

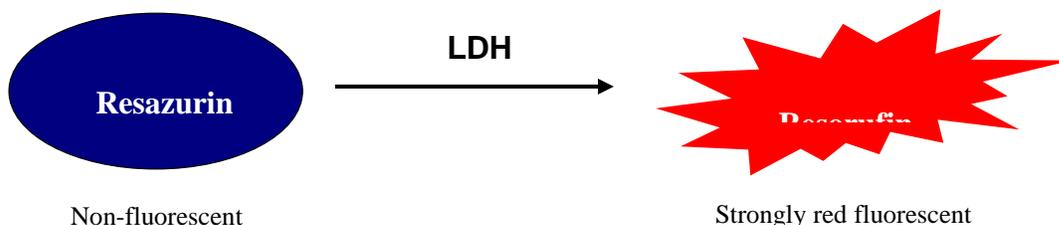
## Cell Meter™ Fluorimetric Cell Cytotoxicity Assay Kit

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 22781 (1,000 assays), 22781-B (5,000 assays)	Keep in freezer and protect from light	Fluorescence microplate readers

### Introduction

The measurement of mitochondrial dehydrogenases (e.g. LDH) activity is a well-accepted assay to quantify cell numbers and monitor cell viability. Our Cell Meter™ Fluorimetric Cell Cytotoxicity Assay Kit provides a fast, simple, accurate and homogeneous assay for the colorimetric or fluorimetric detection of viable cells. This assay is based on the observation that oxidized non-fluorescent blue resazurin is reduced to a red fluorescent dye (resorufin) by accepting an electron from mitochondrial respiratory chain in live cells. The amount of resorufin produced is directly proportional to the number of living cells.

Our Cell Meter™ Fluorimetric Cell Cytotoxicity Assay Kit is more sensitive for cell proliferation and cytotoxicity than other assays such as MTT. The kit components are quite stable with minimal cytotoxicity, thus a longer incubation time (such as 24 to 48 hours) is possible if required. The characteristics of its high sensitivity (<100 CHO cells), non-radioactivity and no-wash method make the kit suitable for high throughput screening of cell proliferation or cytotoxicity against a variety of compounds. The assay can be performed in a convenient 96-well and 384-well microtiter-plate format with a filter set of Ex/Em = ~540/590 nm, cutoff 570 nm.



### Kit Key Features

<i>Non-Radioactive:</i>	No special requirements for waste treatment.
<i>Continuous:</i>	Easily adapted to automation without mixing or separation.
<i>Convenient:</i>	Formulated to have minimal hands-on time.
<i>Wide Applications:</i>	Cell proliferation and cytotoxicity.
<i>Sensitive And Accurate:</i>	As low as 100 cells can be accurately quantified.
<i>Enhanced Value:</i>	Less expensive than the sum of individual components.

### Kit Component

Component	Amount	
	#22781	#22781-B
	1,000 assays (96-well)	5,000 assays (96-well)
	2,000 assays (384-well)	10,000 assays (384-well)
Component A: Assay Solution	20 mL	100 mL

### Materials Required (but not provided)

- 96 or 384-well microplates: Tissue culture microplates with black wall and clear bottom are recommended.
- A fluorescence microplate reader: Capable of detecting emission at 590 nm with excitation at 530-560 nm.

## Assay Protocol for One 96-Well Plate

### Brief Summary

**Prepare cells with test compounds (100 µL/well/96-well plate or 50 µL/well/384-well plate)**  
→ Add 1/5 volume of Assay Solution (Component A) → Incubate at 37 °C  
for 1-24 hours → Monitor fluorescence intensity at Ex/Em = 540/590 nm

#### 1. Prepare cells and test compounds:

- 1.1 Plate 100 to 10,000 cells per well in a tissue culture microplate with black wall and clear bottom. Add test compounds into the cells, and incubate for a desired period of time (such as 24, 48 or 96 hours) in a 37 °C, 5% CO<sub>2</sub> incubator. For blank wells (medium without the cells), add the same amount of compound buffer. The suggested total volume is 100 µL for a 96-well plate, and 50 µL for a 384-well plate.
- 1.2 Set up the following controls at the same time.
  - Positive control contains cells and known proliferation or cytotoxicity inducer.
  - Negative control contains cells but no test compounds.
  - Vehicle control contains cells and the vehicle used to deliver test compounds.
  - Non-cell control contains growth medium without cells.  
*Note: LDH in serum will contribute to background fluorescence.*
  - Test compound control contains the vehicle used to deliver test compounds [Hank's balance salt solution (HBSS) or phosphate-buffered saline (PBS)] and test compound. Some test compounds have strong autofluorescence and may give false positive results.  
*Note: Match the total volume of all the controls to 100 µL for a 96-well plate or 50 µL for a 384-well plate with growth medium.*

