

PRODUCT INFORMATION



Norstictic Acid

Item No. 29581

CAS Registry No.: 571-67-5
Formal Name: 1,3-dihydro-1,4,10-trihydroxy-5,8-dimethyl-3,7-dioxo-7H-isobenzofuro[4,5-b][1,4]benzodioxepin-11-carboxaldehyde

MF: C₁₈H₁₂O₉

FW: 372.3

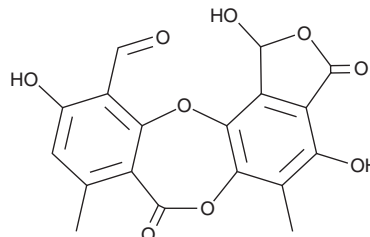
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Lobaria* sp.



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

Laboratory Procedures

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

Description

Norstictic acid is a lichen metabolite that has been found in *R. farinacea* and has antimicrobial and anticancer activities.^{1,2} It is active against *A. hydrophilia*, *B. subtilis*, *L. monocytogenes*, *P. vulgaris*, *S. faecalis*, *C. albicans*, and *C. glabrata* (MICs = 0.1-6.7 mM).¹ Norstictic acid (1-100 μM) inhibits proliferation, migration, and invasion of MDA-MB-231 cells.² *In vivo*, norstictic acid (15 mg/kg, i.p.) reduces tumor growth in an MDA-MB-231/GFP mouse xenograft model.

References

1. Tay, T., Türk, A.Ö., Yilmaz, M., *et al.* Evaluation of the antimicrobial activity of the acetone extract of the lichen *Ramalina farinacea* and its (+)-usnic acid, norstictic acid, and protocetraric acid constituents. *Z. Naturforsch. C J. Biosci.* **59(5-6)**, 384-388 (2004).
2. Ebrahim, H.Y., Elsayed, H.E., Mohyeldin, M.M., *et al.* Norstictic acid inhibits breast cancer cell proliferation, migration, invasion, and *in vivo* invasive growth through targeting c-Met. *Phytother. Res.* **30(4)**, 557-566 (2016).

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, 08/15/2023

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