

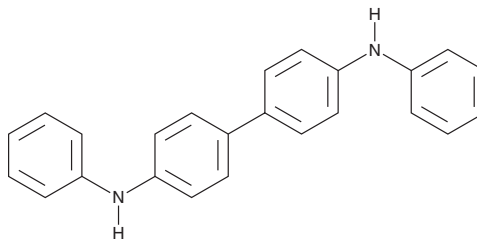
PRODUCT INFORMATION



DPB

Item No. 45199

CAS Registry No.: 531-91-9
Formal Name: N⁴,N⁴-diphenyl-[1,1'-biphenyl]-4,4'-diamine
Synonyms: Diphenylbenzidine, N,N'-Diphenylbenzidine
MF: C₂₄H₂₀N₂
FW: 336.4
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

DPB is supplied as a solid. A stock solution may be made by dissolving the DPB in the solvent of choice, which should be purged with an inert gas. DPB is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1) and DMSO.

Description

DPB is an aromatic diamine that has been used as a synthetic intermediate in the synthesis of soluble and electroactive polymers.¹

Reference

1. Niu, J., Lee, J.Y., Neudeck, A., *et al.* Electrochemical and optical characteristics of the hole conducting material N,N'-diphenylbenzidine during oxidation in DMF. *Synth. Met.* **99(2)**, 133-142 (1999).

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, 02/29/2026

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