

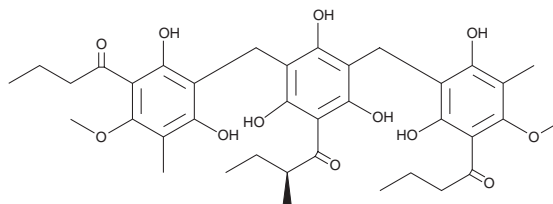
PRODUCT INFORMATION



Agrimol B

Item No. 34996

CAS Registry No.: 55576-66-4
Formal Name: (2S)-1-[3,5-bis[[2,6-dihydroxy-4-methoxy-3-methyl-5-(1-oxobutyl)phenyl]methyl]-2,4,6-trihydroxyphenyl]-2-methyl-1-butanone
MF: $C_{37}H_{46}O_{12}$
FW: 682.8
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 230, 288 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Agrimonia pilosa*



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

Laboratory Procedures

Agrimol B is supplied as a solid. A stock solution may be made by dissolving the agrimol B in the solvent of choice, which should be purged with an inert gas. Agrimol B is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Agrimol B is a polyphenol that has been found in *A. pilosa* with adipogenesis inhibitory and anticancer activities.^{1,2} It induces nuclear translocation of sirtuin 1 (SIRT1) and reduces expression of peroxisome proliferator-activated receptor γ (PPAR γ), C/EBP α , and fatty acid synthase (FASN) in 3T3-L1 adipocytes when used at a concentration of 10 μM . Agrimol B reduces lipid droplet accumulation in 3T3-L1 adipocytes ($\text{IC}_{50} = 3.35 \mu\text{M}$). It is cytotoxic to PC3 prostate and A549 lung cancer cells *in vitro* and reduces tumor volume in a PC3 mouse xenograft model when administered at a dose of 10 mg/kg.²

References

1. Wang, S., Zhang, Q., Zhang, Y., *et al.* Agrimol B suppresses adipogenesis through modulation of SIRT1-PPAR gamma signal pathway. *Biochem. Biophys. Res. Commun.* **477**(3), 454-460 (2016).
2. Hnit, S.S.T., Ding, R., Bi, L., *et al.* Agrimol B present in *Agrimonia pilosa* Ledeb impedes cell cycle progression of cancer cells through G₀ state arrest. *Biomed. Pharmacother.* **141**, 111795 (2021).

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, 11/23/2022

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