

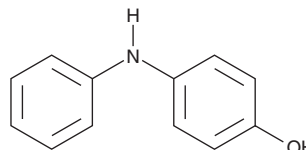
PRODUCT INFORMATION



4-Hydroxydiphenylamine

Item No. 40730

CAS Registry No.: 122-37-2
Formal Name: 4-(phenylamino)-phenol
Synonyms: 4-HDPA, *p*-Hydroxydiphenylamine, *para*-Hydroxydiphenylamine, NSC 1543
MF: C₁₂H₁₁NO
FW: 185.2
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

4-Hydroxydiphenylamine (4-HDPA) is supplied as a solid. A stock solution may be made by dissolving the 4-HDPA in the solvent of choice, which should be purged with an inert gas. 4-HDPA is soluble (≥10 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 4-HDPA can be prepared by directly dissolving the solid in aqueous buffers. 4-HDPA is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

4-HDPA is an aqueous transformation product of the substituted *p*-phenylenediamine and antiozonant 6-PPD (Item No. 38246).¹ It is also a biotransformation product of 6-PPD in zebrafish embryos and an active metabolite of the antioxidant and pesticide diphenylamine (Item No. 24250).^{2,3} 4-HDPA scavenges carbon radicals in an AIBN-induced methyl methacrylate polymerization assay in a concentration-dependent manner.⁴ It has been found in tire wear particles, roadway runoff, and roadway-impacted creeks.¹

References

1. Zhao, H.N., Hu, X., Tian, Z., *et al.* Transformation products of tire rubber antioxidant 6PPD in heterogeneous gas-phase ozonation: Identification and environmental occurrence. *Environ. Sci. Technol.* **57**(14), 5621-5632 (2023).
2. Grasse, N., Seiwert, B., Massei, R., *et al.* Uptake and biotransformation of the tire rubber-derived contaminants 6-PPD and 6-PPD quinone in the zebrafish embryo (*Danio rerio*). *Environ. Sci. Technol.* **57**(41), 15598-15607 (2023).
3. Alexander, W.E., Ryan, A.J., and Wright, S.E. Metabolism of diphenylamine in the rat, rabbit and man. *Food Cosmet. Toxicol.* **3**(4), 571-579 (1965).
4. Fujisawa, S., Ishihara, M., and Kadoma, Y. Antioxidant activity of diphenylamine-related compounds as scavengers of carbon radicals. *J. Mol. Des.* **4**, 711-720 (2005).

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, 04/16/2024

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