

Plasma Membrane Isolation Kit

To isolate intact plasma membrane from mammalian cells.

The Plasma Membrane proteins are a group of specialized proteins involved in transport, cell signaling and cell adhesion, recognition and communication. The low abundance of these proteins makes their study challenging. AkrivisBio's Plasma Membrane Isolation Kit provides optimized reagents in a two phase extraction system for the effective isolation of the plasma membrane from the other cellular membranes based on differential affinity towards polyethylene glycol and dextran. The plasma membrane selectively enters the upper PEG rich phase while the other membranes are found in the lower dextran rich phase.

Ordering Information

[Order Online »](#)

AKR-PI-0104	50preparations
-------------	----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Species independent

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