

# Glycogen Assay

A plate based colorimetric/fluorometric assay to quantify the amount of glycogen present in samples.

Glycogen, the main form of glucose storage in animals is synthesized and stored primarily in liver and muscle. Glycogen is formed on a core of glycogenin which links the first few molecules of glucose after which other enzymes continue adding glucose monomers to the growing polymer with  $\alpha$ -1,4 linkages extending linear chains, and  $\alpha$ -1,6 linkages giving a branched structure. A number of disease states are related to the inability to store or utilize stored glycogen properly. AkrivisBio's Glycogen Assay is a simple sensitive method to measure glycogen levels in biological samples. In the assay, glycogen is hydrolyzed, and the glucose formed is oxidized with the formation of H<sub>2</sub>O<sub>2</sub> which is utilized by peroxidase to oxidize ADHP resulting in the formation of intense color ( $\lambda_{\text{max}}$  570 nm) and fluorescence (Excitation/Emission = 535/587 nm). The assay can determine glycogen content from the low nanogram to high microgram level.

## Ordering Information

[Order Online »](#)

AKR-MA-0112	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)