

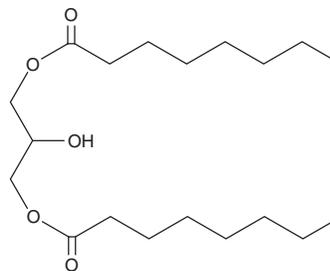
PRODUCT INFORMATION



1,3-Dioctanoyl Glycerol

Item No. 27012

CAS Registry No.: 1429-66-9
Formal Name: octanoic acid, 1,1'-(2-hydroxy-1,3-propanediyl) ester
Synonyms: DG(8:0/0:0/8:0), 1,3-Dioctanoin
MF: C₁₉H₃₆O₅
FW: 344.5
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,3-Dioctanoyl glycerol is supplied as a neat oil. A stock solution may be made by dissolving the 1,3-dioctanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dioctanoyl glycerol is soluble in organic solvents such as ethanol, chloroform, DMSO, and dimethyl formamide (DMF). The solubility of 1,3-dioctanoyl glycerol in ethanol and chloroform is approximately 10 mg/ml. The solubility of 1,3-dioctanoyl glycerol in DMSO and DMF is approximately 1 and 30 mg/ml, respectively.

1,3-Dioctanoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,3-dioctanoyl glycerol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 1,3-Dioctanoyl glycerol has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,3-Dioctanoyl glycerol is a diacylglycerol that contains octanoic acid at the *sn*-1 and *sn*-3 positions.

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, 11/28/2022

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