

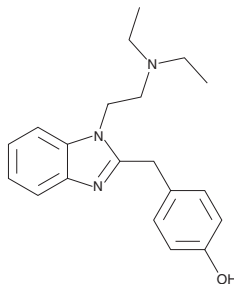
PRODUCT INFORMATION



4'-hydroxy Desnitazene

Item No. 39321

CAS Registry No.: 110151-18-3
Formal Name: 4-[[1-[2-(diethylamino)ethyl]-1H-benzimidazol-2-yl]methyl]-phenol
Synonym: O-desethyl Etodesnitazene
MF: C₂₀H₂₅N₃O
FW: 323.4
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

4'-hydroxy Desnitazene (Item No. 39321) is an analytical reference standard categorized as an opioid metabolite.^{1,2} 4'-hydroxy Desnitazene is a metabolite of etodesnitazene (Item Nos. 29916 | 39892). This product is intended for research and forensic applications.

References

1. Grigoryev, A. Kavanagh, P. Dowling, G., *et al.* Tentative identification of etazene (etodesnitazene) metabolites in rat serum and urine by gas chromatography-mass spectrometry and accurate mass liquid chromatography-mass spectrometry. *J. Anal. Toxicol.* **46(9)**, 1032-1037 (2023)
2. Chaturvedi, K., Hewamanna, I., Pandey, P., *et al.* Identification of the putative binding site of a benzimidazole opioid (etazene) and its metabolites at μ -opioid receptor: A human liver microsomal assay and systematic computational study. *Molecules* **28(4)**, 1601 (2023).

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, 04/07/2025

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