

Cytotoxicity

LDH Assay Kit-WST

FEATURES

- ◇ Colorimetric Assay System
- ◇ Utilize for both Homo & Non-homogeneous Assays
- ◇ Very Stable Reagent



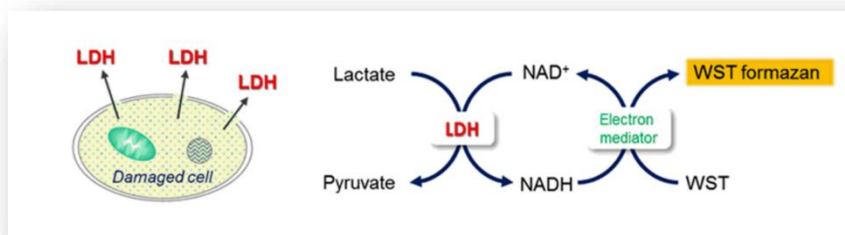
CK12-01 100tests
 CK12-05 500tests
 CK12-20 2000tests



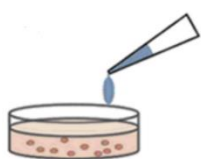
PRODUCT DESCRIPTION

Cytotoxicity LDH Assay Kit-WST is a kit for determination of cytotoxicity by measuring a lactate dehydrogenase (LDH) 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. Since Cytotoxicity LDH Assay kit-WST neither reflects the activity of living cells nor is harmful to cells, cytotoxicity can be measured with living cells (homogeneous assay). In addition, non-homogeneous assay that is performed by using the cell culture supernatant is also possible. Unlike competitive products, the reconstituted Working Solution is stable under refrigerated condition and the Working Solution can be used for long periods as ready-to-use solution after the preparation.

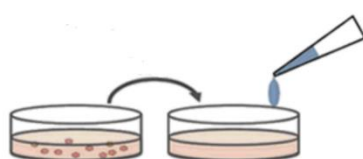
ASSAY PRINCIPLE



NO NEED TO CHOOSE THE TYPE OF A LDH ASSAY KIT



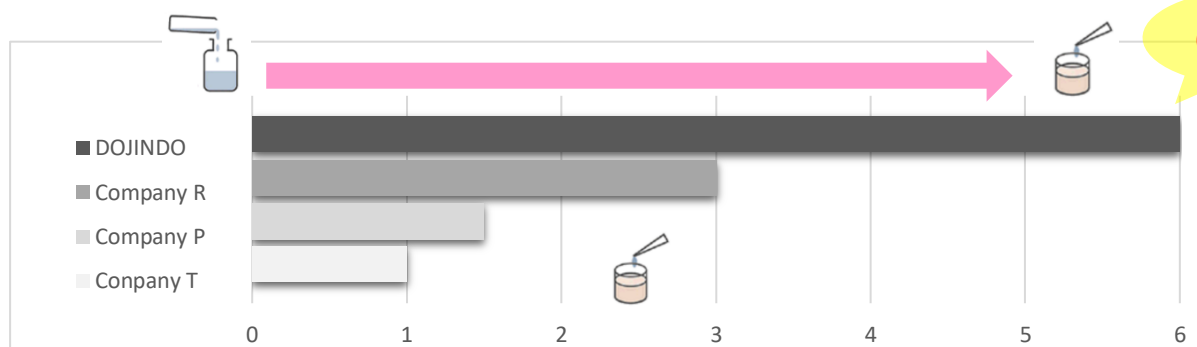
Homogeneous Assay



Non-homogeneous Assay

Whereas other LDH kit suppliers offer individual kit for either homogeneous or non-homogeneous assay, Dojindo' LDH kit allows both assay by using a single product.

