

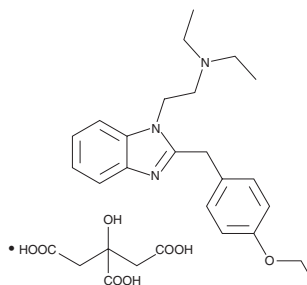
PRODUCT INFORMATION



Etodesnitazene (citrate)

Item No. 29916

Formal Name: 2-[(4-ethoxyphenyl)methyl]-N,N-diethyl-1H-benzimidazole-1-ethanamine, 2-hydroxy-1,2,3-propanetricarboxylate
Synonyms: Desnitroetonitazene, Etazene, Etazone
MF: C₂₂H₂₉N₃O • C₆H₈O₇
FW: 543.6
Purity: ≥98%
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

Etodesnitazene (citrate) (Item No. 29916) is an analytical reference standard categorized as an opioid.¹ Etodesnitazene has been found in parcels seized by law enforcement.² Etodesnitazene is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

- Vandeputte, M.M., Van Uytvanghe, K., Layle, N.K., *et al.* Synthesis, chemical characterization, and μ -opioid receptor activity assessment of the emerging group of “nitazene” 2-benzylbenzimidazole synthetic opioids. *ACS Chem. Neurosci.* **12(7)**, 1241-1251 (2021).
- Siczek, M., Zawadzki, M., Siczek, M., *et al.* Etazene (N,N diethyl 2-[(4-ethoxyphenyl)methyl] 1H benzimidazol 1 yl) ethan 1 amine (dihydrochloride): A novel benzimidazole opioid NPS identified in seized material: Crystal structure and spectroscopic characterization. *Forensic Toxicol.* **39(1)**, 146-155 (2021).

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, 07/07/2025

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