
Product Manual

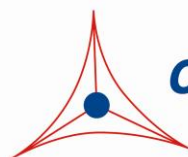
CytoSelect™ LDH Cytotoxicity Assay Kit

Catalog Number

CBA-241

960 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



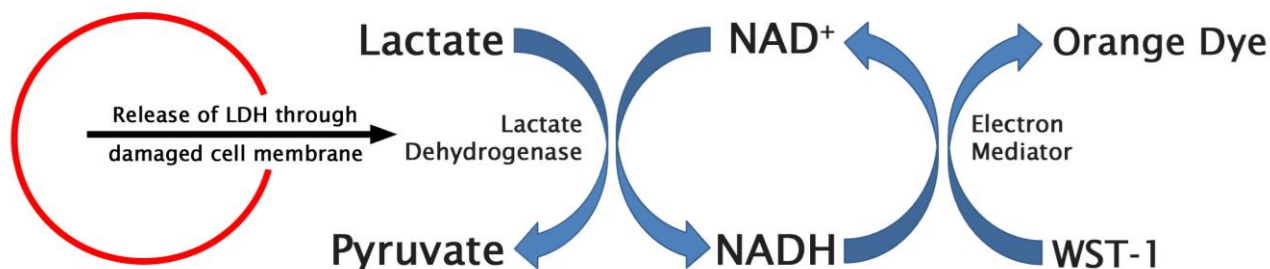
CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

The measurement and monitoring of cell cytotoxicity is an essential technique in any laboratory focused on cell-based research. This skill allows for the optimization of cell culture conditions. More importantly, the cytotoxic nature of anticancer compounds in toxicology testing, the toxicity of therapeutic chemicals in drug screening, and cell-mediated cytotoxicity can all be assessed through this assay-based approach.

Cell cytotoxicity characteristics include loss of cellular metabolic activity or cell membrane integrity. One method for measuring metabolic activity is to incubate the cells with a tetrazolium salt such as MTT, which is cleaved into a colored formazan product by metabolically active cells. Similarly, the green fluorescent dye Calcein AM can measure intracellular esterase activity in proliferating live cells, while dyes such as trypan blue or propidium iodide can enter and stain cells that have lost membrane integrity.

Cell Biolabs' CytoSelect™ LDH Cytotoxicity Assay Kit provides a colorimetric format for measuring and monitoring cell cytotoxicity. The kit contains sufficient reagents for the evaluation of 960 assays in 96-well plates. Cells can be plated and then treated with compounds or agents that affect cell viability. Upon cell death, lactate dehydrogenase (LDH), a soluble enzyme found in the cytoplasm, is released into the growth media. The growth media is then transferred to another plate and the released LDH is then detected with a cytotoxicity reagent. In the presence of lactate substrate (included in the LDH Cytotoxicity Reagent) LDH converts lactate to pyruvate and generates nicotinamide adenine dinucleotide (NADH). The WST-1 molecule, also present in the LDH Cytotoxicity Reagent, is converted from WST-1 to the orange formazan form. An increase in cell cytotoxicity is accompanied by increased LDH release and increased colorimetric signal. The assay principles are basic and can be applied to most eukaryotic cell lines, including adherent and non-adherent cells and certain tissues, depending on LDH expression levels. The LDH Cytotoxicity Reagent can be used to detect cytotoxicity in mammalian cells.



Related Products

1. CBA-232: Quantitative Cellular Senescence Assay (SA β -Gal)
2. CBA-240: Cell Viability and Cytotoxicity Assay
3. CBA-251: CytoSelect™ BrdU Cell Proliferation ELISA Kit
4. CBA-252: CytoSelect™ MTT Cell Proliferation Assay
5. CBA-253: CytoSelect™ WST-1 Cell Proliferation Assay Reagent

Kit Components (shipped on blue ice)

1. LDH Cytotoxicity Assay Reagent (Part No. 124101): One 10 mL bottle of reagent
2. Triton X-100 Solution (Part No. 124102): One 10 mL bottle of 10% Triton X-100

Materials Not Supplied

1. Cells for measuring cytotoxicity
2. Cytotoxicity mediating compound to be tested
3. Cell culture medium
4. 96-well clear cell culture plates
5. Microtiter plate reader capable measuring absorbance at 450 nm

Storage

The LDH Cytotoxicity Assay Reagent is a clear, slightly red, ready-to-use solution. Aliquot as needed to avoid repeated freeze-thaw cycles and store at -20°C protected from light. If precipitates or turbidity are observed upon thawing, warm the solution to 37°C for 5-10 minutes and agitate to dissolve the precipitates.

