

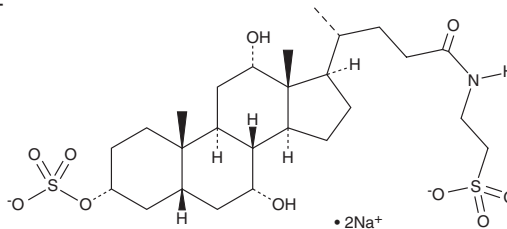
# PRODUCT INFORMATION



## Taurocholic Acid 3-sulfate (sodium salt)

Item No. 9002538

**CAS Registry No.:** 71781-33-4  
**Formal Name:** 2-[[[(3 $\alpha$ ,5 $\beta$ ,7 $\alpha$ ,12 $\alpha$ )-7,12-dihydroxy-24-oxo-3-(sulfooxy)cholan-24-yl]amino]ethanesulfonic acid, disodium salt  
**Synonyms:** 3-Sulfocholyl Taurine, TCA3S  
**MF:** C<sub>26</sub>H<sub>43</sub>NO<sub>10</sub>S<sub>2</sub> • 2Na  
**FW:** 639.7  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Taurocholic acid 3-sulfate (TCA3S) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the TCA3S (sodium salt) in the solvent of choice, which should be purged with an inert gas. TCA3S (sodium salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TCA3S (sodium salt) in these solvents is approximately 0.3 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of TCA3S (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of TCA3S (sodium salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

TCA3S is a metabolite of the conjugated bile acid taurocholic acid (Item No. 16215).<sup>1</sup> Plasma levels of TCA3S are elevated in wild-type and *Sortilin 1* (*Sort1*) knockout mice at 6 hours following bile duct ligation (BDL) and are further elevated in *Sort1* knockout mice at 24 hours post-BDL.<sup>2</sup>

### References

1. Lefebvre, P., Cariou, B., Lien, F., *et al.* Role of bile acids and bile acid receptors in metabolic regulation. *Physiol. Rev.* **89**(1), 147-191 (2009).
2. Li, J., Woolbright, B.L., Zhao, W., *et al.* Sortilin 1 loss-of-function protects against cholestatic liver injury by attenuating hepatic bile acid accumulation in bile duct ligated mice. *Toxicol. Sci.* **161**(1), 34-47 (2018).

#### 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, 03/08/2021

#### 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