

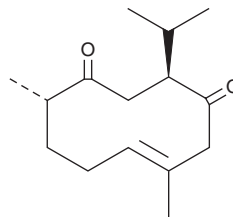
PRODUCT INFORMATION



Curdione

Item No. 30365

CAS Registry No.: 13657-68-6
Formal Name: (3S,6E,10S)-6,10-dimethyl-3-(1-methylethyl)-6-cyclodecene-1,4-dione
Synonym: (+)-Curdione
MF: C₁₅H₂₄O₂
FW: 236.3
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Curcuma zedoaria*



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

Laboratory Procedures

Curdione is supplied as a crystalline solid. A stock solution may be made by dissolving the curdione in the solvent of choice, which should be purged with an inert gas. Curdione is soluble in ethanol, DMSO, and dimethyl formamide. The solubility of curdione in these solvents is approximately 30 mg/ml.

Description

Curdione is a sesquiterpene that has been found in *C. wenyujin* and has diverse biological activities. It reduces the cytopathic effect of influenza A in MDCK cells (IC₅₀ = 28.32 μM).¹ Curdione (100 μM) inhibits thrombin-induced aggregation, AMP-activated protein kinase (AMPK) and integrin αIIbβ3 phosphorylation, and increases in talin and vinculin levels in washed platelets.² It inhibits prostaglandin E₂ (PGE₂; Item No. 14010) production in LPS-stimulated RAW 264.7 macrophages (IC₅₀ = 1.12 μM).³

References

1. Dong, J.-Y., Ma, X.-Y., Cai, X.-Q., *et al.* Sesquiterpenoids from *Curcuma wenyujin* with anti-influenza viral activities. *Phytochemistry* **85**, 122-128 (2013).
2. Hui, F., Beibei, G., Yingli, Z., *et al.* Curdione inhibits thrombin-induced platelet aggregation via regulating the AMP-activated protein kinase-vinculin/talin-integrin αIIbβ3 sign pathway. *Phytomedicine* **61**, 152859 (2019).
3. Oh, O.-J., Min, H.-Y., and Lee, S.K. Inhibition of inducible prostaglandin E₂ production and cyclooxygenase-2 expression by curdione from *Curcuma zedoaria* *Arch. Pharm. Res.* **30(10)**, 1236-1239 (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, 05/07/2020

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