

PRODUCT INFORMATION

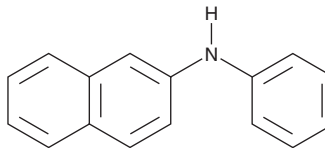


N-Phenyl-2-naphthylamine

Item No. 42580

CAS Registry No.: 135-88-6
Formal Name: N-phenyl-2-naphthalenamine
Synonyms: Aceto PBN, 2-Anilinonaphthalene, Antioxidant 116, Antioxidant D, Antioxidant PBN, 2-Naphthylphenylamine, β -Naphthylphenylamine, NSC 37151, P2NA, PBNA, PNA

MF: $C_{16}H_{13}N$
FW: 219.3
Purity: $\geq 95\%$
Supplied as: A 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

N-Phenyl-2-naphthylamine (P2NA) is supplied as a solid. A stock solution may be made by dissolving the P2NA in the solvent of choice, which should be purged with an inert gas. P2NA is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

P2NA is an aromatic amine antioxidant.^{1,2} It is non-carcinogenic, however, it can be dephenylated *in vivo* to 2-naphthylamine (2NA), a carcinogen associated with bladder cancer. P2NA was previously used as a common additive in rubber.

References

1. Weiss, T., Bolt, H.M., Schlüter, G., *et al.* Metabolic dephenylation of the rubber antioxidant N-phenyl-2-naphthylamine to carcinogenic 2-naphthylamine in rats. *Arch. Toxicol.* **87**(7), 1265-1272 (2013).
2. Marek, E.M., Koslitz, S., Weiss, T., *et al.* Quantification of N-phenyl-2-naphthylamine by gas chromatography and isotope-dilution mass spectrometry and its percutaneous absorption *ex vivo* under workplace conditions. *Arch. Toxicol.* **91**(11), 3587-3596 (2017).

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, 03/25/2025

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