

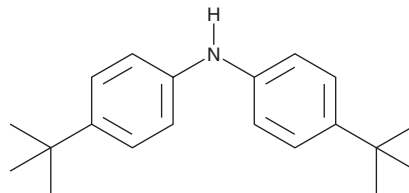
PRODUCT INFORMATION



DBDPA

Item No. 45197

CAS Registry No.: 4627-22-9
Formal Name: 4-(1,1-dimethylethyl)-N-[4-(1,1-dimethylethyl) phenyl]-benzenamine
Synonym: 4,4'-di-*tert*-butyl Diphenylamine
MF: C₂₀H₂₇N
FW: 281.4
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

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

Description

DBDPA is an aromatic amine antioxidant.¹ It has been found in children's car seats.

Reference

1. Wu, Y. and Venier, M. High levels of synthetic antioxidants and ultraviolet filters in children's car seats. *Sci. Total Environ.* **855**, 158637 (2023).

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, 02/12/2026

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