

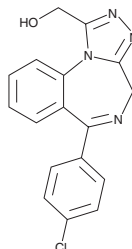
PRODUCT INFORMATION



α -hydroxy-4'-chloro Deschloroalprazolam

Item No. 40819

Formal Name: (6-(4-chlorophenyl)-4H-benzo[f][1,2,4]triazolo[4,3-a][1,4]diazepin-1-yl)methanol
MF: C₁₇H₁₃ClN₄O
FW: 324.8
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Description

α -hydroxy-4'-chloro Deschloroalprazolam (Item No. 40819) is an analytical reference standard categorized as a benzodiazepine. α -hydroxy-4'-chloro Deschloroalprazolam is a potential metabolite of 4'-chloro deschloroalprazolam (Item No. 36057) based on the published metabolism of triazolam (Item Nos. 39683 | 22577).¹ At the time α -hydroxy-4'-chloro deschloroalprazolam (Item No. 40819) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

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 (2000).

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, 04/18/2024

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