



Product Manual

LDH Cytotoxicity WST Assay

Catalog #: ENZ-KIT157

ASSAY FOR THE DETERMINATION OF CYTOTOXICITY BY LACTATE
DEHYDROGENASE ACTIVITY

500 Tests

5 x 500 Tests



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

CONTENTS



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

Introduction	2
Principle of the Assay.....	2
Materials Supplied.....	2
Additional Materials and Equipment Needed.....	3
Precautions	3
Reagents and Sample Preparation	3
Assay Protocol - Homogeneous Assay	3
Optimization of Cell Concentration	3
Cytotoxicity Assay	4
Calculation of Cytotoxicity.....	5
Assay Protocol - Non-Homogeneous Assay	5
Optimization of Cell Concentration	5
Cytotoxicity Assay	6
Cytotoxicity of Mitomycin C Using HeLa Cells.....	7
Contact Information.....	10

INTRODUCTION

LDH Cytotoxicity WST Assay is a kit for the determination of cytotoxicity by measuring 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. LDH Cytotoxicity WST Assay can be used to measure the released LDH according to the following scheme. LDH catalyzes dehydrogenation of lactate to pyruvate thereby 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 the formazan dye thus formed is proportional to that of released LDH into the medium, which is an indication of cytotoxicity.

Since LDH Cytotoxicity WST Assay neither reflects the activity of living cells nor is harmful to cells, cytotoxicity can be measured with the 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.

PRINCIPLE OF THE ASSAY

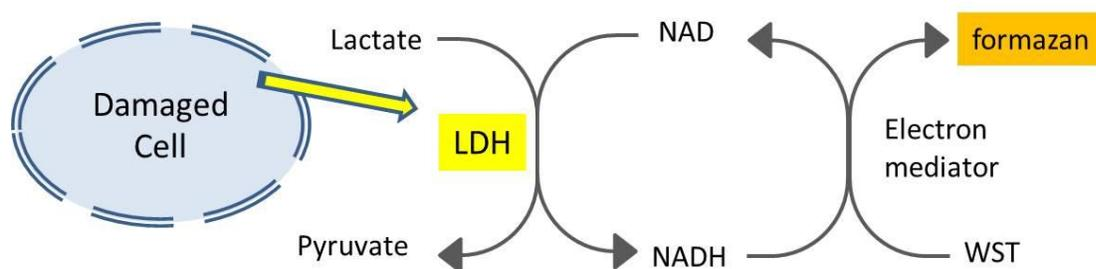


Fig.1 Principle of cytotoxicity measurement

MATERIALS SUPPLIED

	500 tests	5 x 500 tests
Dye Mixture	Bottle x 1	Bottle x 5
Assay Buffer	55 mL x 1	55 mL x 5
Lysis Buffer	5.5 mL x 1	5.5 mL x 5
Stop Solution	27.5 mL x 1	27.5 mL x 5

ADDITIONAL MATERIALS AND EQUIPMENT NEEDED

- CO₂ incubator
- Microplate reader (490 nm filter)
- 96-well tissue culture plate (flat-bottomed). For suspension cells in non-homogeneous assay: Round or V-bottomed plate
- 20 and 100-200 µL multichannel pipettes
- 96-well optically clear plate (flat-bottomed) for non-homogeneous assay

PRECAUTIONS



Handle with care

- This kit contains a glass bottle with an aluminum cap. Use protective gloves and be cautious in handling.
- The amount of LDH is dependent on cell type. We recommend carrying out a preliminary experiment to optimize the cell concentration which is determined by maximum absorbance difference at 490 nm between low and high controls.

REAGENTS AND SAMPLE PREPARATION

1. Add 5 mL of Assay Buffer to the Dye Mixture vial. Close the cap and dissolve the contents completely.
2. Add the whole volume of the mixture prepared in Step 1 to Assay Buffer bottle.
3. Store the Working Solution at 0°C - 5°C and protect it from light. It is stable for 2 months.



Storage temp

ASSAY PROTOCOL – HOMOGENEOUS ASSAY

Optimization of Cell Concentration

1. Collect cells and wash them with the assay medium. Prepare cell suspension to 5.0×10^5 cells/mL in the assay medium.
2. Add 100 µL of the assay medium to each well of a flat-bottom 96-well tissue culture plate.
3. Prepare 2-fold serial dilution of each well in triplicate set of wells for the high-control, low-control and background control (medium only) (Refer to Fig.2 for the plate arrangement).

