

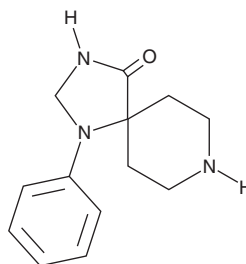
PRODUCT INFORMATION



nor-Spiroorphine

Item No. 45135

CAS Registry No.: 1021-25-6
Formal Name: 1-phenyl-1,3,8-triazaspiro[4.5]
decan-4-one
Synonym: NSC 96918
MF: C₁₃H₁₇N₃O
FW: 231.3
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-Spiroorphine (Item No. 45135) is an analytical reference standard categorized as a precursor in the synthesis of synthetic opioids, including spirochlorphine and spirobrorphine (Item No. 42301). nor-Spiroorphine is also a potential metabolite of spirochlorphine and spirobrorphine based on the published metabolism of brorphine (Item No. 27816).¹ At the time nor-spiroorphine (Item No. 45135) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

Reference

1. Krotulski, A.J., Papsun, D.M., Noble, C., *et al.* Brorphine-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).

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, 02/24/2026

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