

Phosphofructokinase Activity Assay

A plate based colorimetric assay to quantify Phosphofructokinase (PFK) activity in a variety of biological samples with a sensitivity in the microunit range.

Phosphofructokinase (PFK) is a key glycolytic enzyme and plays a major regulatory role in glycolysis. PFK catalyzes the conversion of fructose-6-phosphate to fructose-1,6-bisphosphate. It is allosterically regulated by several chemicals including ATP, ADP and citrate. When energy levels are low the ADP to ATP ratio is high which turns on PFK, accelerating glycolysis. Conversely when energy levels are high ATP and citrate work to turn off PFK. Dysregulation of PFK leads to metabolic disorders. PFK has been targeted in cancer therapy since cancer cells display enhanced glycolytic activity. AkrivisBio's Phosphofructokinase Activity Assay is a simple sensitive means of quantifying PFK activity in a variety of biological samples with a sensitivity in the microunit range.

Ordering Information

[Order Online »](#)

AKR-MA-0165	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