Serial Dilution Procedure:

- a. Add 100 μL of the cell suspension (5.0×10^5 cells/mL) to the first well [A] and mix by pipetting. This well contains the maximum number of cells (2.5×10^4 cells/well).
- b. Transfer 100 μL from the first well to the next well [B] and mix by pipetting.
- c. Repeat this procedure and discard 100 μL from the last dilution.

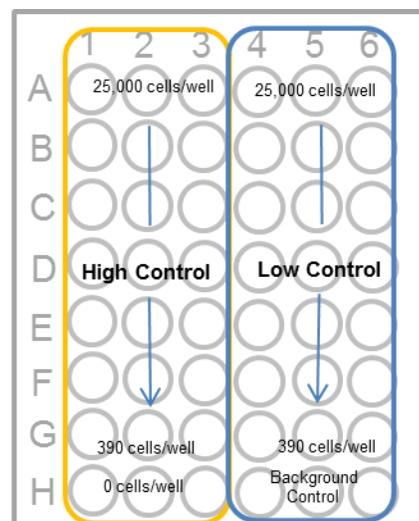


Fig.2 Plate arrangement

4. Incubate the plate at 37°C for an appropriate time in a CO₂ incubator (use the same incubation time in the cytotoxicity assay).
5. Add 10 μL of the Lysis Buffer to each well of the high control.
6. Incubate the plate at 37°C for 30 minutes in a CO₂ incubator.
7. Add 100 μL of the Working Solution to each well. Protect the plate from light and incubate it at room temperature for 30 minutes.
8. Add 50 μL of the Stop Solution to each well.
9. Measure the absorbance at 490 nm by a microplate reader.



Protect from light

Cytotoxicity Assay

1. Add 50 μL of cell suspension to each well of a flat-bottom 96-well tissue culture plate.
For adherent cells: Incubate the plate at 37°C overnight in a CO₂ incubator to allow the cells to adhere tightly. Remove the medium and add 50 μL of fresh assay medium.
2. Add 50 μL of assay medium containing test substance that adjusted to the desired concentration (Refer to Table 1).

Table 1: Amount of each solution (Homogeneous Assay)

	Test Substance	High Control	Low Control	Background Control
Assay medium	-	50 μL	50 μL	100 μL
Cell suspension	50 μL	50 μL	50 μL	-
Test substance in culture medium	50 μL	-	-	-
Lysis Buffer	-	10 μL	-	-

* The difference of total volume of test substance and high control does not affect the result.



Protect from
light

3. Incubate the plate at 37°C for an appropriate time period in a CO₂ incubator.
4. Add 10 µL of the Lysis Buffer to each well of the high control. Incubate the plate at 37°C for 30 minutes in a CO₂ incubator.
5. Add 100 µL of the Working Solution to each well. Protect the plate from light and incubate it at room temperature for 30 minutes.
6. Add 50 µL of the Stop Solution to each well.
7. Measure the absorbance at 490 nm using a microplate reader.

Calculation of Cytotoxicity

Calculate the average absorbance from each triplicate set of wells and subtract the background control value from each absorbance value. Calculate the percent cytotoxicity by the following equation:

$$\text{Cytotoxicity(\%)} = \frac{(A-C)}{(B-C)} \times 100$$

A: Test substance
B: High control
C: Low control

ASSAY PROTOCOL – NON-HOMOGENEOUS ASSAY

Optimization of Cell Concentration

1. Collect cells and wash them with the assay medium. Prepare cell suspension to 5.0×10⁵ cells/mL in the assay medium.
2. Add 100 µL of the assay medium to each well of a 96-well tissue culture plate. Use round or v-bottomed plate for suspension cells, flat-bottomed plate for adherent cells.
3. Prepare 2-fold serial dilution of each well in triplicate set of wells for the high-control, low-control and background control (medium only) (Refer to Fig. 2 for the plate arrangement).

Serial Dilution Procedure:

- a. Add 100 µL of the cell suspension (5.0×10⁵ cells/mL) to the first well [A] and mix by pipetting. This well contains the maximum number of cells (2.5×10⁴ cells/well).
- b. Transfer 100 µL from the first well to the next well [B] and mix by pipetting.

- c. Repeat this procedure and discard 100 μ L from the last dilution.
4. Add 100 μ L of the assay medium to each well.
5. Incubate the plate at 37°C for an appropriate time in a CO₂ incubator (use the same incubation time in the cytotoxicity assay).
6. Add 20 μ L of the Lysis Buffer to each well of the high control.
7. Incubate the plate at 37°C for 30 minutes in a CO₂ incubator.
8. Centrifuge the plate at 250xg for 2 minutes to precipitate the cells (for suspension cells).
9. Transfer 100 μ L of the supernatant from each well to an optically clear 96-well plate.
10. Add 100 μ L of the Working Solution to each well. Protect the plate from light and incubate it at room temperature for 30 minutes.
11. Add 50 μ L of the Stop Solution to each well.
12. Measure the absorbance at 490 nm by a microplate reader.



