

PRODUCT INFORMATION

Hyodeoxycholic Acid MaxSpec® Standard

Item No. 31606

CAS Registry No.: 83-49-8

Formal Name: (3 α ,5 β ,6 α)-3,6-dihydroxy-cholan-24-oic acid

Synonyms: α -Hyodeoxycholic Acid, HDCA, NSC 60672

MF: C₂₄H₄₀O₄

FW: 392.6

Purity: $\geq 95\%$

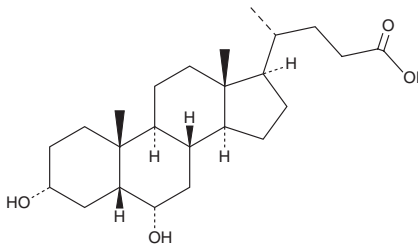
Supplied as: A solution in methanol; in a deactivated glass ampule

Concentration: 1 mg/ml (nominal); see certificate of analysis for verified concentration

Storage: -20°C

Stability: ≥ 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

Hyodeoxycholic acid (HDCA) is a secondary bile acid.¹ It is produced from lithocholic acid (Item No. 20253) by gut bacteria.¹⁻³ Dietary administration of HDCA (1.25% w/w) decreases plasma VLDL and LDL cholesterol levels and reduces fasting glucose levels and atherosclerotic lesion size in LDL receptor knockout mice fed a Western diet.⁴ Serum levels of HDCA are increased in patients with Crohn's disease or ulcerative colitis.⁵

Hyodeoxycholic acid MaxSpec® standard is a quantitative grade standard of HDCA (Item No. 20294) 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 hyodeoxycholic acid 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. Einarsson, K. On the formation of hyodeoxycholic acid in the rat. Bile acids and steroids 154. *J. Biol. Chem.* **241**(3), 534-539 (1966).
2. Madsen, D., Beaver, M., Chang, L., et al. Analysis of bile acids in conventional and germfree rats. *J. Lipid. Res.* **17**(2), 107-111 (1976).
3. Sacquet, E., Parquet, M., Riottot, M., et al. Intestinal absorption, excretion, and biotransformation of hyodeoxycholic acid in man. *J. Lipid. Res.* **24**(5), 604-613 (1983).
4. Shih, D.M., Shaposhnik, Z., Meng, Y., et al. Hyodeoxycholic acid improves HDL function and inhibits atherosclerotic lesion formation in LDLR-knockout mice. *FASEB J.* **27**(9), 3805-3817 (2013).
5. Gnewuch, C., Liebisch, G., Langmann, T., et al. Serum bile acid profiling reflects enterohepatic detoxification state and intestinal barrier function in inflammatory bowel disease. *World J. Gastroenterol.* **15**(25), 3134-3141 (2009).

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