

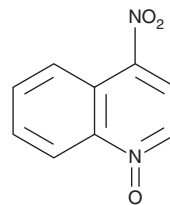
PRODUCT INFORMATION



4-NDQ

Item No. 42994

CAS Registry No.: 56-57-5
Formal Name: 4-nitroquinoline 1-oxide
Synonyms: Nitrochin, 4-Nitroquinoline N-oxide, 4-NQO, NSC 19645
MF: C₉H₆N₂O₃
FW: 190.2
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

4-NDQ is supplied as a solid. A stock solution may be made by dissolving the 4-NDQ in the solvent of choice, which should be purged with an inert gas. 4-NDQ is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

4-NDQ is a carcinogen.¹ It forms adducts with adenine and guanine, which can be removed by base excision repair, in *E. coli* DNA. 4-NDQ (100 µg/ml in the drinking water) induces tongue cancer in mice.²

References

1. Ikenaga, M., Ichikawa-Ryo, H., and Kondo, S. The major cause of inactivation and mutation by 4-nitroquinoline 1-oxide in *Escherichia coli*: Excisable 4NQO-purine adducts. *J. Mol. Biol.* **92(2)**, 341-356 (1975).
2. Tang, X.-H., Knudsen, B., Bemis, D., et al. Oral cavity and esophageal carcinogenesis modeled in carcinogen-treated mice. *Clin. Cancer Res.* **10(Pt 1)**, 301-313 (2004).

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, 03/27/2025

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