

Oil Red O Lipid Staining Kit

To qualitatively visualize lipid accumulations in cells grown in cell culture under a microscope.

Cellular lipids are usually found in the membranes of the cell and its organelles. In those cells where lipid is stored for energy production or as source for more complex structures, it is most often seen as droplets consisting primarily of triglycerides although other neutral lipids may be present. Differentiation of adipocytes into mature cells involves a dramatic accumulation of lipid droplets. There are a number of lipid storage diseases (Gaucher's, Niemann-Pick, Fabry's, etc.) where the accumulation of excess lipids can cause severe pathology. Oil Red O is a lipid-soluble azo dye which preferentially stains neutral lipids red so is ideal for visualization of neutral lipid accumulations. The AkrisBio Oil Red O Lipid Staining Kit which also includes hematoxylin to counterstain nuclei blue, is an easy, inexpensive way to qualitatively visualize lipid accumulations in cells grown in cell culture under a microscope. The kit provides enough materials to stain two plates (6- 12-, 24- or 96-well) or five 100 mm culture dishes.

Ordering Information

[Order Online »](#)

AKR-COS-0101	1Kit
--------------	------

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

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