Store the Triton X-100 Solution at Room Temperature.

Assay Protocol

Note: Both a negative and a positive control should be run alongside experimental samples. Each experimental sample and control should be assayed in duplicate.

1. Prepare a cell suspension containing $0.1\text{--}1.0 \times 10^6$ cells/mL in medium.
2. Add the cell suspension to a 96-well cell culture plate for each experimental sample, negative and positive control well according to Table 1 below. Include the compound to be tested in experimental wells only.

	Experimental Sample	Negative Control	Positive Control
Cell Suspension	150 μ L	150 μ L	150 μ L
Cytotoxicity-mediating compound to be tested	+	-	-

Table 1. Preparation of Experimental and Control Wells.

3. Culture the cells for 24-96 hours at 37°C and 5% CO₂ in a humidified incubator.
4. Add sterile water or the provided Triton X-100 Solution to each well according to Table 2 below.

	Experimental Sample	Negative Control	Positive Control
Sterile Water	15 μ L	15 μ L	-
Triton X-100 Solution	-	-	15 μ L

Table 2. Lysis of Positive Control Cells.

- Incubate 5-10 minutes at room temperature.
- Transfer 90 μ L of medium from each well to a clean 96-well plate suitable for a plate reader.
- Add 10 μ L of LDH Cytotoxicity Assay Reagent.
- Incubate plate at 37°C and 5% CO₂ for 0.5-4 hours.
- Read OD using 450 nm as the primary wave length.

Example of Results

The following figure demonstrates typical results with the LDH Cytotoxicity Assay Kit. One should use the data below for reference only. This data should not be used to interpret actual results.

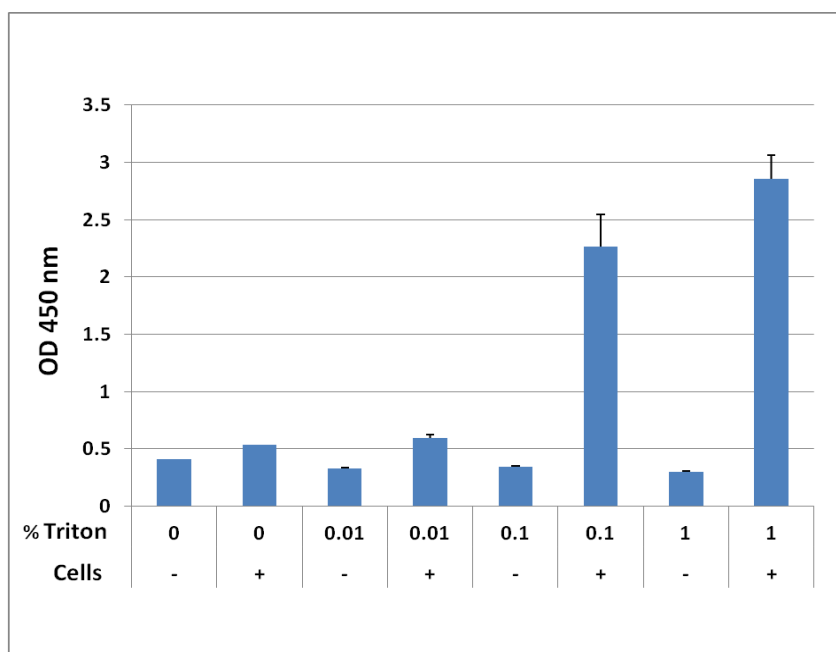


Figure 1. LDH Release from Human HEK 293 Cells. HEK 293 cells were seeded at 20000 cells per well and allowed to grow for 24 hours. After adding various concentrations of Triton X-100, LDH Cytotoxicity Assay Reagent was added. Cells were incubated for 30 minutes at 37°C and 5% CO₂, and then read on a colorimetric plate reader.

