

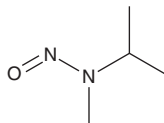
PRODUCT INFORMATION



N-Nitrosomethylisopropylamine

Item No. 36220

CAS Registry No.: 30533-08-5
Formal Name: N-methyl-N-nitroso-2-propanamine
Synonym: NMIPA
MF: $C_4H_{10}N_2O$
FW: 102.1
Purity: $\geq 90\%$
UV/Vis.: λ_{max} : 227 nm
Supplied as: A liquid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

N-Nitrosomethylisopropylamine is an N-nitrosamine.¹ It induces tumor formation in mice.

Reference

1. Dunn, W.J.I. and Wold, S. The carcinogenicity of N-nitroso compounds: A SIMCA pattern recognition study. *Bioor. Chem.* **10(1)**, 29-45 (1981).

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, 03/18/2026

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