

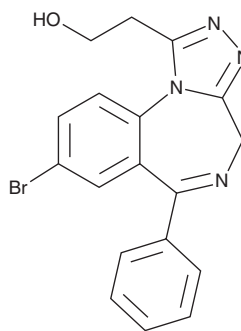
PRODUCT INFORMATION



α -hydroxy Ethylbromazolam

Item No. 45237

Formal Name: 2-(8-bromo-6-phenyl-4H-benzo[f][1,2,4]triazolo[4,3-a][1,4]diazepin-1-yl)ethan-1-ol
MF: C₁₈H₁₅BrN₄O
FW: 383.2
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 Ethylbromazolam (Item No. 45237) is an analytical reference standard categorized as a benzodiazepine. α -hydroxy Ethylbromazolam is a potential metabolite of ethylbromazolam (Item No. 42723) similar to the published metabolism of flubromazolam (Item Nos. 26352 | 16435).¹ At the time α -hydroxy ethylbromazolam (Item No. 45237) 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. El Balkhi, S., Chaslot, M., Picard, N., *et al.* Characterization and identification of eight designer benzodiazepine metabolites by incubation with human liver microsomes and analysis by a triple quadrupole mass spectrometer. *Int. J. Legal Med.* **131**(4), 979-988 (2017).

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, 03/04/2026

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