

GAPDH Activity Assay

A plate based colorimetric assay to measure GAPDH activity in a variety of samples with a sensitivity into the low $\mu\text{U/ml}$.

GAPDH catalyzes the conversion of glyceraldehyde-3-phosphate to 1, 3-bisphosphoglycerate (1,3-BPG), creating a high-energy phosphate bond while simultaneously facilitating the formation of NADH. This is a key step in glycolysis. 1,3-BPG goes on to result in ATP production while the NADH powers the electron transport chain, synthesizing more ATP. GAPDH is an essential constitutive enzyme whose activity is regulated by several factors, including cellular energy demand, redox state and post-translational modifications. Dysregulation of GAPDH is associated with cancer and neurodegenerative disorders. AkrisBio's GAPDH Assay is a simple, sensitive means of measuring GAPDH activity in a variety of samples with a sensitivity into the low $\mu\text{U/ml}$.

Ordering Information

[Order Online »](#)

AKR-MA-0153	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