

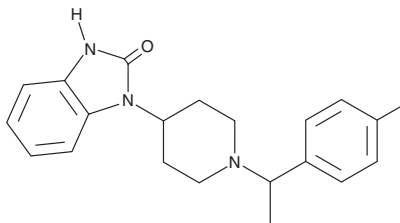
PRODUCT INFORMATION



Iodorphine

Item No. 37062

Formal Name: 1-(1-(1-(4-iodophenyl)ethyl)piperidin-4-yl)-1,3-dihydro-2H-benzo[d]imidazol-2-one
MF: C₂₀H₂₂IN₃O
FW: 447.3
Purity: ≥98%
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

Iodorphine (Item No. 37062) is an analytical reference standard categorized as an opioid.¹ Iodorphine induces antinociception and respiratory depression in mice, which can be reversed by naloxone (Item Nos. ISO60191 | 15594). This product is intended for research and forensic applications.

Reference

1. Vandeputte, M.M., Bilel, S., Tirri, M., *et al.* Elucidating the harm potential of brorphine analogues as new synthetic opioids: Synthesis, *in vitro*, and *in vivo* characterization. *Neuropharmacology* **260**, 110113 (2024).

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, 12/04/2024

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