

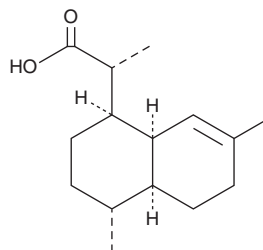
PRODUCT INFORMATION



Dihydroartemisinin Acid

Item No. 25142

CAS Registry No.: 85031-59-0
Formal Name: (α R,1R,4R,4aS,8aS)-1,2,3,4,4a,5,6,8a-octahydro- α ,4,7-trimethyl-1-naphthaleneacetic acid
MF: C₁₅H₂₄O₂
FW: 236.4
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Dihydroartemisinin acid is supplied as a crystalline solid. A stock solution may be made by dissolving the dihydroartemisinin acid in the solvent of choice, which should be purged with an inert gas. Dihydroartemisinin acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of dihydroartemisinin acid in these solvents is approximately 16, 10, and 20 mg/ml, respectively.

Dihydroartemisinin acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dihydroartemisinin acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Dihydroartemisinin acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dihydroartemisinin acid is a biosynthetic precursor to the antimalarial agent artemisinin (Item No. 11816) and an intermediate in the synthesis of artemisinin from artemisinic acid (Item No. 25059).^{1,2}

References

1. Tian, N., Li, J., Liu, S., *et al.* Simultaneous isolation of artemisinin and its precursors from *Artemisia annua* L. by preparative RP-HPLC. *Biomed. Chromatogr.* **26(6)**, 708-713 (2012).
2. Roth, R.J. and Acton, N. A simple conversion of artemisinic acid into artemisinin. *J. Nat. Prod.* **52(5)**, 1183-1185 (1989).

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, 12/01/2022

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