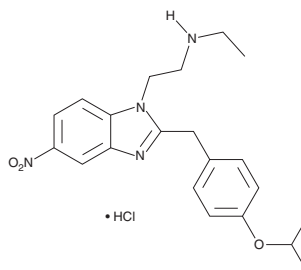


PRODUCT INFORMATION

N-desethyl Isotonitazene (hydrochloride)

Item No. 30216

Formal Name: N-ethyl-2-[[4-(1-methylethoxy)phenyl]methyl]-5-nitro-1H-benzimidazole-1-ethanamine, monohydrochloride
MF: C₂₁H₂₆N₄O₃ • HCl
FW: 418.9
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 301 nm
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

N-desethyl Isotonitazene (hydrochloride) (Item No. 30216) is an analytical reference standard categorized as an opioid.^{1,2} N-desethyl Isotonitazene is a metabolite of isotonitazene (Item Nos. 30880 | 27255). N-desethyl Isotonitazene is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Krotulski, A.J., Papsun, D.M., Kacinko, S.L., *et al.* Isotonitazene quantitation and metabolite discovery in authentic forensic casework. *J. Anal. Toxicol.* **44**(6), 521-530 (2020).
2. Vandeputte, M.M., Van Uytanghe, 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).

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, 07/29/2024

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