

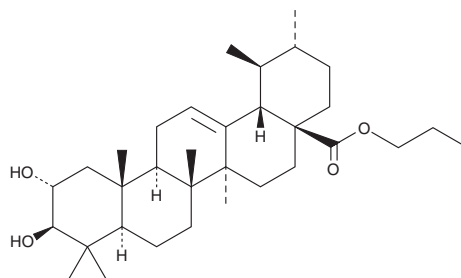
PRODUCT INFORMATION



Corosolic Acid β

Item No. 42738

CAS Registry No.: 958237-73-5
Formal Name: 2 α ,3 β -dihydroxy-urs-12-en-28-oic acid, propyl ester
Synonym: CA β
MF: C₃₃H₅₄O₄
FW: 514.8
Purity: $\geq$ 98%
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: -20°C
Stability: $\geq$ 1 year



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

Description

Corosolic acid β is a synthetic pentacyclic triterpenoid and derivative of corosolic acid (Item No. 27887) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA and siRNA *in vitro* and *in vivo*.¹ It also inhibits glycogen phosphorylase (IC_{50} = 82.5 μ M for the rabbit muscle enzyme).² LNPs containing corosolic acid β and encapsulating an mRNA reporter show greater transfection efficiency than those containing cholesterol (Item Nos. 9003100 | 39088) in 4T1, B16/F10, and HeLa cells.¹ Corosolic acid β -containing LNPs selectively accumulate in tumors and do not accumulate in the liver compared with cholesterol-containing LNPs, which accumulate in both, following intratumoral injection. LNPs containing corosolic acid β and encapsulating siRNA targeting mRNA encoding Bcl-2 decrease tumor volume and lung metastasis and increase survival in a 4T1 murine mammary carcinoma model.

References

1. Liu, Y., Zhang, R., Yang, Y., *et al.* Corosolic acid derivative-based lipid nanoparticles for efficient RNA delivery. *J. Control. Release* **378**, 1-17 (2024).
2. Wen, X., Xia, J., Cheng, K., *et al.* Pentacyclic triterpenes. Part 5: Synthesis and SAR study of corosolic acid derivatives as inhibitors of glycogen phosphorylases. *Bioorg. Med. Chem. Lett.* **17**(21), 5777-5782 (2007).

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, 11/24/2025

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