

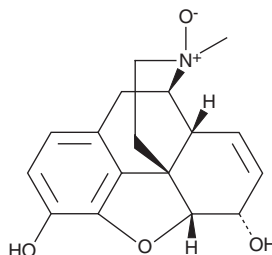
PRODUCT INFORMATION



Morphine N-oxide

Item No. 44701

CAS Registry No.: 639-46-3
Formal Name: 7,8-didehydro-4,5 α -epoxy-17-methyl-morphinan-3,6 α -diol, 17-oxide
Synonyms: Genomorphine, Morphine Oxide
MF: C₁₇H₁₉NO₄
FW: 301.3
Purity: $\geq 98\%$
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

Morphine N-oxide (Item No. 44701) is an analytical reference standard categorized as an opioid.¹ Morphine N-oxide is a degradation product and an active metabolite of morphine (Item Nos. ISO60147 | 15464 | 25955) that induces analgesia in mice and rats to a lesser extent than morphine.^{1,2} This product is intended for research and forensic applications.

References

1. Fennessy, M.R. The analgesic action of morphine-N-oxide. *Br. J. Pharmacol.* **34**(2), 337-344 (1968).
2. Yeh, S.-Y. and Lach, J.L. Stability of morphine in aqueous solution. II. Separation of morphine from its degradation products. *J. Pharm. Sci.* **50**, 30-34 (1961).

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, 06/29/2026

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