

LDH Activity Assay

A plate based colorimetric/fluorometric assay to quantify Alanine aminotransferase activity in a variety of biological samples with a detection limit below 10 mU per well.

Lactate Dehydrogenase (LDH) is an enzyme found in various tissues and is involved in the conversion of lactate to pyruvate and vice versa. LDH is commonly used to assess cell viability, disease progression etc. in clinical and research settings. AkrivisBio's LDH Activity Assay is a simple, and sensitive assay to measure LDH activity in a wide variety of samples with a sensitivity to accurately quantify as little as 0.02 mU of LDH activity. In this assay, L-lactate is converted to pyruvate in the presence of LDH with the formation of NADH. NADH then reduces WST, a readily reducible nearly colorless tetrazolium to a highly colored formazan measured at OD 450 nm.

Ordering Information

[Order Online »](#)

AKR-MA-0117	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)