

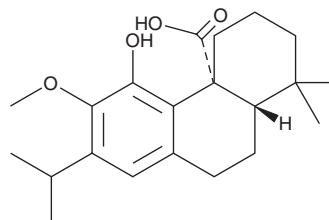
PRODUCT INFORMATION



12-O-Methylcarnosic Acid

Item No. 36175

CAS Registry No.: 62201-71-2
Formal Name: (4aR,10aS)-1,3,4,9,10,10a-hexahydro-5-hydroxy-6-methoxy-1,1-dimethyl-7-(1-methylethyl)-4a(2H)-phenanthrenecarboxylic acid
Synonyms: Carnosic acid 12-methyl ether, 12-Methoxycarnosic Acid
MF: C₂₁H₃₀O₄
FW: 346.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Rosmarinus officinalis*



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

Laboratory Procedures

12-O-Methylcarnosic acid is supplied as a solid. A stock solution may be made by dissolving the 12-O-methylcarnosic acid in the solvent of choice, which should be purged with an inert gas. 12-O-Methylcarnosic acid is soluble in chloroform.

Description

12-O-Methylcarnosic acid is a diterpene that has been found in *S. repens* and has antileishmanial and gastroprotective activities.^{1,2} It induces mortality in *L. donovani* amastigotes (IC₅₀ = 0.75 μM).¹ 12-O-Methylcarnosic acid (10 mg/kg) reduces lesion area in a mouse model of hydrochloric acid and ethanol-induced gastric ulcers.²

References

1. Mokoka, T.A., Peter, X.K., Fouche, G., *et al.* Antileishmanial activity of 12-methoxycarnosic acid from *Salvia repens* burch. ex benth. (lamiaceae). *S. Afr. J. Bot.* **90**, 93-95 (2014).
2. Theodulz, C., Pertino, M.W., Rodríguez, J.A., *et al.* Gastroprotective effect and cytotoxicity of carnosic acid derivatives. *Planta Med.* **77**(9), 882-887 (2011).

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, 12/19/2022

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