

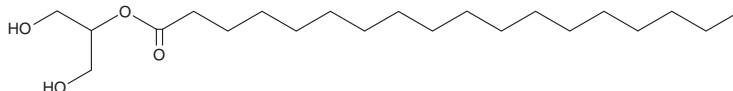
PRODUCT INFORMATION



2-Stearoyl-*rac*-glycerol

Item No. 27282

CAS Registry No.: 621-61-4
Formal Name: octadecanoic acid, 2-hydroxy-1-(hydroxymethyl)ethyl ester
Synonyms: MG(0:0/18:0/0:0), 18:0-MG, 2-Monostearin
MF: C₂₁H₄₂O₄
FW: 358.6
Purity: ≥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

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

Description

2-Stearoyl-*rac*-glycerol is a monoacylglycerol containing stearic acid (Item No. 10011298) at the *sn*-2 position. 2-Stearoyl-*rac*-glycerol levels are increased in mouse plasma following exposure to low-dose ionizing radiation, as well as in rat kidney in a model of ischemia-reperfusion injury.^{1,2}

References

1. Lee, D.Y., Bowen, B.P., Nguyen, D.H., *et al.* Low-dose ionizing radiation-induced blood plasma metabolic response in a diverse genetic mouse population. *Radiat. Res.* **178**(6), 551-555 (2012).
2. Huang, H., van Dulleman, L.F.A., Akhtar, M.Z., *et al.* Proteo-metabolomics reveals compensation between ischemic and non-injured contralateral kidneys after reperfusion. *Sci. Rep.* **8**(1):8539 (2018).

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, 12/22/2022

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