

Phosphate Assay

A plate-based colorimetric assay to measure as low as 1 μM of Phosphate levels in samples.

Phosphate is found virtually everywhere in the biological system. It is used as a building block for several crucial substances including those used by the cell for energy, cell membranes, and DNA. There are over 500 kinases and 200 phosphatases in the human genome alone, which function by adding or removing phosphate groups to/from hundreds or thousands of molecules including proteins, sugars, nucleic acids, lipids and a plethora of other small molecules. AkrivisBio's Phosphate Assay provides a simple, quick and precise method for measuring phosphate levels in various samples. In this assay, phosphate ions in samples react with malachite green and ammonium molybdate to form a chromogenic complex measured at OD 650 nm. The intensity of the color of the chromogenic complex is directly proportional to the phosphate concentration in the sample.

Ordering Information

[Order Online »](#)

AKR-MA-0105	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

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