

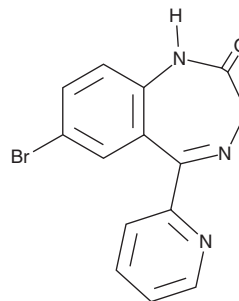
PRODUCT INFORMATION



Bromazepam

Item No. 14260

CAS Registry No.: 1812-30-2
Formal Name: 7-bromo-1,3-dihydro-5-(2-pyridinyl)-2H-1,4-benzodiazepin-2-one
Synonyms: Compendium, Lexotanil, NSC 140669, Ro 5-3350
MF: C₁₄H₁₀BrN₃O
FW: 316.2
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 322 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Bromazepam is a psychoactive benzodiazepine derivative and prescription drug that is commonly abused and often used in 'drug-facilitated crimes'.^{1,2} This product is intended for forensic and research applications.

References

1. Yakushiji, T., Fukuda, T., Oyama, Y., *et al.* Effects of benzodiazepines and non-benzodiazepine compounds on the GABA-induced response in frog isolated sensory neurones. *Br. J. Pharmacol.* **98**(3), 735-740 (1989).
2. Lendoiro, E., Quintela, O., de Castro, A., *et al.* Target screening and confirmation of 35 licit and illicit drugs and metabolites in hair by LC-MSMS. *Forensic Sci. Int.* **217**(1-3), 207-215 (2012).

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, 07/14/2023

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