

L-Lactate Assay II

A plate based colorimetric assay to measure L(+)-Lactate concentration in various samples.

L(+)-Lactate is formed in animals by the action of the enzyme Lactate Dehydrogenase. L-Lactate is the major lactate stereoisomer formed in human metabolism and D-Lactate is present at much lower levels (~1-5% of L-Lactate). L-Lactate plays an important role in many biological processes. High lactate concentrations are associated with pathological conditions such as diabetes and lactic acidosis. AkrisBio's L-Lactate Assay is a simple, colorimetric assay to measure Lactate in various biological samples. In this assay, L-lactate is oxidized by L-Lactate Dehydrogenase to form pyruvate and NADH. NADH then reduces WST, a readily reducible nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm. This assay is useful for detecting L(+)-Lactate at concentrations of 0.05 to 20 mM. There is no need for pretreatment or purification of samples.

Ordering Information

[Order Online »](#)

AKR-MA-0123	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