

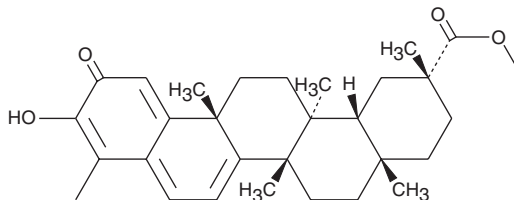
PRODUCT INFORMATION



Pristimerin

Item No. 13621

CAS Registry No.: 1258-84-0
Formal Name: 3-hydroxy-9 β ,13 α -dimethyl-2-oxo-24,25,26-trinoroleana-1(10),3,5,7-tetraen-29-oic acid, methyl ester
Synonyms: Celastrol methyl ester, NSC 99281
MF: C₃₀H₄₀O₄
FW: 464.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 203, 419 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Pristimerin is supplied as a crystalline solid. A stock solution may be made by dissolving the pristimerin in the solvent of choice, which should be purged with an inert gas. Pristimerin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pristimerin in ethanol is approximately 3 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Monoacylglycerol lipase (MAGL) hydrolyzes the endogenous cannabinoid 2-arachidonoyl glycerol (2-AG), terminating its capacity to activate cannabinoid receptors. Pristimerin is a naturally occurring terpenoid that potently inhibits MAGL (IC₅₀ = 93 nM).¹ Its actions are rapid, reversible, and noncompetitive.¹ Pristimerin (1 μ M) significantly increases 2-AG levels in isolated rat neurons, indicating that it inhibits endogenous MAGL in cultured cells.¹ Moreover, it does not increase levels of palmitoyl ethanolamide, suggesting that pristimerin does not affect the activity of fatty acid amide hydrolase (FAAH).

Reference

1. King, A.R., Dotsey, E.Y., Lodola, A., *et al.* Discovery of potent and reversible monoacylglycerol lipase inhibitors. *Chem. Biol.* **16**(10), 1045-1052 (2009).

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, 10/19/2022

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