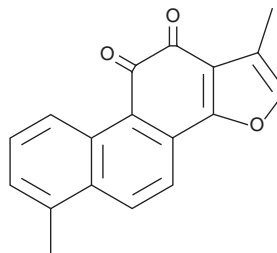


# PRODUCT INFORMATION



## Tanshinone I Item No. 26852

**CAS Registry No.:** 568-73-0  
**Formal Name:** 1,6-dimethyl-phenanthro[1,2-b]furan-10,11-dione  
**Synonym:** Tanshinone A  
**MF:** C<sub>18</sub>H<sub>12</sub>O<sub>3</sub>  
**FW:** 276.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 245, 420 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Salviae Miltiorrhizae Radix et Rhizoma*



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

### Laboratory Procedures

Tanshinone I is supplied as a crystalline solid. A stock solution may be made by dissolving the tanshinone I in the solvent of choice. Tanshinone I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of tanshinone I in these solvents is approximately 0.1, 0.3, and 1 mg/ml, respectively.

### Description

Tanshinone I is a diterpene that has been found in *S. miltiorrhiza* and has diverse biological activities, including anticancer, antidiabetic, and neuroprotective properties.<sup>1-4</sup> It inhibits proliferation in a variety of cancer cell lines, including KB/VCR, MCF-7/ADR, and K562/A02 multidrug-resistance cell lines.<sup>2</sup> It inhibits topoisomerase I- and topoisomerase II-mediated supercoiled DNA relaxation activity when used at concentrations of 200 and 25 μM, respectively. Tanshinone I also inhibits proliferation of H358-IR and H157-IR radioresistant lung cancer cells (IC<sub>50</sub>s = 10.87 and 6.1 μM, respectively) and sensitizes them to radiation.<sup>3</sup> It reduces plasma levels of total cholesterol, triglycerides, LDL, and nonesterified fatty acids and decreases blood glucose levels in a rat model of type 2 diabetes induced by streptozotocin (Item No. 13104) and a high-fat diet.<sup>5</sup> Tanshinone I reduces the LPS-induced expression of TNF-α, IL-1β, and IL-6 in microglia *in vitro*.<sup>6</sup> It is neuroprotective in a mouse model of Parkinson's disease induced by MPTP, preventing dopaminergic cell death in the striatum and improving motor function when administered at a dose of 10 mg/kg per day for two weeks.

### References

1. Yang, L., Ding, G., Lin, H., *et al.* Transcriptome analysis of medicinal plant *Salvia miltiorrhiza* and identification of genes related to tanshinone biosynthesis. *PLoS One* **8**(11), e80464 (2013).
2. Tian, Q.-T., Ding, C.-Y., Song, S.-S., *et al.* New tanshinone I derivatives S222 and S439 similarly inhibit topoisomerase I/II but reveal different p53-dependency in inducing G2/M arrest and apoptosis. *Biochem. Pharmacol.* **156**, 255-264 (2018).
3. Yan, Y., Su, W., Zeng, S., *et al.* Effect and mechanism of tanshinone I on the radiosensitivity of lung cancer cells. *Mol. Pharm.* **15**(11), 4843-4853 (2018).
4. Wang, S., Jing, H., Yang, H., *et al.* Tanshinone I selectively suppresses pro-inflammatory genes expression in activated microglia and prevents nigrostriatal dopaminergic neurodegeneration in a mouse model of Parkinson's disease. *J. Ethnopharmacol.* **164**, 247-255 (2015).
5. Wei, Y., Gao, J., Qin, L., *et al.* Tanshinone I alleviates insulin resistance in type 2 diabetes mellitus rats through IRS-1 pathway. *Biomed. Pharmacother.* **93**, 352-356 (2017).

#### 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