

# Lipid Extraction Kit

To extract lipids from various sample types.

One of the largest most diverse subset of molecules found in living systems are the lipids. This includes various fats, sterols, mono-, di- and tri-glycerides and phospholipids. While typically viewed as primarily a structural element of the cell (and subcellular) membranes, lipids are also important as a source of energy and in signaling. AkrivisBio's Lipid Extraction Kit provides an easy, efficient method for preparing lipid extracts on a small scale from various sample types. One isolation using 1 mg of sample should recover up to 80  $\mu\text{g}$  of lipid. The assay uses hexane and methanol for extracting almost all lipid types with high efficiency.

## Ordering Information

[Order Online »](#)

AKR-PI-0101	100preparations
-------------	-----------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**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](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)