

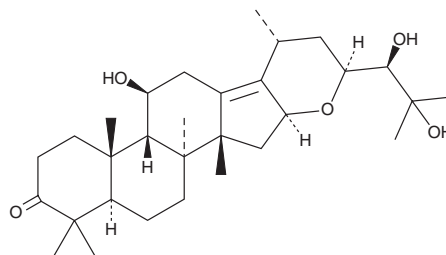
PRODUCT INFORMATION



Alisol F

Item No. 35627

CAS Registry No.: 155521-45-2
Formal Name: (8 α ,9 β ,11 β ,14 β ,16 β ,23S,24R)-16,23-epoxy-11,24,25-trihydroxy-dammar-13(17)-en-3-one
MF: C₃₀H₄₈O₅
FW: 488.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Alisma plantago-aquatica* tubers



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

Laboratory Procedures

Alisol F is supplied as a solid. A stock solution may be made by dissolving the alisol F in the solvent of choice, which should be purged with an inert gas. Alisol F is soluble in chloroform, dichloromethane, ethyl acetate, acetone and DMSO.

Description

Alisol F is a triterpene that has been found in *A. orientale* and has anti-inflammatory and hepatoprotective activities.^{1,2} It reduces levels of inducible nitric oxide synthase (iNOS) and COX-2 and the production of TNF- α , IL-6, and IL-1 β in LPS-stimulated RAW 264.7 cells.² Alisol F (20 mg/kg) reduces plasma alanine transaminase (ALT) and aspartate aminotransferase (AST) levels, hepatic fibrosis, and hepatic levels of phosphorylated ERK in a mouse model of LPS/D-galactosamine-induced acute liver injury.^{1,2}

References

1. Choi, E., Jang, E., and Lee, J.-H. Pharmacological activities of *Alisma orientale* against nonalcoholic fatty liver disease and metabolic syndrome: Literature review. *Evid. Based Complement. Alternat. Med.* 2943162 (2019).
2. Bi, X., Wang, P., Ma, Q., et al. Anti-inflammatory activities and liver protection of alisol F and 25-anhydroalisol F through the inhibition of MAPK, STAT3, and NF- κ B activation in vitro and in vivo. *Molecules* **22**(6), 951 (2017).

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, 12/12/2022

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