

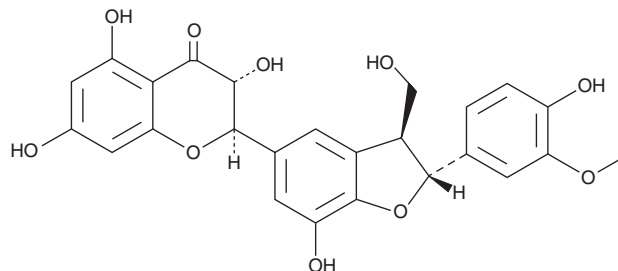
# PRODUCT INFORMATION



## Silychristin

Item No. 34658

**CAS Registry No.:** 33889-69-9  
**Formal Name:** (2R,3R)-2-[(2R,3S)-2,3-dihydro-7-hydroxy-2-(4-hydroxy-3-methoxyphenyl)-3-(hydroxymethyl)-5-benzofuranyl]-2,3,4H-1-benzopyran-4-one  
**Synonyms:** Silychristin A, (+)-Silychristin  
**MF:** C<sub>25</sub>H<sub>22</sub>O<sub>10</sub>  
**FW:** 482.4  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Silybum marianum*



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

### Laboratory Procedures

Silychristin is supplied as a solid. A stock solution may be made by dissolving the silychristin in the solvent of choice, which should be purged with an inert gas. Silychristin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of silychristin in these solvents is approximately 25 mg/ml.

Silychristin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, silychristin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Silychristin has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Silychristin is a flavonolignan that has been found in *S. marianum* and has diverse biological activities.<sup>1-3</sup> It inhibits monocarboxylate transporter 8 (MCT8; IC<sub>50</sub> = ~100 nM) and MCT8-dependent uptake of the thyroid hormone 3,3',5-triiodo-L-thyronine (T3) in MDCK cells and primary mouse astrocytes (IC<sub>50</sub>s = 170 and 50 nM, respectively).<sup>1</sup> Silychristin scavenges DPPH (Item No. 14805) radicals in a cell-free assay (IC<sub>50</sub> = 18.97 μM) and inhibits *tert*-butyl-induced lipid peroxidation in rat liver microsomes (IC<sub>50</sub> = 5.45 μM).<sup>2</sup> It also inhibits P-glycoprotein (P-gp) in P-gp-containing membranes (IC<sub>50</sub> = 21 μM) and LPS-induced production of nitric oxide (NO) in RAW 264.7 cells (IC<sub>50</sub> = 65 μM).<sup>3</sup>

### References

1. Johannes, J., Jayarama-Naidu, R., Meyer, F., *et al.* Silychristin, a flavonolignan derived from the milk thistle, is a potent inhibitor of the thyroid hormone transporter MCT8. *Endocrinology* **157**(4), 1694-1701 (2016).
2. Biedermann, D., Buchta, M., Holečková, V., *et al.* Silychristin: Skeletal alterations and biological activities. *J. Nat. Prod.* **79**(12), 3086-3092 (2016).
3. Viktorová, J., Dobiasová, S., Řehořová, K., *et al.* Antioxidant, anti-inflammatory, and multidrug resistance modulation activity of silychristin derivatives. *Antioxidants* **8**(8), 303 (2019).

**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, 11/30/2022

### 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