

# LDH Based Cell Cytotoxicity Assay

For the detection of LDH released from damaged cells.

Cell death or cytotoxicity is classically evaluated by assessing the degree of plasma membrane damage. Lactate dehydrogenase (LDH) is an intracellular enzyme that is released into the extracellular medium upon cellular damage or lysis, making it a useful indicator of cytotoxic effects. AkrivisBio's LDH Cytotoxicity Assay Kit is used to assess cell membrane integrity and measure cell death or cytotoxicity. The assay utilizes WST, a tetrazolium detection reagent to give a fast, highly sensitive measure of LDH release from damaged cells. The intensity of color correlates directly with the number of damaged cells.

## Ordering Information

[Order Online »](#)

AKR-CPT-0101	500Reactions
--------------	--------------

## 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](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)