

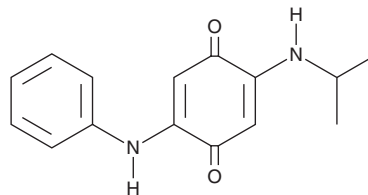
PRODUCT INFORMATION



IPPD-Q

Item No. 40829

CAS Registry No.: 68054-73-9
Formal Name: 2-[(1-methylethyl)amino]-5-(phenylamino)-2,5-cyclohexadiene-1,4-dione
Synonym: N-Isopropyl-N'-phenyl-p-phenylenediamine-quinone
MF: C₁₅H₁₆N₂O₂
FW: 256.3
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

IPPD-Q is supplied as a solid. A stock solution may be made by dissolving the IPPD-Q in the solvent of choice, which should be purged with an inert gas. IPPD-Q is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

IPPD-Q is a *p*-phenylenediamine quinone (PPD-Q) and an oxidized derivative of the antiozonant IPPD (Item No. 40725).¹ It is toxic to the bacterium *V. fischeri* (EC₅₀ = 2.97 mg/L). It has been found in fine particulate matter 2.5 (PM_{2.5}) samples collected at the roadside in China.²

References

1. Wang, W., Chen, Y., Fang, J., *et al.* Toxicity of substituted *p*-phenylenediamine antioxidants and their derived novel quinones on aquatic bacterium: Acute effects and mechanistic insights. *J. Hazard. Mater.* **469:133900**, (2024).
2. 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, 05/20/2024

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