

Pyruvate Kinase Activity Assay

A plate-based colorimetric/fluorometric assay to measure pyruvate kinase activity in samples below 50 μ U.

Pyruvate kinase (PK) is a key enzyme in the glycolytic pathway catalyzing the conversion of phosphoenolpyruvate to pyruvate forming ATP. This reaction serves as a key control step regulating the overall rate of glycolysis and determining the fate of glucose metabolism. Pyruvate kinase plays a central role providing energy for cellular processes, as pyruvate is a metabolite common to several pathways. Pyruvate assay helps in understanding glycolytic flux, assessing enzyme regulation, and evaluating metabolic alterations in diseases including cancer, diabetes, or genetic disorders. AkrivisBio's Pyruvate Kinase Activity Assay is a simple, sensitive assay for determining PK activity in a variety of sample types below 50 μ U. In the assay, pyruvate kinase converts Phosphoenolpyruvate and ADP to generate pyruvate and ATP. Pyruvate is then oxidized by pyruvate oxidase, forming hydrogen peroxide. Hydrogen peroxide is utilized by peroxidase to oxidize ADHP resulting in the formation of intense color/fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

Ordering Information

[Order Online »](#)

AKR-MA-0107	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