

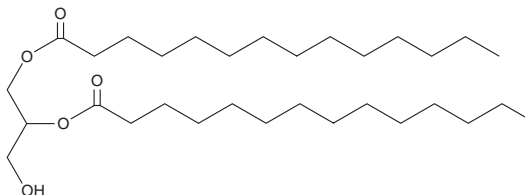
PRODUCT INFORMATION



1,2-Dimyristoyl-*rac*-glycerol

Item No. 28171

CAS Registry No.: 20255-94-1
Formal Name: tetradecanoic acid, 1,1'-[1-(hydroxymethyl)-1,2-ethanediyl] ester
Synonyms: 1,2-Dimyristin, TG(14:0/14:0/0:0)
MF: C₃₁H₆₀O₅
FW: 512.8
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

1,2-Dimyristoyl-*rac*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2-dimyristoyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1,2-Dimyristoyl-*rac*-glycerol is soluble in chloroform.

Description

1,2-Dimyristoyl-*rac*-glycerol is a diacylglycerol that contains myristic acid (Item No. 13351) at the *sn*-1 and *sn*-2 positions. It increases phosphorylation of the insulin receptor in IM-9 cells when used at a concentration of 100 µg/ml.¹ *In vivo*, 1,2-dimyristoyl-*rac*-glycerol (0.5 mg/kg) reduces tumor weight in an LY5178Y mouse xenograft model.²

References

1. Grunberger, G. and Levy, J. Diacylglycerols modulate phosphorylation of the insulin receptor from human mononuclear cells. *Eur. J. Biochem.* **187**(1), 191-198 (1990).
2. Fukusen, N., Kido, H., Kato, Y., *et al.* Diacylglycerols enhance the anti-tumor effect of glucocorticoid on L5178Y lymphoblasts *in vivo*. *Jpn. J. Cancer Res.* **80**(10), 963-967 (1989).

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, 10/31/2022

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