

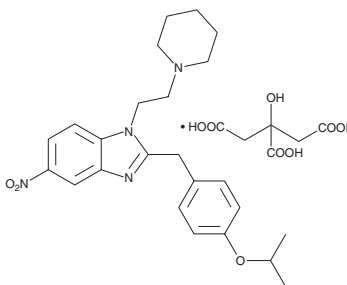
PRODUCT INFORMATION



N-Piperidinyl Isotonitazene (citrate)

Item No. 35106

Formal Name: 2-(4-isopropoxybenzyl)-5-nitro-1-(2-(piperidin-1-yl)ethyl)-1H-benzo[d]imidazole, monocation
MF: $C_{24}H_{30}N_4O_3 \cdot C_6H_8O_7$
FW: 614.7
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 239 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-Piperidinyl isotonitazene (citrate) (Item No. 34909) is an analytical reference standard that is categorized as an opioid.¹ This product is intended for research and forensic applications.

Reference

1. Kozell, L.B., Eshleman, A.J., Wolfrum, K.M., *et al.* Pharmacology of newly identified nitazene variants reveals structural determinants of affinity, potency, selectivity for mu opioid receptors. *Neuropharmacology* **276**, 110512 (2025).

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, 05/12/2026

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