

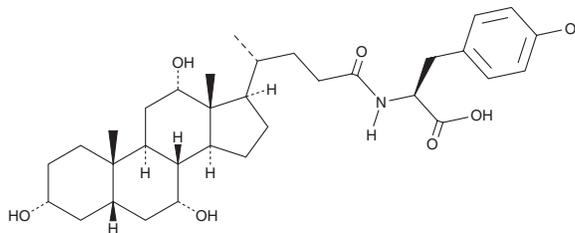
PRODUCT INFORMATION



N-Cholyl-L-Tyrosine

Item No. 44858

CAS Registry No.: 18485-51-3
Formal Name: N-[5 β -3 α ,7 α ,12 α -trihydroxy-24-oxocholan-24-yl]-L-tyrosine
Synonyms: N-Cholyl-Tyr, TyrCA
MF: C₃₃H₄₉NO₇
FW: 571.8
Purity: \geq 98% (mixture of epimers)
Supplied as: A crystalline 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

N-Cholyl-L-tyrosine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-cholyl-L-tyrosine in the solvent of choice, which should be purged with an inert gas. N-Cholyl-L-tyrosine is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

N-Cholyl-L-tyrosine is a bile acid conjugate composed of cholic acid (Item Nos. 20250 | 44008) and L-tyrosine (Item No. 36333).

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, 02/02/2026

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