

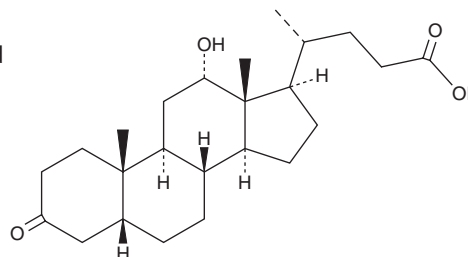
# PRODUCT INFORMATION



## 3-oxo Deoxycholic Acid

Item No. 35227

CAS Registry No.: 4185-01-7  
Formal Name: (5 $\beta$ )-12 $\alpha$ -hydroxy-3-oxo-cholan-24-oic acid  
Synonym: 3-oxo DCA  
MF: C<sub>24</sub>H<sub>38</sub>O<sub>4</sub>  
FW: 390.6  
Purity:  $\geq$ 95%  
Supplied as: A solid  
Storage: -20°C  
Stability:  $\geq$ 4 years



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

### Laboratory Procedures

3-oxo Deoxycholic acid is supplied as a solid. A stock solution may be made by dissolving the 3-oxo deoxycholic acid in the solvent of choice, which should be purged with an inert gas. 3-oxo Deoxycholic acid is slightly soluble in chloroform and methanol.

### Description

3-oxo Deoxycholic acid is a secondary bile acid and a metabolite of deoxycholic acid (Item Nos. 20756 | 18231) in *C. perfringens*, a bacteria found in the intestines of mammals and in the environment.<sup>1</sup>

### Reference

1. Hirano, S., Masuda, N., Oda, H., *et al.* Transformation of bile acids by *Clostridium perfringens*. *Appl. Environ. Microbiol.* **42(3)**, 394-399 (1981).

#### 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, 12/02/2022

#### 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