

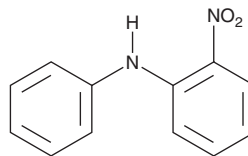
PRODUCT INFORMATION



2-Nitrodiphenylamine

Item No. 45198

CAS Registry No.: 119-75-5
Formal Name: 2-nitro-N-phenyl-benzenamine
Synonyms: C.I. 10335, NDPA, 2-nDPA, 2-nitroDPA, 2-NO₂-DPA, NSC 4754, NSC 105613, NSC 629274
MF: C₁₂H₁₀N₂O₂
FW: 214.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

2-Nitrodiphenylamine is supplied as a solid. A stock solution may be made by dissolving the 2-nitrodiphenylamine in the solvent of choice, which should be purged with an inert gas. 2-Nitrodiphenylamine is soluble (≥10 mg/ml) in DMSO. 2-Nitrodiphenylamine is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

2-Nitrodiphenylamine is a non-*p*-phenylenediamine (non-PPD) aromatic amine antioxidant.¹ It has been found in soil, ground water, and marine sediment.^{1,2} Formulations containing 2-nitrodiphenylamine have been used as stabilizing additives in torpedo propellants and as antioxidant additives in rubber.

References

1. Drzyzga, O. Diphenylamine and derivatives in the environment: A review. *Chemosphere* **53**(8), 809-818 (2003).
2. Jin, R., Li, Y., Saito, Y., *et al.* Amino accelerators and antioxidants in sediments from the Dong Nai River System, Vietnam: Distribution and influential factors. *Environ. Chem.* **227**, 115712 (2023).

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, 04/06/2026

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