

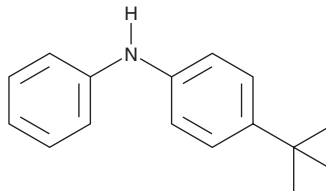
PRODUCT INFORMATION



BDPA

Item No. 45202

CAS Registry No.: 4496-49-5
Formal Name: 4-(1,1-dimethylethyl)-N-phenylbenzenamine
Synonyms: 4-*tert*-butyl-N-Phenylaniline, 4-*tert*-Butyldiphenylamine
MF: C₁₆H₁₉N
FW: 225.3
Purity: ≥95%
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

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

Description

BDPA is a substituted diphenylamine (SDPA) amino accelerator and antioxidant (AAL/O), which is a class of compounds used in tire rubber, plastics, and lubricants.¹ It has been found in road dust and river estuary sediment.^{1,2} BDPA has also been found in human plasma.³

References

1. Jin, R., Wu, Y., He, Q., *et al.* Ubiquity of amino accelerators and antioxidants in road dust from multiple land types: Targeted and nontargeted analysis. *Environ. Sci. Technol.* **57**(28), 10361-10372 (2023).
2. Jin, R., Li, B., Wu, Y., *et al.* Unpuzzling spatio-vertical and multi-media patterns of aniline accelerators/antioxidants in an urban estuary. *Water Res.* **266**, 122427 (2024).
3. Liang, B., Deng, Q., Pan, Z., *et al.* Solid-phase extraction based on Captiva EMR-lipid for determination of 19 aromatic amine antioxidants and two p-phenylenediamine quinones in human plasma. *Toxics* **14**(3), 187 (2026).

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, 03/18/2026

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