Calculation of Results

Positive control wells represent maximal LDH release, while negative control wells represent background LDH release. The OD for negative controls is subtracted from both experimental and positive control OD values, and results are reported as a relative cytotoxicity percentage:

OD experimental sample — OD negative control

OD positive control — OD negative control

x 100 = % Relative Cytotoxicity

References

1. Jacobsen MD, Weil M, Raff MC. (1996) *J Cell Biol* **133**, 1041.
2. Liu X, Cui X, Shan N, Li Y, Fang X, Ding M, Wang X. (2013) *Cancer Cell Int.* **13**, 112
3. Yamaori S, Ishii H, Chiba K, Yamamoto I, Watanabe K (2013) *Toxicology* **314**, 251
4. Wang Y, Qu L, Gong L, Sun L, Gong R, Si J (2013) *Cancer Biother Radiopharm.* **28**, 623
5. Xue TY, He YZ, Zhang JJ, Zhang XY, Xin YJ, Lu Y, Li YH (2013) *Clin Lab* **59**, 613

Recent Product Citations

1. Dutta, B. et al. (2023). Plant-derived cell-penetrating microprotein α -astratide aM1 targets Akt signaling and alleviates insulin resistance. *Cell Mol Life Sci.* **80**(10):293. doi: 10.1007/s00018-023-04937-y.
2. Konaka, H. et al. (2023). Secretion of mitochondrial DNA via exosomes promotes inflammation in Behçet's syndrome. *EMBO J.* doi: 10.15252/emj.2022112573.
3. Haba, D. et al. (2023). Investigation of ultrasound and low-frequency vibrations for glycometabolism promotion in 3T3-L1 adipocytes. *Japanese Journal of Electrophysical Agents.* doi: 10.57337/jjeapt.2023-001.
4. Eller, K.A. et al. (2023). Photoactivated antibiotics to treat intracellular infection of bacteria. *Nanoscale Adv.* **5**(7):1910-1918. doi: 10.1039/d2na00378c.
5. Yu, Z. et al. (2022). UVA induces retinal photoreceptor cell death via receptor interacting protein 3 kinase mediated necroptosis. *Cell Death Discov.* **8**(1):489. doi: 10.1038/s41420-022-01273-1.
6. Faist, A. et al. (2022). Inhibition of p38 signaling curtails the SARS-CoV-2 induced inflammatory response but retains the IFN-dependent antiviral defense of the lung epithelial barrier. *Antiviral Res.* doi: 10.1016/j.antiviral.2022.105475.
7. Tanas, A. et al. (2022). In Vitro and In Vivo Neuroprotective Effects of Sarcosine. *Biomed Res Int.* doi: 10.1155/2022/5467498.
8. Pardo-Mora, D.P. et al. (2022). Microarray analysis of canine osteosarcoma cells exposed to Colombian propolis. *Phytomedicine Plus.* doi: 10.1016/j.phyplu.2022.100356.
9. Tang, Y. et al. (2022). Cross-seeding between A β and SEVI indicates a pathogenic link and gender difference between alzheimer diseases and AIDS. *Commun Biol.* **5**(1):417. doi: 10.1038/s42003-022-03343-7.
10. Schreiber, A. et al. (2022). The MEK1/2-inhibitor ATR-002 efficiently blocks SARS-CoV-2 propagation and alleviates pro-inflammatory cytokine/chemokine responses. *Cell Mol Life Sci.* **79**(1):65. doi: 10.1007/s00018-021-04085-1.
11. Griffiths, A. et al. (2021). Nicotinamide N-methyltransferase upregulation via the mTORC1-ATF4 pathway activation contributes to palmitate-induced lipotoxicity in hepatocytes. *Am J Physiol Cell Physiol.* **321**(3):C585-C595. doi: 10.1152/ajpcell.00195.2021.
12. Hiu, J.J. & Yap, M.K.K. (2021). The effects of Naja sumatrana venom cytotoxin, sumaCTX on alteration of the secretome in MCF-7 breast cancer cells following membrane permeabilization. *Int J Biol Macromol.* **184**:776-786. doi: 10.1016/j.ijbiomac.2021.06.145.

