

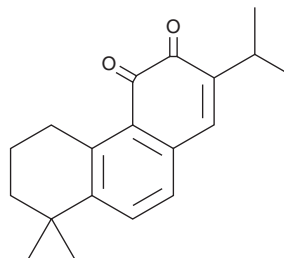
# PRODUCT INFORMATION



## Miltirone

Item No. 36663

**CAS Registry No.:** 27210-57-7  
**Formal Name:** 5,6,7,8-tetrahydro-8,8-dimethyl-2-(1-methylethyl)-3,4-phenanthrenedione  
**Synonyms:** NSC 639662, Rosmariquinone  
**MF:** C<sub>19</sub>H<sub>22</sub>O<sub>2</sub>  
**FW:** 282.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 258 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

Miltirone is supplied as a solid. A stock solution may be made by dissolving the miltirone in the solvent of choice, which should be purged with an inert gas. Miltirone is soluble in chloroform and ethyl acetate.

### Description

Miltirone is a tanshinone that has been found in *S. miltiorrhiza* and has diverse biological activities.<sup>1-5</sup> It protects lard from thermal oxidation when used at a concentration of 0.02%.<sup>1</sup> Miltirone is a partial agonist of the benzodiazepine receptor (IC<sub>50</sub> = 0.3 μM).<sup>2</sup> It inhibits carboxylesterase 2 (CES2; K<sub>i</sub> = 0.04 μM) as well as severe acute respiratory syndrome coronavirus (SARS-CoV) main protease (M<sup>pro</sup>), also known as 3C-like protease (3CL<sup>pro</sup>), and papain-like protease (PL<sup>pro</sup>; IC<sub>50</sub>s = 21.1 and 30 μM, respectively).<sup>3,4</sup> Miltirone is cytotoxic to HCT116 colon cancer cells (IC<sub>50</sub> = 4.27 μM) and induces mitochondrial dysfunction and the production of reactive oxygen species (ROS) in the same cells.<sup>5</sup> *In vivo*, miltirone (10-60 mg/kg) increases the number of punished crossings in the four-plate test, indicating tranquilizing activity, in mice.<sup>2</sup>

### References

1. Houlihan, C.M., Ho, C.-T., and Chang, S.S. The structure of rosmariquinone - A new antioxidant isolated from *Rosmarinus officinalis* L. *JAOCS* **62**(1), 96-98 (1985).
2. Lee, C.-M., Wong, H.N.C., Chui, K.-Y., et al. Miltirone, a central benzodiazepine receptor partial agonist from a Chinese medicinal herb *Salvia miltiorrhiza*. *Neurosci. Lett.* **127**(2), 237-241 (1991).
3. Hatfield, J.M., Tsurkan, L.G., Hyatt, J.L., et al. Modulation of esterified drug metabolism by tanshinones from *Salvia miltiorrhiza* ("Danshen"). *J. Nat. Prod.* **76**(1), 36-44 (2012).
4. Park, J.-Y., Kim, J.H., Kim, Y.M., et al. Tanshinones as selective and slow-binding inhibitors for SARS-CoV cysteine proteases. *Bioorg. Med. Chem.* **20**(19), 5928-5935 (2012).
5. Wang, L., Hu, T., Shen, J., et al. Miltirone induced mitochondrial dysfunction and ROS-dependent apoptosis in colon cancer cells. *Life Sci.* **151**, 224-234 (2016).

#### 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, 12/20/2022

#### 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