

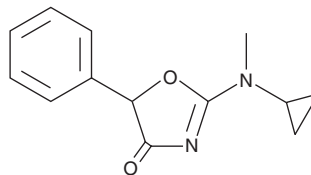
PRODUCT INFORMATION



N-methyl Cyclazodone

Item No. 27947

CAS Registry No.: 14461-92-8
Formal Name: 2-(cyclopropylmethylamino)-5-phenyl-4(5H)-oxazolone
Synonym: LD 4202
MF: C₁₃H₁₄N₂O₂
FW: 230.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

N-methyl Cyclazodone is supplied as a solid. A stock solution may be made by dissolving the N-methyl cyclazodone in the solvent of choice, which should be purged with an inert gas. N-methyl Cyclazodone is soluble in acetonitrile, methanol, and DMSO.

Description

N-methyl Cyclazodone (Item No. 27947) is an analytical reference standard that is structurally similar to known stimulants.¹ N-methyl Cyclazodone has been abused recreationally. This product is intended for research and forensic applications.

References

1. Holborn, T.J., Page, R., Schifano, F., *et al.* Self-medication with novel psychoactive substances (NPS): A systematic review. *Int. J. Ment. Health Ad.* (2023).

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, 01/30/2024

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