

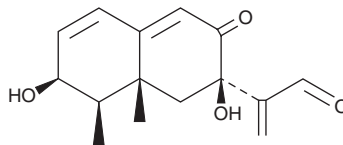
PRODUCT INFORMATION



Dihydrobipolaroxin

Item No. 40916

CAS Registry No.: 100665-30-3
Formal Name: 1,2R,3,7S,8R,8aR-hexahydro-2,7-dihydroxy-8,8a-dimethyl- α -methylene-3-oxo-2-naphthaleneacetaldehyde
MF: $C_{15}H_{18}O_4$
FW: 262.3
Purity: $\geq 70\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Curvularia lunata*



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

Laboratory Procedures

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

Description

Dihydrobipolaroxin is a sesquiterpene that has been found in *B. cynodontis*.¹

Reference

1. Sugawara, F., Strobel, G., Fisher, L.E., et al. Bipolaroxin, a selective phytotoxin produced by *Bipolaris cynodontis*. *Proc. Nat. Acad. Sci. USA* **82**(24), 8291-8294 (1985).

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, 05/20/2024

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