

Glycogen Assay II

A plate based colorimetric assay to measure glycogen levels in biological samples.

Glycogen is a polysaccharide that serves as a storage form of glucose in cells, particularly in liver and muscle tissues. Metabolic abnormalities of glycogen metabolism, referred to as glycogen storage diseases, include Von Gierke disease, Pompe disease and Cori disease. AkrisBio's Glycogen Assay II provides a simple, sensitive method of measuring glycogen levels in biological materials. In the assay, glycogen is hydrolyzed to glucose, which is oxidized to form NADH. The NADH formed then reduces a nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm. The Glycogen Assay II is valuable in various research areas, including glycogen metabolism, energy balance, and glycogen-related disorders.

Ordering Information

[Order Online »](#)

AKR-MA-0136	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

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