

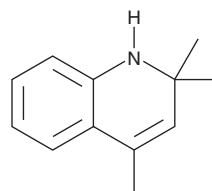
PRODUCT INFORMATION



TMQ

Item No. 42736

CAS Registry No.: 26780-96-1
Formal Name: 1,2-dihydro-2,2,4-trimethyl-quinoline, homopolymer
Synonym: Poly(1,2-dihydro-2,2,4-trimethylquinoline)
MF: (C₁₂H₁₅N)_x
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

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

Description

TMQ is an antioxidant.¹ It induces acanthosis at the site of administration but does not induce cancer in mice when administered at a dose of 100 mg/kg per day for two years.² TMQ has been used as an antioxidant and antiozonant in tires.

References

1. Ferradino, A.G. Antioxidant selection for peroxide cure elastomer applications. *Rubber Chemistry and Technology* **76**(3), 694-718 (2003).
2. Irwin, R.D., French, J.E., Elwell, M., *et al.* Carcinogenicity of dermally administered 1,2-dihydro-2,2,4-trimethylquinoline monomer in F344 rats and B6C3F₁ mice. *J. Appl. Toxicol.* **19**(2), 125-132 (1999).

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, 05/20/2025

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