

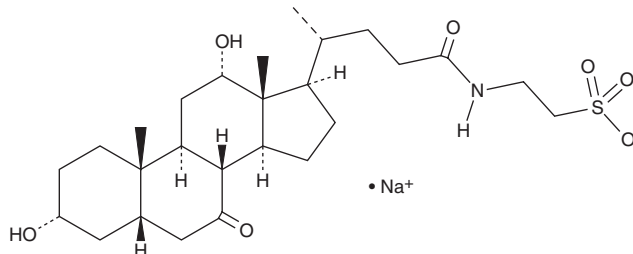
PRODUCT INFORMATION



Tauro-7-keto Deoxycholic Acid (sodium salt)

Item No. 9005003

CAS Registry No.: 2674426-87-8
Formal Name: 2-[[[(5 β)-3 α ,12 α -dihydroxy-7,24-dioxocholan-24-yl]amino]-ethanesulfonic acid, monosodium salt
Synonyms: Tauro-7-keto DCA, Tauro-7- oxo DCA, Tauro-7-oxo Deoxycholic Acid
MF: C₂₆H₄₂NO₇S • Na
FW: 535.7
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Tauro-7-keto deoxycholic acid (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the tauro-7-keto deoxycholic acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. Tauro-7-keto deoxycholic acid (sodium salt) is soluble in methanol.

Description

Tauro-7-keto deoxycholic acid is a taurine-conjugated form of 7-keto deoxycholic acid (Item No. 35264), a metabolite of the bile acid cholic acid (Item No. 20250).

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, 05/07/2026

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