

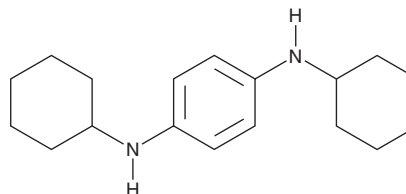
PRODUCT INFORMATION



CCPD

Item No. 42610

CAS Registry No.: 4175-38-6
Formal Name: N¹,N⁴-dicyclohexyl-1,4-benzenediamine
Synonym: N,N'-Dicyclohexyl-*p*-phenylenediamine
MF: C₁₈H₂₈N₂
FW: 272.4
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

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

Description

CCPD is a *p*-phenylenediamine (PPD).¹ It induces microcephaly and toxicity in chicken egg embryos when used at a concentration of 4.4 μM/egg. CCPD has been used as an antiozonant in rubber.²

References

1. Korhonen, A., Hemminki, K., and Vainio, H. Embryotoxicity of sixteen industrial amines to the chicken embryo. *J. Appl. Toxicol.* **3**(2), 112-117 (1983).
2. Dunn, J.R. The oxidative aging in ultraviolet light of dicumyl peroxide vulcanizates of natural rubber in presence of various antioxidants. *J. Appl. Polym. Sci.* **4**(11), 151-158 (1960).

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, 01/30/2025

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