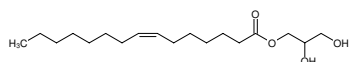


**7.8 MAG**

1-(7Z-pentadecenoyl)-rac-glycerol

Cat. No.	Amount
X-LCP-106	100 mg



Structural formula of 7.8 MAG

**For general laboratory use.****Shipping:** shipped on dry ice**Storage Conditions:** store at -20 °C**Shelf Life:** 6 months**Molecular Formula:** C<sub>18</sub>H<sub>34</sub>O<sub>4</sub>**Molecular Weight:** 314.46**CAS#:** 1309764-71-3**Purity:** > 99 %**Description:**

Host Lipid for protein crystallization in the Lipidic Cubic Phase (LCP) [1-3].

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

**Related Products:**

Monoolein, 9.9 MAG, #X-LCP-101

Monopalmitolein, 9.7 MAG, #X-LCP-102

Monovaccenin, 11.7 MAG, #X-LCP-103

Monoeicosenoin, 11.9 MAG, #X-LCP-104

7.7 MAG, #X-LCP-105

7.9 MAG, #X-LCP-107

**Selected References:**[1] 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.[2] Fowler *et al.* (2015) Gating Topology of the Proton-Coupled Oligopeptide Symporters. *Structure* **23**:290.[3] Li *et al.* (2013) Crystal structure of the integral membrane diacylglycerol kinase. *Nature* **497**:521.