

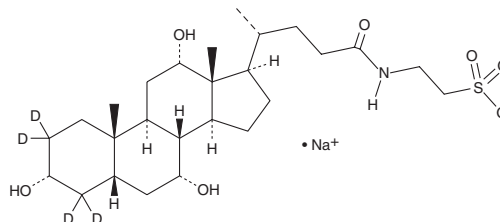
PRODUCT INFORMATION



Taurocholic Acid-d₄ (sodium salt)

Item No. 21891

CAS Registry No.: 2410279-93-3
Formal Name: 2-[[[(3 α ,5 β ,7 α ,12 α)-3,7,12-trihydroxy-24-oxocholan-24-yl]-2,2,4,4-d₄]amino]ethanesulfonic acid, monosodium salt
Synonym: Ethanesulfonic Acid-d₄, TCA-d₄
MF: C₂₆H₄₀D₄NO₇S • Na
FW: 541.7
Chemical Purity: ≥98% (Taurocholic Acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A solid
Storage: -20°C; Store in the dark in a tightly sealed container
Stability: ≥4 years
Special Conditions: Light sensitive



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

Laboratory Procedures

Taurocholic acid-d₄ (TCA-d₄) (sodium salt) is intended for use as an internal standard for the quantification of TCA (Item No. 16215) 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).

Description

TCA is a taurine-conjugated form of the primary bile acid cholic acid (Item No. 20250).¹ Serum levels of TCA are decreased in patients with Crohn's disease and those with ulcerative colitis with no extraintestinal manifestations, but are increased in patients with ulcerative colitis accompanied by hepatobiliary disease.²

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. 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, 11/11/2025

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