

# PRODUCT INFORMATION



## Glycodeoxycholic Acid MaxSpec® Standard

Item No. 31599

CAS Registry No.: 360-65-6

Formal Name: N-[(3 $\alpha$ ,5 $\beta$ ,12 $\alpha$ )-3,12-dihydroxy-24-oxocholan-24-yl]-glycine

Synonym: GDCA

MF: C<sub>26</sub>H<sub>43</sub>NO<sub>5</sub>

FW: 449.6

Purity:  $\geq 95\%$

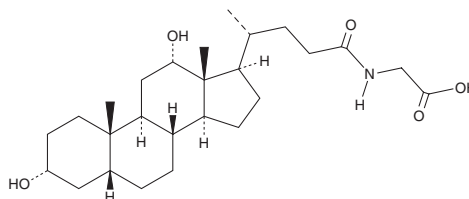
Supplied as: A solution in methanol; in a deactivated glass ampule

Concentration: 100  $\mu\text{g/ml}$  (nominal); see certificate of analysis for verified concentration

Storage: -20°C

Stability:  $\geq 5$  years; Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.

Special Conditions: Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.



### Description

Glycodeoxycholic acid (GDCA) is a glycine-conjugated form of the secondary bile acid deoxycholic acid (Item Nos. 18231 | 20756).<sup>1</sup> It induces a reversible, concentration-dependent reduction in myogenic tone in rats and decreases expression of the gene encoding the cytochrome P450 (CYP) isoform 7A1 (CYP7A1) in rabbits.<sup>2,3</sup> Serum levels of GDCA are elevated in non-surviving patients with acetaminophen-induced acute liver failure (AALF) compared with survivors.<sup>4</sup> GDCA levels are also increased in the plasma of patients with asthma.<sup>5</sup>

GDCA MaxSpec® standard is a quantitative grade standard of GDCA (Item No. 20274) that has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. The verified concentration is provided on the certificate of analysis. This GDCA MaxSpec® standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product. **Note:** The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.

### 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. Khurana, S., Raina, H., Pappas, V., *et al.* Effects of deoxycholyglycine, a conjugated secondary bile acid, on myogenic tone and agonist-induced contraction in rat resistance arteries. *PLoS One* **7**(2), e32006 (2012).
3. Shang, Q., Guo, G.L., Honda, A., *et al.* Bile acid flux through portal but not peripheral veins inhibits CYP7A1 expression without involvement of ileal FGF19 in rabbits. *Am. J. Physiol. Gastrointest. Liver Physiol.* **307**(4), G479-G486 (2014).
4. Woolbright, B.L., McGill, M.R., Staggs, V.S., *et al.* Glycodeoxycholic acid levels as prognostic biomarker in acetaminophen-induced acute liver failure patients. *Toxicol. Sci.* **142**(2), 436-444 (2014).
5. Comhair, S.A.A., McDunn, J., Bennett, C., *et al.* Metabolomic endotype of asthma. *J. Immunol.* **195**(2), 643-650 (2015).

#### 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, 07/10/2023

#### 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