

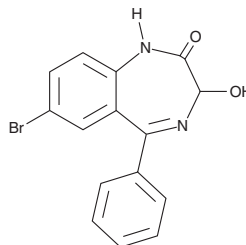
PRODUCT INFORMATION



3-hydroxy Desalkylgidazepam

Item No. 42258

CAS Registry No.: 37891-18-2
Formal Name: 7-bromo-1,3-dihydro-3-hydroxy-5-phenyl-2H-1,4-benzodiazepin-2-one
Synonyms: 3-hydroxy Bromonordiazepam
MF: $C_{15}H_{11}BrN_2O_2$
FW: 331.2
Purity: $\geq 98\%$
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 Desalkylgidazepam (Item No. 42258) is an analytical reference standard categorized as a benzodiazepine.¹ 3-hydroxy Desalkylgidazepam is an active metabolite of gidazepam (Item No. 36421) and desalkylgidazepam (Item Nos. 40621 | 36608).^{1,2} It has been used as a precursor in the synthesis of substituted benzodiazepinones.³ This product is intended for research and forensic applications.

References

1. Maskell, P.D., Wilson, G., and Manchester, K.R. Designer benzodiazepines gidazepam and desalkylgidazepam (bromonordiazepam): What do we know? *J. Anal. Toxicol.* **47**(6), e53 (2023).
2. Kolyvanov, G.B., Zherdev, V.P., Chirkov, A.M., et al. [Gidazepam biotransformation and pharmacokinetics in different species of animals and man]. *Eksp. Klin. Farmakol.* **56**(3), 48-50 (1993).
3. Pavlovskii, V.I., Ushakov, I.Y., Kabanova, T.A., et al. Synthesis and analgesic activity of 3-arylamino-1,2-dihydro-3H-1,4-benzodiazepin-2-ones. *Pharm. Chem. J.* **19**(9), 592-597 (2015).

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, 10/28/2024

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