

PRODUCT INFORMATION



1'-hydroxy Midazolam

Item No. 10385

CAS Registry No.: 59468-90-5
Formal Name: 8-chloro-6-(2-fluorophenyl)-4H-imidazo[1,5-a][1,4]benzodiazepine-1-methanol

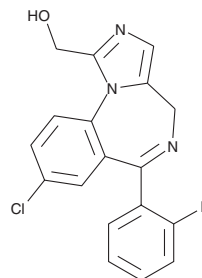
Synonym: Ro 21-6347
MF: C₁₈H₁₃ClFN₃O
FW: 341.8

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Description

1'-hydroxy Midazolam is the major metabolite of the anesthetic, midazolam. It is produced rapidly by the actions of hepatic cytochrome P450 3A and has been shown to be equipotent to midazolam.¹

Reference

1. Perloff, M.D., Von Moltke, L.L., Court, M.H., *et al.* Midazolam and triazolam biotransformation in mouse and human liver microsomes: Relative contribution of CYP3A and CYP2C isoforms. *J. Pharmacol. Exp. Ther.* **292**(2), 618-628 (2016).

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, 09/01/2016

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