

Lactate Assay

A plate based colorimetric/fluorometric assay to measure L-lactate in a variety of biological samples.

Lactate (L-lactate) is the primary end product of anaerobic metabolism. It accumulates when its consumption via the respiratory chain is slower than glycolysis, driven by the immediate need for energy. The Lactate pool can either be oxidized to pyruvate and further processed via normal mitochondrial oxidative phosphorylation generating ATP or it can be converted back to glucose via gluconeogenesis. In addition to being a metabolic fuel, lactate has been discovered to be a metabolic feedback regulator and a signaling molecule. Abnormally high lactate levels can be caused either by enhanced production or impaired utilization in pathological conditions thereby providing valuable information about cellular metabolism or diseased conditions. AkrivisBio's Lactate Assay is a simple, specific and sensitive method to measure L-lactate levels in a variety of biological samples. In this assay, lactate is oxidized by L-lactate oxidase thereby forming pyruvate and hydrogen peroxide. Hydrogen peroxide is then utilized by the peroxidase to oxidize ADHP thereby resulting in the formation of intense color or fluorescence measured at OD 570 nm or at Ex/Em = 535/587 nm respectively.

Ordering Information

[Order Online »](#)

AKR-MA-0115	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