

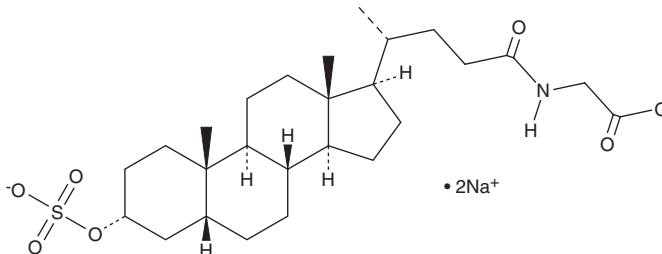
# PRODUCT INFORMATION



## Glycolithocholic Acid 3-sulfate (sodium salt)

Item No. 35298

**CAS Registry No.:** 64936-82-9  
**Formal Name:** N-[(3 $\alpha$ ,5 $\beta$ )-24-oxo-3-(sulfooxy)cholan-24-yl]-glycine, disodium salt  
**Synonyms:** GLCA 3-sulfate, Lithocholylglycine 3-sulfate, 3-Sulfoglycolithocholic Acid, Sulfolithocholylglycine  
**MF:** C<sub>26</sub>H<sub>41</sub>NO<sub>7</sub>S • 2Na  
**FW:** 557.7  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Glycolithocholic acid 3-sulfate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the glycolithocholic acid 3-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Glycolithocholic acid 3-sulfate (sodium salt) is slightly soluble in methanol.

Glycolithocholic acid 3-sulfate (sodium salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Glycolithocholic acid 3-sulfate is a metabolite of glycolithocholic acid (Item Nos. 21723 | 20273).<sup>1</sup> It is formed from glycolithocholic acid by hepatocytes.<sup>2</sup> Glycolithocholic acid 3-sulfate (240  $\mu$ mol/kg) induces cholestasis and increases biliary cholesterol secretion in rats.<sup>3</sup> Serum glycolithocholic acid 3-sulfate levels are increased in patients with cirrhosis, hepatitis, or cholestasis.<sup>4</sup>

### References

1. Palmer, R.H. The formation of bile acid sulfates: A new pathway of bile acid metabolism in humans. *Proc. Natl. Acad. Sci. USA* **58**(3), 1047-1050 (1967).
2. Kirkpatrick, R.B. and Belsaas, R.A. Formation and secretion of glycolithocholate-3-sulfate in primary hepatocyte cultures. *J. Lipid Res.* **26**(12), 1431-1437 (1985).
3. Yousef, I.M., Tuchweber, B., Vonk, R.J., et al. Lithocholate cholestasis-sulfated glycolithocholate-induced intrahepatic cholestasis in rats. *Gastroenterology* **80**(2), 233-241 (1981).
4. Demers, L.M. and Hepner, G.W. Levels of immunoreactive glycine-conjugated bile acids in health and hepatobiliary disease. *Am. J. Clin. Pathol.* **66**(5), 831-839 (1976).

#### 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, 03/17/2025

#### 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](http://WWW.CAYMANCHEM.COM)