

# Monopalmitolein

Host Lipid for protein crystallization in the Lipidic Cubic Phase (LCP). 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.

## Ordering Information

[Order Online »](#)

JBS-X-LCP-102	1g
---------------	----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name 9.7 MAG, 1-(9Z-hexadecenoyl)-rac-glycerol

CAS 37515-61-0

Formula  $C_{19}H_{36}O_4$

MW 328.43

Purity > 99 %

Solubility Soluble in Hexane

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)