

Alkaline Phosphatase Activity Assay

A plate based colorimetric assay to measure the presence and activity of alkaline phosphatase enzyme in various sample types.

Alkaline phosphatase (ALP) is an enzyme that catalyzes the hydrolysis of phosphate esters under alkaline conditions with the production of free phosphate. AkrivisBio's ALP Activity Assay is a simple, sensitive assay to measure ALP activity in a variety of biological samples. In this assay the specific alkaline phosphatase substrate, p-nitrophenyl phosphate is hydrolyzed to phosphate ion and p-nitrophenol. P-nitrophenol formation is then monitored using a spectrophotometer at λ_{\max} 405 nm. The absorbance or color intensity is directly proportional to the alkaline phosphatase activity in the sample.

Ordering Information

[Order Online »](#)

AKR-MA-0109	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°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