

Glucose Assay

A plate based colorimetric/fluorometric assay to measure glucose concentration in a wide range of biological samples.

Glucose is the primary source of energy in all organisms. In plants, glucose is stored as starch and as glycogen in animals. Blood glucose concentration is highly regulated and a variety of metabolic disorders is reflected in a significant deviation in glucose concentration.

AkrivisBio's Glucose Assay is a simple, sensitive way to measure Glucose in a wide range of biological samples. In this assay, Glucose is oxidized by glucose oxidase forming gluconolactone and hydrogen peroxide. Hydrogen peroxide is utilized by the peroxidase to oxidize ADHP thereby resulting in the formation of intense color or fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

Ordering Information

[Order Online »](#)

AKR-MA-0122	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