

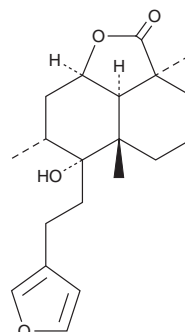
PRODUCT INFORMATION



Marrubiin

Item No. 33876

CAS Registry No.: 465-92-9
Formal Name: (2aS,5aS,6R,7R,8aR,8bR)-6-[2-(3-furanyl)ethyl]decahydro-6-hydroxy-2a,5a,7-trimethyl-2H-naphtho[1,8-bc]furan-2-one
Synonyms: (+)-Marrubiin, NSC 36693
MF: C₂₀H₂₈O₄
FW: 332.4
Purity: ≥98%
UV/Vis.: λ_{max}: 213 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Marrubium vulgare*



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

Laboratory Procedures

Marrubiin is supplied as a solid. A stock solution may be made by dissolving the marrubiin in the solvent of choice, which should be purged with an inert gas. Marrubiin is soluble in ethanol and chloroform.

Description

Marrubiin is a labdane diterpenoid lactone originally isolated from *M. vulgare* that has diverse biological activities.¹⁻³ It inhibits potassium chloride-induced contractions of rat aortic rings (IC₅₀ = 24 μM).¹ Marrubiin (100 mg/kg) inhibits ovalbumin-induced allergic ear edema in mice.² It inhibits acetic acid-induced writhing and capsaicin-induced paw licking in mice with ID₅₀ values of 2.2 and 28.8 μmol/kg, respectively.³

References

1. El Bardai, S., Morel, N., Wibo, M., *et al.* The vasorelaxant activity of marrubenol and marrubiin from *Marrubium vulgare*. *Planta Med.* **69(1)**, 75-77 (2003).
2. Stulzer, H.K., Tagliari, M.P., Zampirolo, J.A., *et al.* Antioedematogenic effect of marrubiin obtained from *Marrubium vulgare*. *J. Ethnopharmacol.* **108(3)**, 379-384 (2006).
3. De Jesus, R.A., Cechinel-Filho, V., Oliveira, A.E., *et al.* Analysis of the antinociceptive properties of marrubiin isolated from *Marrubium vulgare*. *Phytomedicine* **7(2)**, 111-115 (2000).

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, 11/28/2022

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