

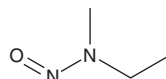
PRODUCT INFORMATION



N-Nitrosomethylethylamine

Item No. 36074

CAS Registry No.: 10595-95-6
Formal Name: N-methyl-N-nitroso-ethanamine
Synonym: NEMA
MF: $C_3H_8N_2O$
FW: 88.1
Purity: $\geq 95\%$
UV/Vis.: λ_{\max} : 229 nm
Supplied as: A liquid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

N-Nitrosomethylethylamine is supplied as a liquid. A stock solution may be made by dissolving the N-nitrosomethylethylamine in the solvent of choice, which should be purged with an inert gas. N-Nitrosomethylethylamine is slightly soluble in chloroform, ethyl acetate, and methanol.

Description

N-Nitrosomethylethylamine is an N-nitrosamine.¹ It has been found at low levels in some processed meat products.²

References

1. IARC monographs on the evaluation of the carcinogenic risk of chemicals to humans: Some N-nitroso compounds. (1978).
2. Niklas, A.A., Herrmann, S.S., Pedersen, M., *et al.* The occurrence of volatile and non-volatile N-nitrosamines in cured meat products from the Danish market. *Food Chem.* **378**, 132046 (2022).

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, 06/14/2022

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