

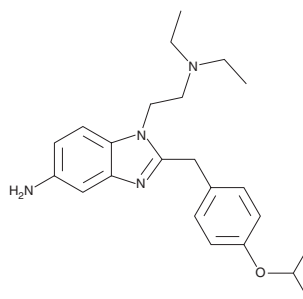
# PRODUCT INFORMATION



## 5-Aminoisotonitazene

Item No. 29318

**CAS Registry No.:** 2732926-25-7  
**Formal Name:** 5-amino-N,N-diethyl-2-[[4-(1-methylethoxy)phenyl]methyl]-1H-benzimidazole-1-ethanamine  
**MF:** C<sub>23</sub>H<sub>32</sub>N<sub>4</sub>O  
**FW:** 380.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221 nm  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

5-Aminoisotonitazene (Item No. 29318) is an analytical reference standard categorized as an opioid.<sup>1,2</sup> It is a metabolite of isotonitazene (Item Nos. 30880 | 27255). This product is intended for research and forensic applications.

### Reference

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 Uytfanghe, K., Layle, N.K., *et al.* Synthesis, chemical characterization, and  $\mu$ -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, 07/26/2023

#### 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