

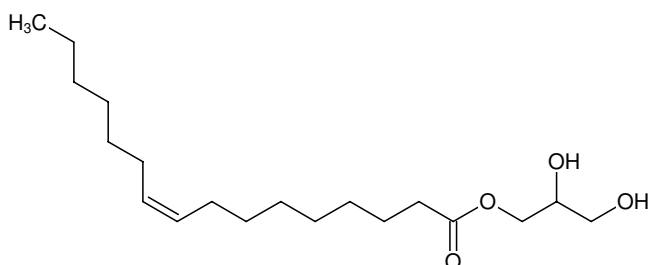


Monopalmitolein

9.7 MAG

1-(9Z-hexadecenoyl)-rac-glycerol

Cat. No.	Amount
X-LCP-102	1 g



Structural formula of Monopalmitolein

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months

Molecular Formula: C₁₉H₃₆O₄

Molecular Weight: 328.43

CAS#: 37515-61-0

Purity: > 99 %

Solubility: Soluble in Hexane

Description:

Host Lipid for protein crystallization in the Lipidic Cubic Phase (LCP)^[1,2]. Used in combination with DSPG to create thermodynamically stable ultraswollen bicontinuous cubic phases with water channels five times larger than traditional lipidic mesophases, suitable for the crystallization of membrane proteins with large extracellular domains^[1].

Once opened, use within 6 months. Purge the vial with nitrogen or argon after usage and store at -20°C.

Related Products:

JBScreen LCP, #CS-340, #CS-213L

LCP Mixer Kit, #X-LCP-M

Monoolein, 9.9 MAG, #X-LCP-101

Monovaccenin, 11.7 MAG, #X-LCP-103

Monoeicosenoin, 11.9 MAG, #X-LCP-104

7.7 MAG, #X-LCP-105

7.8 MAG, #X-LCP-106

7.9 MAG, #X-LCP-107

DSPG, #X-LCP-108

Selected References:

[1] Zabara *et al.* (2018) Design of ultra-swollen lipidic mesophases for the crystallization of membrane proteins with large extracellular domains. *Nat. Commun.* **9**:544.

[2] Caffrey (2015) A comprehensive review of the lipid cubic phase or *in meso* method for crystallizing membrane and soluble proteins and complexes. *Acta Cryst F* **71**:3.