

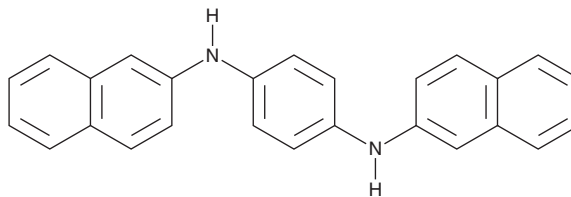
PRODUCT INFORMATION



N,N'-Di-2-naphthyl-*p*-phenylenediamine

Item No. 40728

CAS Registry No.: 93-46-9
Formal Name: N¹,N⁴-di-2-naphthalenyl-1,4-benzenediamine
Synonyms: DNPd, NSC 3410
MF: C₂₆H₂₀N₂
FW: 360.5
Purity: ≥90%
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

N,N'-Di-2-naphthyl-*p*-phenylenediamine (DNPd) is supplied as a solid. A stock solution may be made by dissolving the DNPd in the solvent of choice, which should be purged with an inert gas. DNPd is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DNPd is a *p*-phenylenediamine (PPD).¹ It has been found in fine particulate matter 2.5 (PM_{2.5}) samples collected at the roadside in China.

Reference

1. Jiang, N., Hao, X., Wang, Z., *et al.* *p*-Phenylenediamine antioxidants in PM_{2.5}: New markers for traffic in positive matrix factorization source apportionment. *J. Hazard Mater.* **476**, 135122 (2024).

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, 02/04/2026

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