

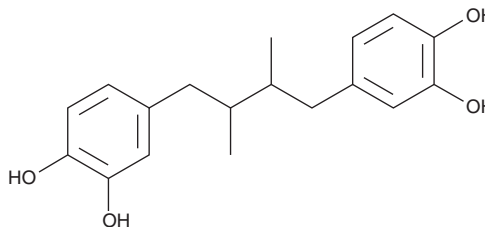
Product Information



Nordihydroguaiaretic Acid

Catalog No. 70300

CAS Registry No.: 500-38-9
Formal Name: 4,4'-(2,3-dimethyl-1,4-butanediyl)bis-1,2-benzenediol
Synonym: NDGA
MF: C₁₈H₂₂O₄
FW: 302.4
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that Nordihydroguaiaretic Acid (NDGA) be stored as supplied at -20°C. It should be stable for at least one year.

Concentrated stock solutions of NDGA can be prepared by dissolving the crystalline solid in an organic solvent such as ethanol, methanol, acetone, DMSO, or acetonitrile. The solubility of NDGA in these solvents is approximately 100 mg/ml. Stock solutions prepared in oxygen-free solvents will be stable for at least one week if protected from light and stored at 0-4°C.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Also, ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

NDGA is a non-selective lipoxygenase inhibitor which blocks cysteinyl-leukotriene synthesis. NDGA inhibits A23187-induced cysteinyl-leukotriene biosynthesis in rat peritoneal cells with an IC₅₀ value of 5-7 μM.¹ The IC₅₀ values are 3.0-5.0 μM for human platelet 12-lipoxygenase and 0.91 μM for rabbit reticulocyte 15-lipoxygenase.^{1,2}

References

1. Hope, W.C., Welton, A.F., Fiedler-Nagy, C., *et al.* *In vitro* inhibition of the biosynthesis of slow reacting substance of anaphylaxis (SRS-A) and lipoxygenase activity by quercetin. *Biochem. Pharmacol.* **32**, 367-371 (1983).
2. Pavani, M., Fones, E., Oksenberg, D., *et al.* Inhibition of tumoral cell respiration and growth by nordihydroguaiaretic acid. *Biochem. Pharmacol.* **48**, 1935-1942 (1994).

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 01/14/2010