Protect from light

Cytotoxicity Assay

1. Add 100 μ L of cell suspension to each well of a flat-bottom 96-well tissue culture plate.

For adherent cells: Incubate the plate at 37°C overnight in a CO₂ incubator to allow the cells to adhere tightly. Remove the medium and add 100 μ l of fresh assay medium.
2. Add 100 μ L of the assay medium containing test substance that adjusted to the desired concentration (Refer to Table 2).

Table 2: Amount of each solution (Non-Homogeneous Assay)

	Test Substance	High Control	Low Control	Background Control
Assay medium	20 μ L	100 μ L	120 μ L	220 μ L
Cell suspension	100 μ L	100 μ L	100 μ L	-
Test substance in culture medium	100 μ L	-	-	-
Lysis Buffer	-	20 μ L	-	-

3. Incubate the plate at 37°C for an appropriate time period in a CO₂ incubator.
4. Add 20 μ L of the Lysis Buffer to each well of the high control. Incubate the plate at 37°C for 30 minutes in a CO₂ incubator.

5. Centrifuge the plate at 250xg for 2 minutes to precipitate the cells (for suspension cells).
6. Transfer 100 μ L of the supernatant from each well to each well of a new optically clear 96-well plate.
7. Add 100 μ L of the Working Solution to each well. Protect the plate from light and incubate it at room temperature for 30 minutes.
8. Add 50 μ L of the Stop Solution to each well.
9. Measure the absorbance at 490 nm by a microplate reader.



Protect from light

10.

Cytotoxicity of Mitomycin C Using HeLa Cells

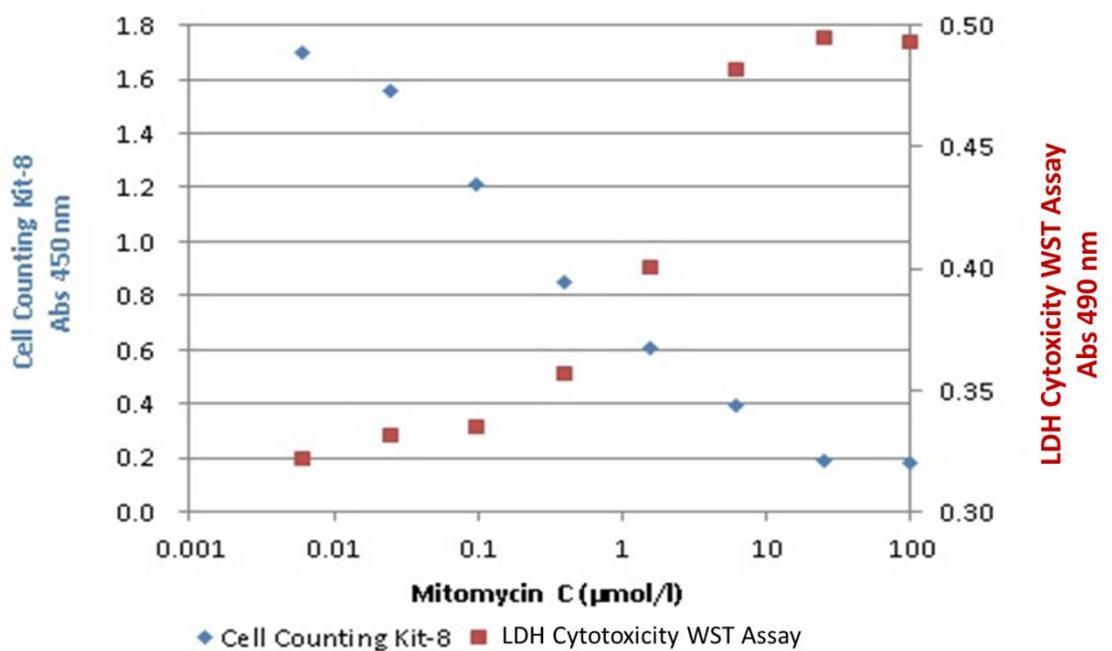


Fig. 3 Cytotoxicity of Mitomycin C Using HeLa cells. Test substance: Mitomycin C, Cell line: HeLa, Culture medium: MEM, 10% FBS, Incubation: 37°C, 5% CO₂, 48 hours



Product Manual

NOTES



Product Manual

NOTES



Product Manual

NOTES



Product Manual

NOTES



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com