

# Free Fatty Acid Assay

A plate based colorimetric/fluorometric assay to measure the concentration of free fatty acids in samples.

Fatty acids are an important source of energy, as structural components and as precursors to a variety of bioactive chemicals collectively referred to as oxylipins with a diverse set of biological actions, including pro- and anti-inflammatory, vasoactive, autocrine and paracrine functions. High levels of free fatty acids have been associated with conditions including obesity, diabetes, liver diseases etc. AkrivisBio's Fatty Acid Assay is a simple and sensitive assay to measure free fatty acids in various sample types. The assay is based upon the conversion of fatty acids to CoA derivatives, which are subsequently oxidized to form color/fluorescence that is quantified at OD 570 nm or at Ex/Em = 535/587 nm respectively. The intensity of color/fluorescence is proportional to the concentration of free fatty acid in samples.

## Ordering Information

[Order Online »](#)

AKR-MA-0111	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°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)