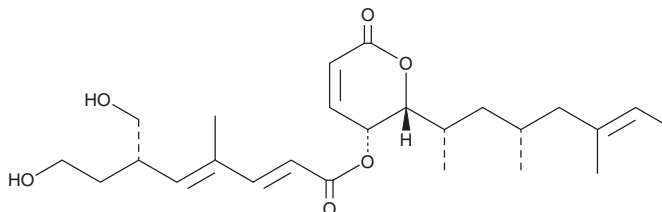


PRODUCT INFORMATION



(-)-Rasfonin Item No. 27581

CAS Registry No.: 303156-68-5
Formal Name: (2E,4E,6S)-8-hydroxy-6-(hydroxymethyl)-4-methyl-2,4-octadienoic acid, (2R,3R)-3,6-dihydro-6-oxo-2-[(1S,3R,5E)-1,3,5-trimethyl-5-hepten-1-yl]-2H-pyran-3-yl ester
Synonym: TT-1
MF: C₂₅H₃₈O₆
FW: 434.6
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years
Item Origin: Fungus/*Trichurus* sp.



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

Laboratory Procedures

(-)-Rasfonin is supplied as a liquid. A stock solution may be made by dissolving the (-)-rasfonin in the solvent of choice. (-)-Rasfonin is soluble in organic solvents such as DMSO, methanol, and ethanol, which should be purged with an inert gas. The solubility of (-)-rasfonin in DMSO is approximately 1 mg/ml.

Description

(-)-Rasfonin is a fungal metabolite that has been found in *T. terrophilus*.¹ It inhibits proliferation of mouse splenocytes induced by concanavalin A (Item No. 14951) and LPS (IC₅₀s = 0.7 and 0.5 µg/ml, respectively).

Reference

1. Fujimoto, H., Okamoto, Y., Sone, E., *et al.* Eleven new 2-pyrones from a fungi imperfecti, *Trichurus terrophilus*, found in a screening study guided by immunomodulatory activity. *Chem. Pharm. Bull. (Tokyo)* 53(8), 923-929 (2005).

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, 08/09/2023

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