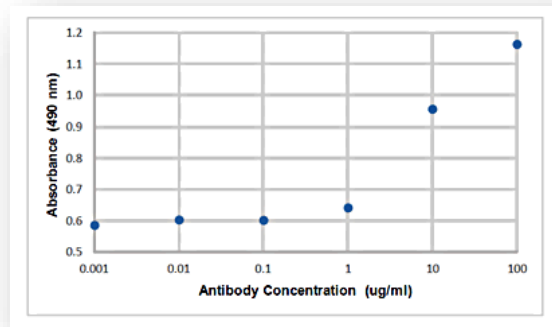
LONGEST STABILITY OF REAGENT AMONG COMMERCIALY AVAILABLE



STATE OF THE ART LIFE SCIENCE TECHNOLOGIES

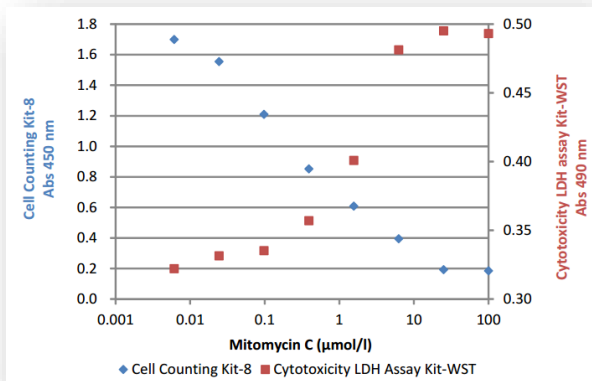
FOR CDC (Complement-Dependent Cytotoxicity) ASSAY

DOJIND's LDH Assay Kit-WST can also be used for CDC assay. Please contact with our technical support at info@dojindo.eu.com to get your assay protocol.



RELATED PRODUCT

Recent research trend may require determination of cell viability by different approaches. Dojindo would like to propose Cell Counting Kit-8 as a suitable co-parameter with LDH released assay.



CELL COUNTING KIT-8

Unit sizes;
CK04-11 1000tests
CK04-13 3000tests
CK04-20 10000tests

Cell Counting Kit-8 is the most well accepted colorimetric cell viability/proliferation assay kit by academia. For more detailed info, please visit us online at www.dojindo.eu.com with a searching keyword, CCK-8.

REFERENCES

- 1) S. Watanabe, C. S. Moniaga, S. Nielsen, M. Hara-Chikuma, "Aquaporin-9 facilitates membrane transport of hydrogen peroxide in mammalian cells.", *Biochem Biophys Res Commun.*, **2016**, 471, 191.
- 2) T. Fukami, A. Iida, K. Konishi, M. Nakajima, "Human arylacetamide deacetylase hydrolyzes ketoconazole to trigger hepatocellular toxicity", *Biochem. Pharmacol.*, **2016**, doi:10.1016/j.jphs.2016.07.007.
- 3) Y. Takanezawa, R. Nakamura, Y. Sone, S. Uruguchi and M. Kiyono, "Atg5-dependent autophagy plays a protective role against methylmercury-induced cytotoxicity", *Toxicol. Lett.*, **2016**, 262, 135.
- 4) S. Wakatsuki, A. Furuno, M. Ohshima, and T. Araki, "Oxidative stress-dependent phosphorylation activates ZNRF1 to induce neuronal/axonal degeneration", *J Cell Biol.*, **2015**, 211, (4), 881.
- 5) A. Ishijima, K. Minamihata, S. Yamaguchi, S. Yamahira, R. Ichikawa, E. Kobayashi, M. Iijima, Y. Shibasaki, T. Azuma, T. Nagamune & I. Sakuma, "Selective intracellular vaporisation of antibody-conjugated phase-change nano-droplets in vitro", *Nature.*, **2017**, DOI: 10.1038/srep44077.



EUROPEAN HEADQUARTERS

DOJINDO EU GMBH
Leopoldstr. 254, 80807 Munich, Germany
Phone +49 89 3540-4806
Fax +49 89 3540-4806
email info@dojindo.eu.com
www.dojindo.eu.com

DISTRIBUTED BY