

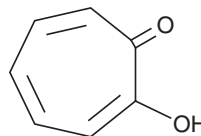
# PRODUCT INFORMATION



## Tropolone

Item No. 29674

**CAS Registry No.:** 533-75-5  
**Formal Name:** 2-hydroxy-2,4,6-cycloheptatrien-1-one  
**Synonyms:** 2-Hydroxytropone,  $\alpha$ -Tropolone, NSC 89303, Purpurocatechol  
**MF:** C<sub>7</sub>H<sub>6</sub>O<sub>2</sub>  
**FW:** 122.1  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 227, 319, 352 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Tropolone is supplied as a solid. A stock solution may be made by dissolving the tropolone in the solvent of choice, which should be purged with an inert gas. Tropolone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tropolone in these solvents is approximately 30, 10, and 5 mg/ml, respectively.

Tropolone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tropolone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Tropolone has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Tropolone is a terpene that has been found in cupressaceous tree bark and has diverse biological activities.<sup>1-5</sup> It inhibits catechol O-methyltransferase (COMT; K<sub>i</sub> = 22  $\mu$ M).<sup>1</sup> Tropolone is bacteriostatic against a panel of Gram-positive and Gram-negative bacteria (MICs = 12-82  $\mu$ M) and bactericidal against *A. hydrophilia* and *C. violaceum* when used at a concentration of 20 mM.<sup>2</sup> It inhibits LPS-induced production of nitric oxide (NO) in RAW 264.7 cells (IC<sub>50</sub> = 12  $\mu$ M).<sup>3</sup> Tropolone (0.25 mmol/animal) reduces thyroid radioiodide uptake in rats.<sup>4</sup> Dietary administration of tropolone induces growth retardation and hyperactivity in mice.

### References

1. Borchardt, R.T. Catechol O-methyltransferase. 1. Kinetics of tropolone inhibition. *J. Med. Chem.* **16**(4), 377-382 (1973).
2. Trust, T.J. Antibacterial activity of tropolone. *Antimicrob. Agents Chemother.* **7**(5), 500-506 (1975).
3. Nishishiro, M., Kurihara, T., Wakabayashi, H., *et al.* Effect of tropolone, azulene and azulenequinone derivatives on prostaglandin E<sub>2</sub> production by activated macrophage-like cells. *Anticancer Res.* **29**(1), 379-383 (2009).
4. Lee, C.P., Hegarty, M.P., and Christie, G.S. Antithyroid and antiperoxidase activity of tropolone and 3-hydroxy-4-pyrone. *Chem. Biol. Interact.* **27**(1), 17-26 (1979).

#### 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, 08/25/2023

#### 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