

Glutathione Peroxidase Activity Assay

A plate based colorimetric assay to measure Glutathione peroxidase (GPx) activity in a

variety of samples with a sensitivity below 0.2 mU per well. Glutathione peroxidase (GPx) is an essential enzyme involved in glutathione metabolism, that helps maintain redox balance and protects against oxidative stress. GPx catalyzes the reduction of H₂O₂ and organic hydroperoxides, using glutathione as a co-substrate preventing accumulation of harmful reactive oxygen species within cells, which cause damage to proteins, lipids, and DNA. GPx is widely distributed throughout the body, particularly abundant in organs with high metabolic activity, (liver, kidney, lung, and red blood cells). Expression levels are modulated in response to environmental factors, including dietary selenium and oxidative stress levels. GPx deficiency is associated with pathological conditions, including neurodegenerative disease, cardiovascular disorders, cancer, and age-related macular degeneration. AkrivisBio's Glutathione Peroxidase Activity Assay is a simple, sensitive means of assessing GPx activity in a variety of samples with a sensitivity below 0.2 mU per well.

Ordering Information

[Order Online »](#)

AKR-MA-0137	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