-1,2,4a,5,6,7,8,8a-octahydro-1,6-dimethyl-2-(1E)-1-propen-1-yl-1-naphthalenyl)methylene]-1-methyl-2,4-pyrrolidinedione
MF: C₂₂H₃₁NO₄
FW: 373.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungi/*Fusarium equiseti*



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

Laboratory Procedures

Epiequisetin is supplied as a solid. A stock solution may be made by dissolving the epiequisetin in the solvent of choice, which should be purged with an inert gas. Epiequisetin is soluble in methanol.

Description

Epiequisetin is a fungal metabolite and an epimer of equisetin (Item No. 18196).¹ It inhibits HIV-1 integrase 3' end-processing and strand transfer activities. Epiequisetin is phytotoxic and inhibits the germination of various seeds and growth of young seedlings.²

Reference

1. Singh, S.B., Zink, D.L., Goetz, M.A., *et al.* Equisetin and a novel opposite stereochemical homolog phomasetin, two fungal metabolites as inhibitors of HIV-1 integrase. *Tetrahedron Lett.* **39(16)**, 2243-2246 (1998).
2. Wheeler, M.H., Stipanovic, R.D., and Puckhaber, L.S. Phytotoxicity of equisetin and epi-equisetin isolated from *Fusarium equiseti* and *F. pallidoroseum*. *Mycol. Res.* **103(8)**, 967-973 (1999).

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, 01/23/2024

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