

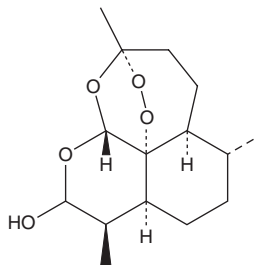
PRODUCT INFORMATION



Dihydroartemisinin (mixed isomers)

Item No. 41908

CAS Registry No.: 131175-87-6
Formal Name: (3R,5aS,6R,8aS,9R,12R,12aR)-decahydro-3,6,9-trimethyl-3,12-epoxy-12H-pyrano[4,3-j]-1,2-benzodioxepin-10-ol
MF: C₁₅H₂₄O₅
FW: 284.4
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Artemisia annua*



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

Laboratory Procedures

Dihydroartemisinin (mixed isomers) is supplied as a crystalline solid. A stock solution may be made by dissolving the dihydroartemisinin (mixed isomers) in the solvent of choice, which should be purged with an inert gas. Dihydroartemisinin (mixed isomers) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile and ethanol.

Description

Dihydroartemisinin (mixed isomers) is a mixture of isomers of dihydroartemisinin, an active metabolite of artemisinin (Item No. 11816).¹ This product contains α-dihydroartemisinin and β-dihydroartemisinin.

Reference

1. Dai, X., Zhang, X., Chen, W., et al. Dihydroartemisinin: A potential natural anticancer drug. *Int. J. Biol. Sci.* **17**(2), 603-622 (2021).

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, 11/08/2024

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