

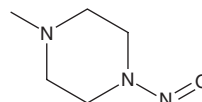
PRODUCT INFORMATION



N-Methyl-N'-nitrosopiperazine

Item No. 40300

CAS Registry No.: 16339-07-4
Formal Name: 1-methyl-4-nitroso-piperazine
Synonyms: MeNP, 1-Methyl-4-nitrosopiperazine, MNNG, MNPaz, NSC 523886
MF: C₅H₁₁N₃O
FW: 129.2
Purity: ≥95%
Supplied as: A neat liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

N-Methyl-N'-nitrosopiperazine is supplied as a neat liquid. A stock solution may be made by dissolving the N-methyl-N'-nitrosopiperazine in the solvent of choice, which should be purged with an inert gas. N-Methyl-N'-nitrosopiperazine is soluble in chloroform.

Description

N-Methyl-N'-nitrosopiperazine is a potential impurity in various pharmaceutical preparations, including sartans, non-steroidal anti-inflammatory drugs (NSAIDs), and thiazide diuretics.^{1,2}

References

1. Kalauz, A., Tiringier, K.V., Horváth, V., *et al.* Simultaneous determination of low molecular weight nitrosamines in pharmaceutical products by fast gas chromatography mass spectrometry. *J. Chromatogr. A* **1708**, 464323 (2023).
2. Schmidtsdorff, S., Neumann, J., Schmidt, A.H., *et al.* Prevalence of nitrosamine contaminants in drug samples: Has the crisis been overcome? *Arch. Pharm. (Weinheim)* **356**(2), e2200484 (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, 03/06/2024

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