

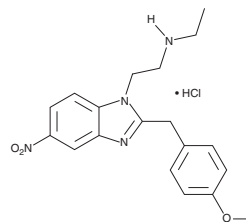
PRODUCT INFORMATION



N-desethyl Metonitazene (hydrochloride)

Item No. 38178

Formal Name: N-ethyl-2-(2-(4-methoxybenzyl)-5-nitro-1H-benzo[d]imidazol-1-yl)ethan-1-amine, monohydrochloride
MF: C₁₉H₂₂N₄O₃ • HCl
FW: 390.9
Purity: ≥98%
UV/Vis.: λ_{max}: 240 nm
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

N-desethyl Metonitazene (Item No. 38178) is an analytical reference standard categorized as an opioid metabolite.¹ N-desethyl Metonitazene is a metabolite of metonitazene (Item No. 26398). This product is intended for research and forensic applications.

Reference

1. Krotulski, A.J., Papsun, D.M., Walton, S.E., *et al.* Metonitazene in the United States-Forensic toxicology assessment of a potent new synthetic opioid using liquid chromatography mass spectrometry. *Drug Test Anal.* **13(10)**, 1697-1711 (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, 07/12/2023

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