

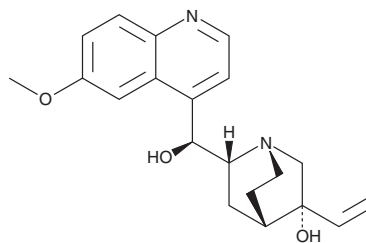
PRODUCT INFORMATION



(3R)-hydroxy Quinidine

Item No. 23410

CAS Registry No.: 60761-51-5
Formal Name: 6'-methoxy-cinchonan-3 α ,9S-diol
Synonym: 3-*epi*-hydroxy Quinidine
MF: C₂₀H₂₄N₂O₃
FW: 340.4
Purity: \geq 98%
Supplied as: A solid
Storage: 4°C
Stability: \geq 4 years



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

Laboratory Procedures

(3R)-hydroxy Quinidine is supplied as a solid. A stock solution may be made by dissolving the (3R)-hydroxy quinidine in the solvent of choice, which should be purged with an inert gas. (3R)-hydroxy Quinidine is soluble in methanol.

Description

(3R)-hydroxy Quinidine is an isomer of (3S)-hydroxy quinidine (Item No. 23411), the major metabolite of quinidine.¹

Reference

1. Leroyer, R., Varoquax, O., Guillebastres, R., *et al.* Study on protein binding of 3-(R)- and 3-(S)-hydroxy derivatives of quinidine and dihydroquinidine. *Therapie* **41(3)**, 159-162 (1986).

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, 12/16/2022

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