

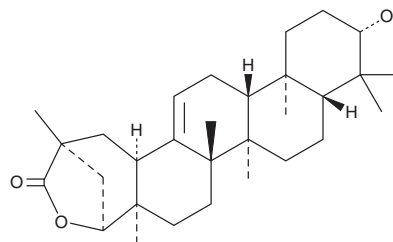
PRODUCT INFORMATION



Wilforlide A

Item No. 34654

CAS Registry No.: 84104-71-2
Formal Name: (3 β ,20 α ,22 α)-3,22-dihydroxy-olean-12-en-29-oic acid, γ -lactone
Synonyms: (-)-Abruslactone, Abruslactone A, Regelide
MF: C₃₀H₄₆O₃
FW: 454.7
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Tripterygium wilfordii*



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

Laboratory Procedures

Wilforlide A is supplied as a solid. A stock solution may be made by dissolving the wilforlide A in the solvent of choice, which should be purged with an inert gas. Wilforlide A is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

Wilforlide A is a triterpene that has been found in *T. wilfordii* and has anti-inflammatory activity.^{1,2} It inhibits xylene-induced ear swelling in mice when administered at doses of 60 or 300 $\mu\text{g/kg}$.¹ Wilforlide A in combination with triptolide reduces inflammation, bone damage, and synovial hyperplasia in the inflamed joints in a rat model of collagen-induced arthritis.²

References

1. Xue, M., Jiang, Z.-Z., Liu, J.-P., *et al.* Comparative study on the anti-inflammatory and immune suppressive effect of Wilforlide A. *Fitoterapia* **81(8)**, 1109-1112 (2010).
2. Zhu, X., Zhang, J., Huo, R., *et al.* Evaluation of the efficacy and safety of different Tripterygium preparations on collagen-induced arthritis in rats. *J. Ethnopharmacol.* **158(Pt A)**, 283-290 (2014).

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, 11/22/2022

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