

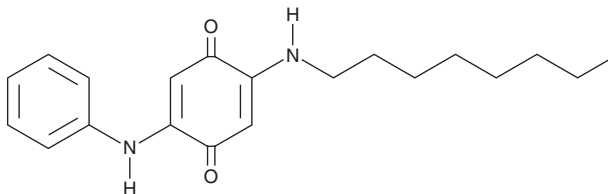
PRODUCT INFORMATION



OPPD-Q

Item No. 40614

CAS Registry No.: 2952779-34-7
Formal Name: 2-(octylamino)-5-(phenylamino)-2,5-cyclohexadiene-1,4-dione
Synonyms: 4-(2-Octylamino)diphenylamine-quinone, OPPD-quinone
MF: C₂₀H₂₆N₂O₂
FW: 326.4
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

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

Description

OPPD-Q is a *p*-phenylenediamine quinone (PPD-Q) and an oxidized derivative of the aromatic amine antioxidant OPPD (Item No. 40732).¹ Unlike 6-PPD-Q (Item No. 38247), OPPD-Q is not toxic to rainbow trout (*O. mykiss*) when used at concentrations up to 50 µg/L for 96 hours.

Reference

1. Nair, P., Sun, J., Xie, L., *et al.* Synthesis and toxicity evaluation of *p*-phenylenediamine-quinones. *Environ. Sci. Technol.* **59**(15), 7485-7494 (2025).

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, 11/14/2025

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