

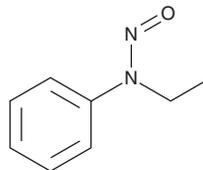
PRODUCT INFORMATION



N-Nitroso-N-ethylaniline

Item No. 21063

CAS Registry No.: 612-64-6
Formal Name: N-ethyl-N-nitroso-benzenamine
Synonyms: N-Ethyl-N-nitrosoaniline, NSC 405070
MF: C₈H₁₀N₂O
FW: 150.2
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 273 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-ethylaniline 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. Solvents such as DMSO purged with an inert gas can be used. The solubility of N-nitroso-N-ethylaniline in DMSO is approximately 5 mg/ml.

Description

N-Nitroso-N-ethylaniline is a member of the N-nitrosamine class of compounds, which are classified as carcinogens by the Environmental Protection Agency.¹ However, N-nitroso-N-ethylaniline is non-carcinogenic in a survival assay using 2FR₄50 Rauscher leukemia virus-infected rat embryo cultures.² It is a synthetic intermediate in the synthesis of 3-amino-4-hydrazine-cyclobut-3-ene-1,2-diones as antagonists of chemokine (C-X-C motif) receptor 2 (CXCR2).³

References

1. Park, J.-e., Seo, J.-e., Lee, J.-y., *et al.* Distribution of seven N-nitrosamines in food. *Toxicol. Res.* **31(3)**, 279-288 (2015).
2. Traul, K., Takayama, F., Kachevsky, V., *et al.* A rapid *in vitro* assay for carcinogenicity of chemical substances in mammalian cells utilizing an attachment-independence endpoint. 2 - Assay validation. *J. Appl. Toxicol.* **1(3)**, 190-195 (1981).
3. Liu, S., Liu, Y., Wang, H., *et al.* Design, synthesis, and evaluation of novel 3-amino-4-hydrazine-cyclobut-3-ene-1,2-diones as potent and selective CXCR2 chemokine receptor antagonists. *Bioorg. Med. Chem. Lett.* **19(19)**, 5741-5745 (2009).

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, 11/01/2022

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