

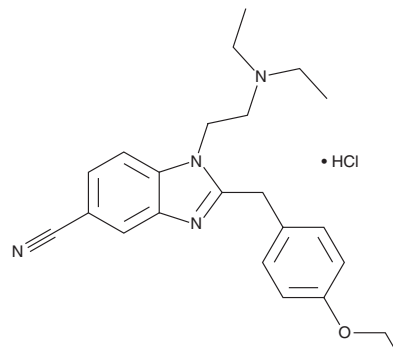
PRODUCT INFORMATION



5-cyano Etodesnitazene (hydrochloride)

Item No. 43960

CAS Registry No.: 15419-87-1
Formal Name: 1-[2-(diethylamino)ethyl]-2-[[4-ethoxyphenyl)methyl]-1H-benzimidazole-5-carbonitrile, monohydrochloride
Synonym: Etocyanazene
MF: C₂₃H₂₈N₄O • HCl
FW: 413.0
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

5-cyano Etodesnitazene (hydrochloride) (Item No. 43960) is an analytical reference standard categorized as an opioid.^{1,2} This product is intended for research and forensic applications.

References

1. Lecolier, S. and Trouiller, G. New benzimidazoles with morphine-like activity. *Laboratories of the Powder Directorate, 12, Quai Henri IV, Paris (4th) 2*, 16-24 (1966).
2. Ujváry, I., Christie, R., Evans-Brown, M., et al. DARK classics in chemical neuroscience: Etonitazene and related benzimidazoles. *ACS Chem. Neurosci.* **12**(7), 1072-1092 (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, 10/10/2025

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