

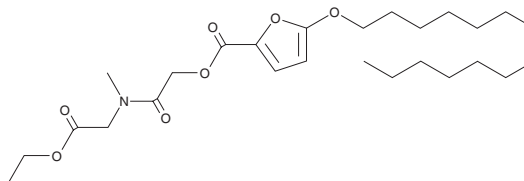
# PRODUCT INFORMATION



## Olumacostat Glasaretil

Item No. 44262

**CAS Registry No.:** 1261491-89-7  
**Formal Name:** 5-(tetradecyloxy)-2-furancarboxylic acid, 2-[(2-ethoxy-2-oxoethyl) methylamino]-2-oxoethyl ester  
**Synonym:** DRM01  
**MF:** C<sub>26</sub>H<sub>43</sub>NO<sub>7</sub>  
**FW:** 481.6  
**Purity:** ≥98%  
**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

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

### Description

Olumacostat glasaretil is an inhibitor of acetyl-CoA carboxylase (ACC).<sup>1</sup> It increases the ratio of acetyl-CoA to free CoA in hamster ear sebaceous glands when applied topically at a concentration of 6% (w/w). Olumacostat glasaretil (0.3, 1, or 3 μM) inhibits triacylglycerol formation induced by T0901317 (Item No. 71810) and insulin in primary human sebocytes. It decreases hamster ear sebaceous gland lipid area when applied topically at a concentration of 6% (w/w) twice per day.

### Reference

1. Hunt, D.W., Winters, G.C., Brownsey, R.W., *et al.* Inhibition of sebum production with the acetyl coenzyme A carboxylase inhibitor olumacostat glasaretil. *J. Invest. Dermatol.* **137(7)**, 1415-1423 (2017).

#### 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, 11/04/2025

#### 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