

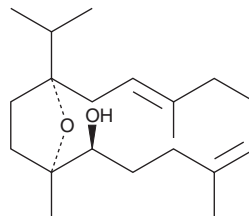
PRODUCT INFORMATION



Incensole

Item No. 22895

CAS Registry No.: 22419-74-5
Formal Name: (1R,2S,5E,9E,12S)-1,5,9-trimethyl-12-(1-methylethyl)-15-oxabicyclo[10.2.1]pentadeca-5,9-dien-2-ol
MF: $C_{20}H_{34}O_2$
FW: 306.5
Purity: $\geq 98\%$
Supplied as: A crystalline 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

Incensole is supplied as a crystalline solid. A stock solution may be made by dissolving the incensole in the solvent of choice, which should be purged with an inert gas. Incensole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of incensole in these solvents is approximately 50 mg/ml.

Incensole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, incensole should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Incensole has a solubility of approximately 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Incensole is a diterpene originally isolated from frankincense produced by *B. carterii*.¹ It inhibits degradation of inhibitor of NF- κ B (I κ B α) in a concentration-dependent manner in TNF- α -stimulated HeLa cells when used at concentrations ranging from 60 to 140 μ M.

Reference

1. Moussaieff, A.S., E., Kashman, Y., Fride, E., *et al.* Incensole acetate, a novel anti-inflammatory compound isolated from *Boswellia* resin, inhibits nuclear factor- κ B activation. *Mol. Pharmacol.* **72**(6), 1657-1664 (2007).

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, 11/03/2022

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