

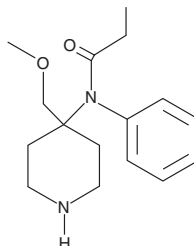
PRODUCT INFORMATION



Norsufentanil

Item No. 9001142

CAS Registry No.: 61086-18-8
Formal Name: N-[4-(methoxymethyl)-4-piperidiny]-N-phenyl-propanamide
Synonym: R 30451
MF: C₁₆H₂₄N₂O₂
FW: 276.4
Purity: ≥95%
Supplied as: A 10 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



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

Description

Sufentanil is a powerful synthetic opioid analgesic that is used in anesthesia and palliative care.¹⁻³ Norsufentanil is a metabolite of sufentanil, produced by oxidative N-dealkylation in the liver by cytochrome P450 isoforms.⁴ The physiological and toxicological properties of this compound are unknown. This product is intended for forensic applications.

References

1. Bovill, J.G., Sebel, P.S., and Stanley, T.H. Opioid analgesics in anesthesia: With special reference to their use in cardiovascular anesthesia. *Anesthesiology* **61**(6), 731-755 (1984).
2. Hodgson, P.S., Neal, J.M., Pollock, J.E., *et al.* The neurotoxicity of drugs given intrathecally (spinal). *Anesth. Analg.* **88**(4), 797-809 (1999).
3. King, S., Forbes, K., Hanks, G.W., *et al.* A systematic review of the use of opioid medication for those with moderate to severe cancer pain and renal impairment: A European Palliative Care Research Collaborative opioid guidelines project. *Palliat. Med.* **25**(5), 525-552 (2011).
4. Tateishi, T., Krivoruk, Y., Ueng, Y.F., *et al.* Identification of human liver cytochrome P-450 3A4 as the enzyme responsible for fentanyl and sufentanil N-dealkylation. *Anesth. Analg.* **82**(1), 167-172 (1996).

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, 10/20/2025

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