

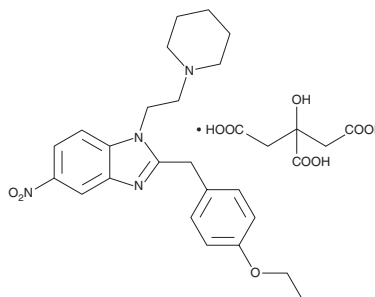
PRODUCT INFORMATION



N-Piperidinyl Etonitazene (citrate) (exempt preparation)

Item No. 45947

Formal Name: 2-[(4-ethoxyphenyl)methyl]-5-nitro-1-[2-(1-piperidinyl)ethyl]-1H-benzimidazole, 2-hydroxy-1,2,3-propanetricarboxylate
Synonyms: Etonitazepipne, N-Piperidyl Etonitazene
MF: $C_{23}H_{28}N_4O_3 \cdot C_6H_8O_7$
FW: 600.6
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 241 nm
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥ 3 years



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

Description

N-Piperidinyl etonitazene (citrate) (exempt preparation) (Item No. 45947) is an analytical reference standard categorized as an opioid.¹ N-Piperidinyl etonitazene induces antinociceptive effects in mice in the hot plate test. N-Piperidinyl etonitazene is regulated as a Schedule I compound in the United States. N-Piperidinyl etonitazene (citrate) (exempt preparation) (Item No. 45947) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Vandeputte, M.M., Verougstraete, N., Walther, D., *et al.* First identification, chemical analysis and pharmacological characterization of N-piperidinyl etonitazene (etonitazepipne), a recent addition to the 2-benzylbenzimidazole opioid subclass. *Arch. Toxicol.* **96**(6), 1865-1880 (2022).

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, 05/12/2026

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