

# Lipid Peroxidation Assay

A plate-based colorimetric assay to measure malondialdehyde (MDA) levels in samples.

Lipids play an important role in human homeostasis. They serve as structural components of cell membranes, organelles and tissues, function as energy storehouses and also act as signaling molecules. Lipids are sensitive to oxidation by reactive oxygen species. Lipid peroxidation is a metabolic process that causes degradation of various lipid molecules by reactive oxygen species thereby resulting in cell damage and eventually, cell death. Lipid peroxidation occurs in both animals and plants and also in vivo during aging and in certain disease states including atherosclerosis, heart failure, Alzheimer's disease, rheumatic arthritis, cancer, and other immunological disorders. Malondialdehyde (MDA) is one of the most known secondary products of lipid peroxidation, and is used as a useful biomarker for the assessment of lipid peroxidation. AkrivisBio's Lipid Peroxidation Assay provides a simple and convenient method for the detection of MDA through its reaction with Thiobarbituric Acid to generate a chromophore. The chromophore is then quantified colorimetrically at 532 nm or fluorometrically at Ex/Em = 532/553 nm.

## Ordering Information

[Order Online »](#)

AKR-MA-0102	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Colorimetric detection

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)