

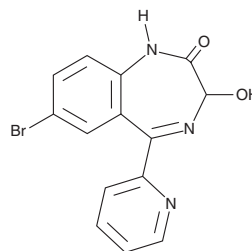
PRODUCT INFORMATION



3-hydroxy Bromazepam

Item No. 44155

CAS Registry No.: 13132-73-5
Formal Name: 7-bromo-1,3-dihydro-3-hydroxy-5-(2-pyridinyl)-2H-1,4-benzodiazepin-2-one
MF: C₁₄H₁₀BrN₃O₂
FW: 332.2
Purity: ≥98%
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

3-hydroxy Bromazepam (Item No. 44155) is an analytical reference standard categorized as a benzodiazepine metabolite.¹ 3-hydroxy Bromazepam is an active metabolite of bromazepam (Item No. 14260).² This product is intended for research and forensic applications.

References

1. Chèze, M., Villain, M., and Pépin, G. Determination of bromazepam, clonazepam and metabolites after a single intake in urine and hair by LC-MS/MS. Application to forensic cases of drug facilitated crimes. *Forensic Sci. Int.* **145(2-3)**, 123-130 (2004).
2. Altamura, A.C., Moliterno, D., Paletta, S., *et al.* Understanding the pharmacokinetics of anxiolytic drugs. *Expert Opin. Drug Metab Toxicol.* **9(4)**, 423-440 (2013).

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, 01/07/2026

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