

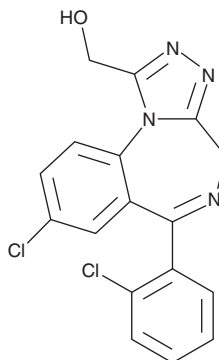
PRODUCT INFORMATION



α -hydroxy Triazolam

Item No. 39390

CAS Registry No.: 37115-45-0
Formal Name: 8-chloro-6-(2-chlorophenyl)-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine-1-methanol
MF: C₁₇H₁₂Cl₂N₄O
FW: 359.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 222 nm
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.

Description

α -hydroxy Triazolam (Item No. 39390) is an analytical reference standard categorized as a benzodiazepine. α -hydroxy Triazolam is an active metabolite of triazolam (Item Nos. 22577 | 39683).^{1,2} It has anticonvulsant and sedative effects in mice.² This product is intended for research and forensic applications.

References

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).
2. Hester, J.B., Jr. and Von Voigtlander, P. 6-Aryl-4H-s-triazolo[4,3-a][1,4]benzodiazepines. Influence of 1-substitution on pharmacological activity. *J. Med. Chem.* **22**(11), 1390-1398 (1979).

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, 01/13/2026

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