

Anionic phospholipid used in combination with monoacylglycerols (MAGs) 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-108	1g
---------------	----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,2-Distearoyl-sn-glycero-3-phospho-rac-glycerol, sodium salt

CAS 200880-42-8

Formula $C_{42}H_{82}NaO_{10}P$

MW 801.07

Purity > 98 %

Solubility Soluble in Chloroform

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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