13. Yu, Z. et al. (2021). Receptor interacting protein 3 kinase, not 1 kinase, through MLKL-mediated necroptosis is involved in UVA-induced corneal endothelium cell death. *Cell Death Discov.* **7**(1):366. doi: 10.1038/s41420-021-00757-w.
14. Haba, D. et al. (2021). Low-frequency vibration promotes AMPK-mediated glucose uptake in 3T3-L1 adipocytes. *Heliyon.* **7**(9):e07897. doi: 10.1016/j.heliyon.2021.e07897.
15. Torrisi, F. et al. (2021). Connexin 43 and Sonic Hedgehog Pathway Interplay in Glioblastoma Cell Proliferation and Migration. *Biology (Basel).* **10**(8):767. doi: 10.3390/biology10080767.
16. Dziwornu, G.A. et al. (2021). Antimalarial Benzimidazole Derivatives Incorporating Phenolic Mannich Base Side Chains Inhibit Microtubule and Hemozoin Formation: Structure-Activity Relationship and In Vivo Oral Efficacy Studies. *J Med Chem.* doi: 10.1021/acs.jmedchem.1c00354.
17. Tang, Y. et al. (2021). Repurposing a Cardiovascular Disease Drug of Cloridarol as hIAPP Inhibitor. *ACS Chem Neurosci.* doi: 10.1021/acscchemneuro.1c00091.
18. Pardo-Mora, D.P. et al. (2021). Apoptosis-related gene expression induced by Colombian propolis samples in canine osteosarcoma cell line. *Vet World.* **14**(4): 964-971. doi: 10.14202/vetworld.2021.964-971.
19. Aydemir Sezer, U. et al. (2020). Zero Valent Iron Nanoparticles Containing Nanofiber Scaffolds for Nerve Tissue Engineering. *J Tissue Eng Regen Med.* doi: 10.1002/term.3137.
20. Probst, A. et al. (2020). Expanding the activity profile of pyrido[1,2-a]benzimidazoles: Synthesis and evaluation of novel N1-1-phenylethanamine derivatives against *Schistosoma mansoni*. *ACS Infect Dis.* doi: 10.1021/acsinfecdis.0c00278.
21. Shun, Q.T. et al. (2020). Naja sumatrana venom cytotoxin, sumaCTX exhibits concentration-dependent cytotoxicity via caspase-activated mitochondrial-mediated apoptosis without transitioning to necrosis. *Toxin Rev.* doi: 10.1080/15569543.2020.1799408.
22. Casanova-Martí, À. et al. (2020). Long Term Exposure to a Grape Seed Proanthocyanidin Extract Enhances L-Cell Differentiation in Intestinal Organoids. *Mol Nutr Food Res.* doi: 10.1002/mnfr.202000303.
23. Fabijańska, M. et al. (2020). Simple Trans-Platinum Complex Bearing 3-Aminoflavone Ligand Could Be a Useful Drug: Structure-Activity Relationship of Platinum Complex in Comparison with Cisplatin. *Int. J. Mol. Sci.* **21**:2116. doi: 10.3390/ijms21062116.
24. Pires, S. et al. (2020). Biological sex influences susceptibility to *Acinetobacter baumannii* pneumonia in mice. *JCI Insight.* pii: 132223. doi: 10.1172/jci.insight.132223.
25. Mavangira, V. et al. (2020). 20-hydroxyeicosatetraenoic acid alters endothelial cell barrier integrity independent of oxidative stress and cell death. *Prostaglandins Other Lipid Mediat.* doi: 10.1016/j.prostaglandins.2020.106425.
26. Pagoni, A. et al. (2019). Novel anti-Alzheimer phenol-lipoyl hybrids: Synthesis, physico-chemical characterization, and biological evaluation. *Eur J Med Chem.* doi: 10.1016/j.ejmech.2019.111880.
27. Vicario, N. et al. (2019). Simultaneous Activation of Mu and Delta Opioid Receptors Reduces Allodynia and Astrocytic Connexin 43 in an Animal Model of Neuropathic Pain. *Mol Neurobiol.* doi: 10.1007/s12035-019-1607-1.
28. Bahmani Jalali, H. et al. (2018). Effective Neural Photostimulation Using Indium-Based Type-II Quantum Dots. *ACS Nano.* **12**(8):8104-8114. doi: 10.1021/acsnano.8b02976.
29. van Krüchten, A. et al. (2018). *Staphylococcus aureus* triggers a shift from influenza virus-induced apoptosis to necrotic cell death. *FASEB J.* **32**(5):2779-2793. doi: 10.1096/fj.201701006R.
30. Geyikoglu, F. et al. (2017). Toxicological evaluation of submerged liquid culture from *Phanerochaete chrysosporium* mycelium on human blood cells: cytotoxicity, genotoxicity and

oxidative damage. *International Journal of Secondary Metabolite*. **4**(2017):319-329. doi: 10.21448/ijsm.373809.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
5628 Copley Drive
San Diego, CA 92111
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2014-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.