

# Triglyceride Assay

**A plate based colorimetric/fluorometric assay to measure 0-10 nanomoles/well (colorimetric) and 1-5 picomoles/well (fluorometric) of Triglyceride levels in samples.**

Triglycerides are the primary constituent of body fat in both animals and plants and are an important energy source. Triglyceride Assay is used to investigate lipid-related disorders in research studies. AkrivisBio's Triglyceride Assay is a simple and sensitive assay to measure triglyceride levels in samples. In this assay, mono, di and triglycerides in samples are hydrolyzed by lipase to generate free fatty acids and free glycerol. Glycerol is then oxidized to generate a product that reacts with the probe to form an intense color or fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

## Ordering Information

[Order Online »](#)

AKR-MA-0110	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)