

PRODUCT INFORMATION



6-PPD-Q-yne

Item No. 45952

Formal Name: 2-((5-methylhex-1-yn-3-yl)amino)-5-(phenylamino)cyclohexa-2,5-diene-1,4-dione

Synonym: 6-PPD-quinone-yne

MF: C₁₉H₂₀N₂O₂

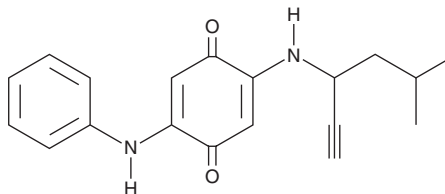
FW: 308.4

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

6-PPD-Q-yne is supplied as a solid. A stock solution may be made by dissolving the 6-PPD-Q-yne in the solvent of choice, which should be purged with an inert gas. 6-PPD-Q-yne is soluble (≥10 mg/ml) in DMSO.

Description

6-PPD-Q-yne is a clickable form of the environmental toxicant 6-PPD-Q (Item Nos. 38247 | 45543). It has been used to study the protein targets and neurotoxicity of 6-PPD-Q in rainbow trout.¹

Reference

1. Cheng, R., Zhang, W., Su, Y., *et al.* Chemoproteomic profiling reveals 6PPDQ-mediated neurotoxicity in rainbow trout through disruption of the glutamate-glutamine cycle. *Environ. Sci. Technol.* **60**(19), 13873-13884 (2026).

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, 07/16/2026

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