

# ATP Assay

A plate based colorimetric/fluorometric assay to quantify the amount of ATP present in various samples. The assay can detect down to 50 pmol (1  $\mu$ M) of ATP.

ATP is the primary energy currency of cells. Virtually all energy dependent processes including metabolism, muscle contraction, signal transduction etc. utilize the chemical energy stored in the phosphate bond of ATP. AkrisBio's ATP assay kit is simple, robust assay to quantify the amount of ATP present in various samples. In this assay, ATP from samples phosphorylate glycerol. Glycerol phosphate is then oxidized with the formation of hydrogen peroxide, which is then utilized by the peroxidase to oxidize ADHP resulting in the formation of intense color/fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm.

## Ordering Information

[Order Online »](#)

AKR-MA-0113	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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)