

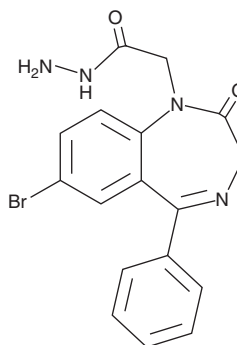
PRODUCT INFORMATION



Gidazepam

Item No. 36421

CAS Registry No.: 129186-29-4
Formal Name: 7-bromo-2,3-dihydro-2-oxo-5-phenyl-1H-1,4-benzodiazepine-1-acetic acid, hydrazide
Synonyms: Hidazepam, Hydazepam
MF: C₁₇H₁₅BrN₄O₂
FW: 387.2
Purity: ≥95%
UV/Vis.: λ_{max}: 229 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Description

Gidazepam (Item No. 36421) is an analytical reference standard categorized as a benzodiazepine. Gidazepam has been abused recreationally and has anxiolytic properties in mice.¹ Gidazepam partially substitutes for phenazepam (Item No. 9000849) in a two-lever drug discrimination test in rats.² This product is intended for research and forensic applications.

References

1. Maskell, P.D., Wilson, G., and Manchester, K.R. Designer benzodiazepines gidazepam and desalkygidazepam (bromonordiazepam): What do we know? *J. Anal. Toxicol.* **47(6)**, e53 (2023).
2. Kalinina, T.S., Garibova, T.L., and Voronina, T.A. Discriminative effects of phenazepam and gidazepam in rats: Comparison with other GABA-related drugs. *Pharmacol. Biochem. Behav.* **64(2)**, 397-401 (1999).

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, 10/17/2023

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