

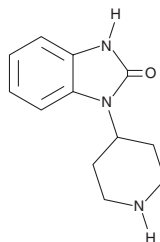
PRODUCT INFORMATION



nor-Orphine

Item No. 37339

CAS Registry No.: 20662-53-7
Formal Name: 1,3-dihydro-1-(4-piperidiny)-2H-benzimidazol-2-one
Synonyms: Orphine N-desalkyl metabolite, N-desalkyl Orphine
MF: C₁₂H₁₅N₃O
FW: 217.3
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

nor-Orphine (Item No. 37339) is an analytical reference standard categorized as an opioid metabolite.¹ nor-Orphine is a metabolite of buprenorphine (Item Nos. 34184 | 27816) and a potential metabolite of various other orphines based on the published metabolism of buprenorphine.^{1,2} It is also a precursor in the synthesis of various orphines. This product is intended for research and forensic applications.

References

1. Krotulski, A.J., Papsun, D.M., Noble, C., *et al.* Buprenorphine-investigation and quantitation of a new potent synthetic opioid in forensic toxicology casework using liquid chromatography-mass spectrometry. *J. Forensic Sci.* **66**(2), 664-676 (2021).
2. Vandeputte, M.M., Bilel, S., Tirri, M., *et al.* Elucidating the harm potential of buprenorphine 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, 01/31/2025

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