

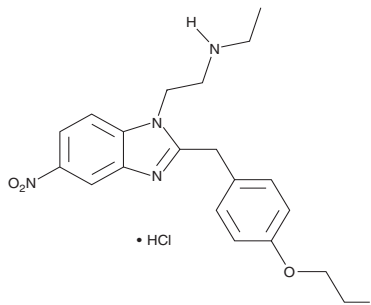
# PRODUCT INFORMATION



## N-desethyl Fluetonitazene (hydrochloride)

Item No. 41296

**Formal Name:** N-ethyl-2-(2-(4-(2-fluoroethoxy)benzyl)-5-nitro-1H-benzo[d]imidazol-1-yl)ethan-1-amine  
**Synonyms:** DFNZ, N-desethyl Fluornitazene  
**MF:** C<sub>20</sub>H<sub>23</sub>FN<sub>4</sub>O<sub>3</sub> • HCl  
**FW:** 422.9  
**Purity:** ≥98%  
**Supplied as:** A 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 Fluetonitazene (DFNZ) (hydrochloride) (Item No. 41296) is an analytical reference standard categorized as an opioid.<sup>1</sup> DFNZ is an active metabolite of fluetonitazene (FNZ; Item No. 40834) that acts as a highly potent  $\mu$ -opioid receptor agonist. It produces analgesia in rodent models of acute and chronic pain without inducing respiratory depression, tolerance, or  $\mu$ -opioid receptor downregulation after repeated dosing. DFNZ also has weaker reinforcing effects and increases dopamine neuron activity more slowly and to a lesser extent than FNZ in rodents. This product is intended for research and forensic applications.

### Reference

1. Gomez, J. L., Ventriglia, E. N., Frangos, Z. J., *et al.* A  $\mu$ -opioid receptor superagonist analgesic with minimal adverse effects. *Nature* **652(8112)**, 1393-1404 (2026).

#### 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, 06/15/2026

#### 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