

PRODUCT INFORMATION



6-PPD-Q-1-OH

Item No. 41591

Formal Name: 2-((1-hydroxy-4-methylpentan-2-yl)amino)-5-(phenylamino)cyclohexa-2,5-diene-1,4-dione

Synonym: 6-PPD-quinone-1-OH

MF: C₁₈H₂₂N₂O₃

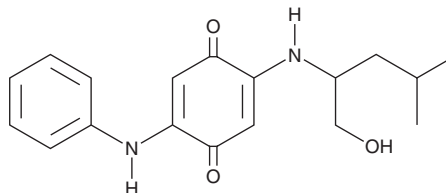
FW: 314.4

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

6-PPD-Q-1-OH is supplied as a solid. A stock solution may be made by dissolving the 6-PPD-Q-1-OH in the solvent of choice, which should be purged with an inert gas. 6-PPD-Q-1-OH is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in chloroform and DMSO.

Description

6-PPD-Q-1-OH is an oxidized derivative of the oxidized *p*-phenylenediamine 6-PPD-Q (Item No. 38247).¹

Reference

1. Ankley, P.J., da Silva, F.C., Jr., Montgomery, D., *et al.* Biotransformation of 6PPD-quinone *in vitro* using RTL-W1 cell line. *Environ. Sci. Tech. Lett.* **11(7)**, 650-769 (2024).

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, 08/08/2025

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