

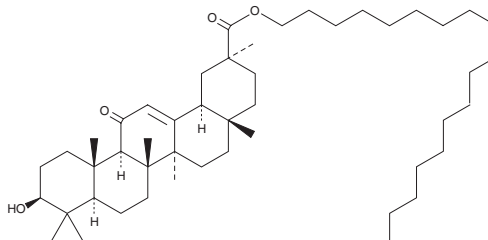
PRODUCT INFORMATION



Stearyl glycyrrhetinate

Item No. 45684

CAS Registry No.: 13832-70-7
Formal Name: (3 β , 20 β)-3-hydroxy-11-oxo-olean-12-en-29-oic acid, octadecyl ester
Synonyms: Glycyrrhetinyl stearate, GR-SA, SG
MF: C₄₈H₈₂O₄
FW: 723.2
Purity: $\geq 95\%$
Supplied as: A 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

Stearyl glycyrrhetinate is supplied as a solid. A stock solution may be made by dissolving the stearyl glycyrrhetinate in the solvent of choice, which should be purged with an inert gas. Stearyl glycyrrhetinate is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Stearyl glycyrrhetinate is a stearyl ester derivative of 18 β -glycyrrhetinic acid (Item No. 11845).¹ It is active against methicillin-susceptible and -resistant *S. aureus* (MRSA; MICs = 64-4,096 and 128-2,048 mg/L, respectively).² Stearyl glycyrrhetinate has been used in the formation of liver-targeting liposomes.^{1,3} Formulations containing stearyl glycyrrhetinate have been used in cosmetics.

References

1. Zhu, J., Zhang, W., Wang, D., *et al.* Preparation and characterization of norcantharidin liposomes modified with stearyl glycyrrhetinate. *Exp. Ther. Med.* **16**(3), 1639-1646 (2018).
2. Oyama, K., Kawada-Matsuo, M., Oogai, Y., *et al.* Antibacterial effects of glycyrrhetinic acid and its derivatives on *Staphylococcus aureus*. *PLoS One* **11**(11), e0165831 (2016).
3. Kim, H., Lee, H.S., Ahn, J.H., *et al.* Targeted nanocomplex delivery for protecting vascular integrity and enhancing anticancer effects in hepatocellular carcinoma. *Small Sci.* **6**(1), e202400616 (2026).

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, 05/22/2026

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