

PRODUCT INFORMATION

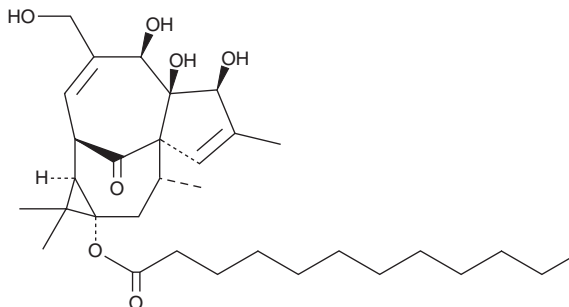


Dodecanoic Acid ingenol ester

Item No. 34213

CAS Registry No.: 54706-70-6
Formal Name: dodecanoic acid,
(1aR,2S,5R,5aR,6S,8aS,9R,10aS)-
1,1a,2,5,5a,6,9,10-octahydro-5,5a,6-
trihydroxy-4-(hydroxymethyl)-1,1,7,9-
tetramethyl-11-oxo-10aH-2,8a-
methanocyclopenta[a]cyclopropa[e]
cyclodecen-10a-yl ester
Synonyms: 13-O-Dodecanoylingenol,
13-Oxyingenol Dodecanoate

MF: C₃₂H₅₀O₇
FW: 546.7
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Dodecanoic acid ingenol ester is supplied as a solid. A stock solution may be made by dissolving the dodecanoic acid ingenol ester in the solvent of choice, which should be purged with an inert gas. Dodecanoic acid ingenol ester is slightly soluble in chloroform and methanol.

Description

Dodecanoic acid ingenol ester is an ingenane diterpene that has been found in *E. kansui* and has antiviral and antinematodal activities.¹ It inhibits replication of HIV-1 in infected MT-4 cells (EC₅₀ = 33.7 nM).² Dodecanoic acid ingenol ester also has antinematodal activity against the plant pathogenic nematode *B. xylophilus* in a *B. cinerea*-based bioassay with a minimum effective dose (MED) of 2.5 µg/cotton ball.¹

References

1. Shi, J., Li, Z., Nitoda, T., *et al.* Antinematodal activities of ingenane diterpenes from *Euphorbia kansui* and their derivatives against the pine wood nematode (*Bursaphelenchus xylophilus*). *Z. Naturforsch. C. J. Biosci.* **63(1-2)**, 59-65 (2008).
2. Liu, Q., Li, W., Huang, L., *et al.* Identification, structural modification, and dichotomous effects on human immunodeficiency virus type 1 (HIV-1) replication of ingenane esters from *Euphorbia kansui*. *Eur. J. Med. Chem.* **156**, 618-627 (2018).

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, 11/18/2022

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