

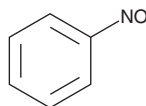
PRODUCT INFORMATION



Nitrosobenzene

Item No. 21009

CAS Registry No.: 586-96-9
Formal Name: nitroso-benzene
Synonym: NSC 66479
MF: C_6H_5NO
FW: 107.1
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 218, 281, 303 nm
Supplied as: A crystalline solid
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

Nitrosobenzene is supplied as a crystalline solid. A stock solution may be made by dissolving the nitrosobenzene in the solvent of choice, which should be purged with an inert gas. Nitrosobenzene is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nitrosobenzene in these solvents is approximately 10 mg/ml.

Nitrosobenzene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nitrosobenzene should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nitrosobenzene has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nitrosobenzene is a spin trap that has been used in the study of oxidative DNA damage and nitroso-compound-induced respiratory burst in neutrophils.¹

Reference

1. Zuman, P. and Shah, B. Addition, reduction, and oxidation reactions of nitrosobenzene. *Chem. Rev.* **94**(6) 1621-1641 (1994).

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