

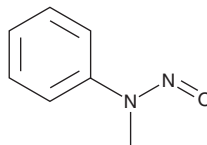
PRODUCT INFORMATION



N-Nitroso-N-methylaniline

Item No. 35563

CAS Registry No.: 614-00-6
Formal Name: N-methyl-N-nitroso-benzenamine
Synonyms: N-Methyl-N-nitrosoaniline, NMA, NSC 137
MF: C₇H₈N₂O
FW: 136.2
Purity: ≥98%
UV/Vis.: λ_{max}: 272 nm
Supplied as: A solution in ethanol
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-Nitroso-N-methylaniline is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. N-Nitroso-N-methylaniline is soluble in methanol, chloroform, and DMSO.

Description

N-Nitroso-N-methylaniline is an N-nitrosamine.¹ It induces the formation of esophageal, pharyngeal, and tongue tumors in rats when administered at a concentration of 200 mg/L in the drinking water or 15 mg/ml via oral gavage.² N-Nitroso-N-methylaniline is toxic to rats with LD₅₀ values of 336 and 225 mg/kg for male and female rats, respectively. It reduces survival of hamsters but only induces a low incidence of tumors in this species when administered at a dose of 5 mg/week for a total dose of 1.8 mmol.¹

References

1. Lijinsky, W. and Kovatch, R.M. Comparative carcinogenesis by nitrosomethylalkylamines in Syrian hamsters. *Cancer Res.* **48(23)**, 6648-6652 (1988).
2. Goodall, C.M., Lijinsky, W., Tomatis, L., et al. Toxicity and oncogenicity of nitrosomethylaniline and nitrosomethylcyclohexylamine. *Toxicol. Appl. Pharmacol.* **17(2)**, 426-432 (1970).

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, 12/12/2022

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