

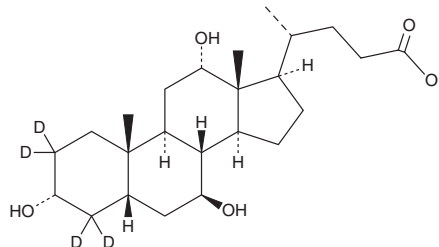
# PRODUCT INFORMATION



## Ursocholic Acid-d<sub>4</sub>

Item No. 35419

<b>Formal Name:</b>	(3 $\alpha$ ,5 $\beta$ ,7 $\beta$ ,12 $\alpha$ )-3,7,12-trihydroxy-cholan-24-oic-2,2,4,4-d <sub>4</sub>
<b>Synonyms:</b>	7-Epichoic Acid-d <sub>4</sub> , 7 $\beta$ -Hydroisocholic Acid-d <sub>4</sub> , 3 $\alpha$ ,7 $\beta$ ,12 $\alpha$ -Trihydroxy-5 $\beta$ -cholanoic Acid-d <sub>4</sub> , UCA-d <sub>4</sub>
<b>MF:</b>	C <sub>24</sub> H <sub>36</sub> D <sub>4</sub> O <sub>5</sub>
<b>FW:</b>	412.6
<b>Chemical Purity:</b>	≥98% (Ursocholic acid)
<b>Deuterium Incorporation:</b>	≥99% deuterated forms (d <sub>1</sub> -d <sub>4</sub> ); ≤1% d <sub>0</sub>
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥2 years



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

### Laboratory Procedures

Ursocholic acid-d<sub>4</sub> (UCA-d<sub>4</sub>) is intended for use as an internal standard for the quantification of UCA (Item No. 29883) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

UCA-d<sub>4</sub> is supplied as a solid. A stock solution may be made by dissolving the UCA-d<sub>4</sub> in the solvent of choice, which should be purged with an inert gas. UCA-d<sub>4</sub> is soluble in DMSO.

### Description

UCA is a bile acid and the 7 $\beta$ -hydroxyepimer of cholic acid (Item No. 20250).<sup>1</sup> Dietary administration of UCA (0.25%) decreases gallbladder cholesterol, phospholipid, and bile acid levels and the occurrence of gallstones in male mice fed a lithogenic diet.

### Reference

1. Uchida, K., Akiyoshi, T., Igmi, H., *et al.* Differential effects of ursodeoxycholic acid and ursocholic acid on the formation of biliary cholesterol crystals in mice. *Lipids* **26**(7), 526-530 (1991).

#### 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, 12/21/2021

#### 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