

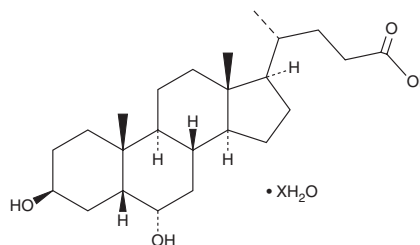
# PRODUCT INFORMATION



## $\beta$ -Hyodeoxycholic Acid (hydrate)

Item No. 35378

**Formal Name:** (5 $\beta$ )-3 $\beta$ ,6 $\alpha$ -dihydroxy-cholan-24-oic acid, hydrate  
**Synonyms:** 3 $\beta$ ,6 $\alpha$ -dihydroxy-5 $\beta$ -Cholan-24-oic Acid, isoHDCA, Isohyodeoxycholic Acid  
**MF:** C<sub>24</sub>H<sub>40</sub>O<sub>4</sub> • XH<sub>2</sub>O  
**FW:** 392.6  
**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

$\beta$ -Hyodeoxycholic acid (hydrate) is supplied as a solid. A stock solution may be made by dissolving the  $\beta$ -hyodeoxycholic acid (hydrate) in the solvent of choice, which should be purged with an inert gas.  $\beta$ -Hyodeoxycholic acid (hydrate) is slightly soluble in ethanol, methanol, and DMSO.

### Description

$\beta$ -Hyodeoxycholic acid is a 3 $\beta$  epimer of the secondary bile acid hyodeoxycholic acid (HDCA; Item No. 20294). Levels of  $\beta$ -hyodeoxycholic acid are increased in the feces of hamsters fed a diet containing HDCA- or hyocholic acid (HCA; Item No. 20293).<sup>1</sup>

### Reference

1. Cohen-Solal, C., Parquet, M., Férézou, J., *et al.* Effects of hyodeoxycholic acid and  $\alpha$ -hyocholic acid, two  $\alpha$ -hydroxylated bile acids, on cholesterol and bile acid metabolism in the hamster. *Biochim. Biophys. Acta* **1257**(2), 189-197 (1995).

#### 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, 10/25/2022

#### 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