

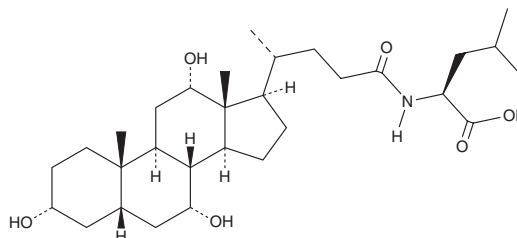
PRODUCT INFORMATION



N-Cholyl-L-Leucine

Item No. 40930

CAS Registry No.: 22154-46-7
Formal Name: N-[(3 α ,5 β ,7 α ,12 α)-3,7,12-trihydroxy-24-oxocholan-24-yl]-L-leucine
Synonyms: N-Cholyl-Leu, Leucocholic Acid
MF: C₃₀H₅₁NO₆
FW: 521.7
Purity: \geq 95%
Supplied as: A 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-leucine is supplied as a solid. A stock solution may be made by dissolving the N-cholyl-L-leucine in the solvent of choice, which should be purged with an inert gas. N-Cholyl-L-leucine is soluble (\geq 10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

N-Cholyl-L-leucine is a bile acid amidate.¹ It has been found in fecal samples from extremely premature infants with mild or severe brain injuries. N-Cholyl-L-leucine has been used as a standard for the detection of endogenous bile acid amidates in human plasma and fecal samples.

Reference

1. Pristner, M., Wasinger, D., Seki, D., *et al.* Neuroactive metabolites and bile acids are altered in extremely premature infants with brain injury. *Cell Rep. Med.* **5**(4), 101480 (2024).

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, 11/03/2025

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