

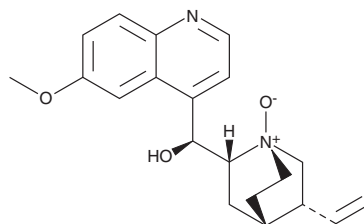
PRODUCT INFORMATION



Quinidine N-oxide

Item No. 22381

CAS Registry No.: 70116-00-6
Formal Name: 6'-methoxy-cinchonan-9S-ol 1-oxide
MF: $C_{20}H_{24}N_2O_3$
FW: 340.4
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Quinidine N-oxide is supplied as a solid. A stock solution may be made by dissolving the quinidine N-oxide in the solvent of choice, which should be purged with an inert gas. Quinidine N-oxide is soluble in DMSO.

Description

Quinidine N-oxide is a pharmacologically inactive quinidine metabolite of quinidine (Item No. 20356).¹ Quinidine, an antiarrhythmic agent, undergoes rapid first-pass metabolism by the cytochrome P450 isoforms CYP3A4, CYP2C9, and CYP2E1, with CYP3A4 being the most active enzyme in quinidine N-oxide formation.²

References

1. Ha, H.-R., Vozeh, S., Uematsu, T., *et al.* Kinetics and dynamics of quinidine-N-oxide in healthy subjects. *Clin. Pharmacol. Ther.* **42**(3), 341-345 (1987).
2. Nielsen, T.L., Rasmussen, B.B., Flinois, J.-P., *et al.* In vitro metabolism of quinidine: the (3S)-3-hydroxylation of quinidine is a specific marker reaction for cytochrome P-4503A4 activity in human liver microsomes. *J. Pharmacol. Exp. Ther.* **289**(1), 31-37 (1999).

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, 11/19/2024

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