

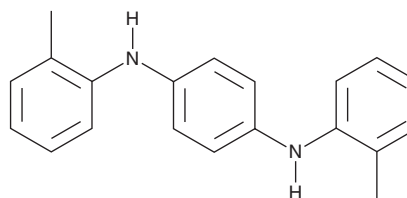
PRODUCT INFORMATION



DTPD

Item No. 40756

CAS Registry No.: 15017-02-4
Formal Name: N¹, N⁴-bis(2-methylphenyl)-1,4-benzenediamine
Synonyms: N¹,N⁴-bis(2-Methylphenyl)benzene-1,4-diamine, N,N'-di(o-tolyl)-p-phenylenediamine
MF: C₂₀H₂₀N₂
FW: 288.4
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

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

Description

DTPD is a substituted *p*-phenylenediamine and non-oxidized version of DTPD-Q (Item No. 40612).¹ It has been found in water treatment plant influent, effluent, biosolids, and landfill leachates, indoor dust samples, and particulate matter 2.5 (PM_{2.5}) in cities and by roadsides.¹⁻³

References

1. Zhang, Z.-F., Zhang, X., Zhang, X., *et al.* Diphenylamine antioxidants in wastewater influent, effluent, biosolids and landfill leachate: Contribution to environmental releases. *Water Res.* **189**, 116602 (2021).
2. Zhu, J., Guo, R., Jiang, S., *et al.* Occurrence of *p*-phenylenediamine antioxidants (PPDs) and PPDs-derived quinones in indoor dust. *Sci. Total Environ.* **912**, 169325 (2024).
3. Wang, W., Cao, G., Zhang, J., *et al.* Beyond substituted *p*-phenylenediamine antioxidants: Prevalence of their quinone derivatives in PM_{2.5}. *Environ. Sci. Technol.* **56(15)**, 10629-10637 (2022).

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, 04/29/2024

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