

LDH Cytotoxicity WST Assay

Colorimetric kit for the determination of cytotoxicity through the measurement of lactate dehydrogenase activity

The LDH cytotoxicity WST assay is a colorimetric assay kit used to determine cytotoxicity by measuring lactate dehydrogenase activity released from damaged cells.

LDH is a stable cytoplasmic enzyme presented in all types of cells and released into the cell culture medium through damaged plasma membrane.

Citations: 7

[View Online »](#)

Ordering Information

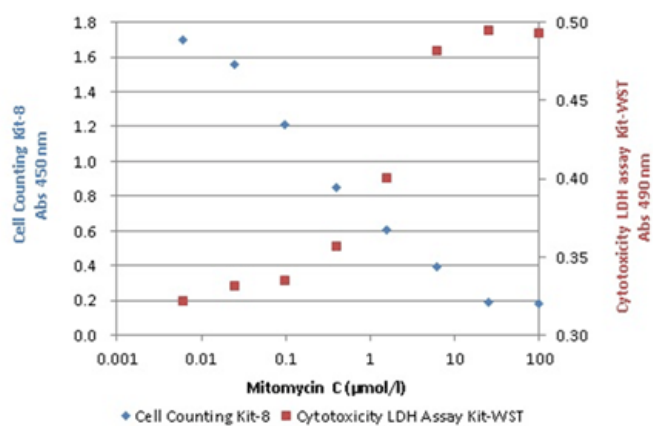
[Order Online »](#)

ENZ-KIT157-0500	500 tests
ENZ-KIT157-2500	5x500 tests

Manuals, SDS & CofA

[View Online »](#)

- Homogenous assay using living cells or non-homogenous assay protocol using cell culture supernatant
- Stable working solution



Difference between the Cell Counting Kit-8 (Prod. No. ALX-850-039) and the LDH Cytotoxicity WST Assay. The Cell Counting Kit-8 determines the number of viable cells in cell proliferation and cytotoxicity assays. On the other hand, the LDH Cytotoxicity WST Assay evaluates the cytotoxic level by measuring the amount of LDH released from damaged cells. Assessing the level of cytotoxicity by combining both the Cell Counting Kit-8 Kit and the LDH Cytotoxicity WST Assay can provide researchers with a complete evaluation of cell health.



Mechanism for LDH Cytotoxicity WST Assay

Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Colorimetric detection
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Quantity	500 tests/5ml 2500 tests/5 x 5ml
Technical Info / Product Notes	<p>Principle: LDH catalyzes dehydrogenation of lactate to pyruvate, reducing NAD to NADH. NADH reduces a water soluble tetrazolium salt (WST) in the presence of an electron mediator to produce an orange formazan dye. The amount of formazan dye produced is proportional to the amount of LDH released into the medium, indicating cytotoxicity.</p>

Last modified: May 29, 2024



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)