#### 2. Assay procedures:

- 2.1 Thaw and warm up the Assay Solution (Component A) to 37 °C, and mix it thoroughly before starting the experiments.
- 2.2 Add 20 µL/well (96-well plate) or 10 µL/well (384-well plate) of Assay Solution (Component A). Mix the reagents by shaking the plate gently for 30 seconds.
- 2.3 Incubate the cells in a 37 °C, 5% CO<sub>2</sub> incubator for 1-24 hours, protected from light.  
*Note 1: The appropriate incubation time depends on the metabolism rate of the individual cell type and cell concentration used. Optimize the incubation time for each experiment.*  
*Note 2: Extremely prolonged incubation time is not recommended since resazurin could be converted to colorless dihydroresorufin.*
- 2.4 Monitor the fluorescence intensity (bottom read) at Ex/Em = 540/590 nm, cutoff 570

#### 3. Perform data analysis:

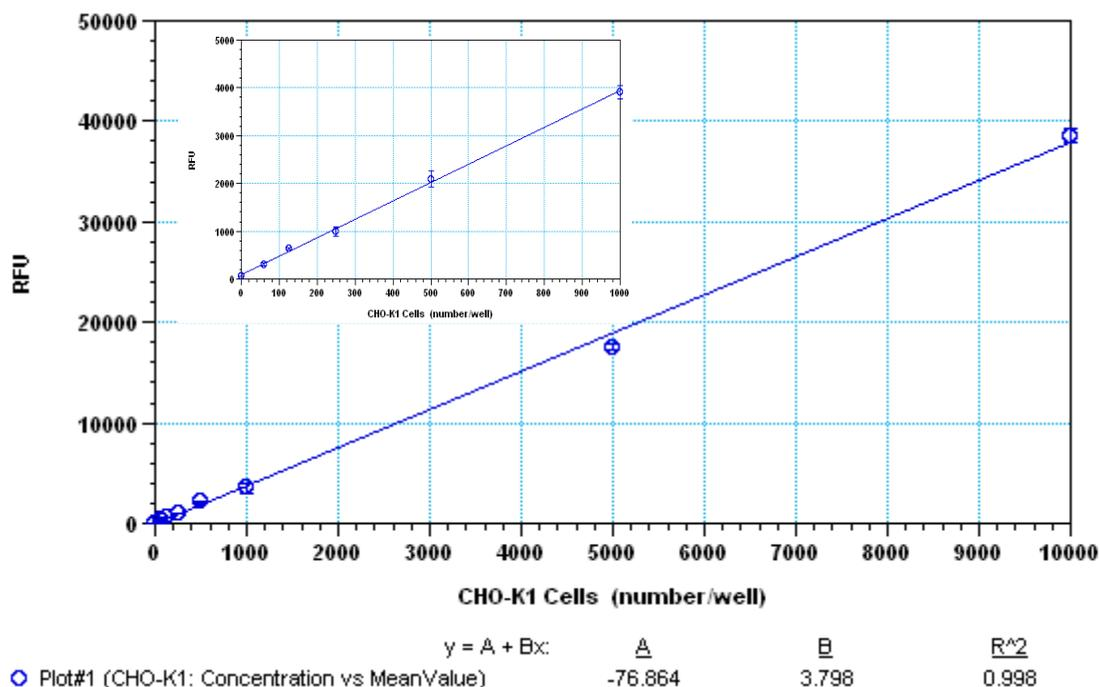
- 3.1 The background fluorescence reading from the non-cell control well is subtracted from the values for those wells containing the cells.  
*Note: The background fluorescence of the blank wells may vary depending on the sources of the growth media or the microtiter plates.*
- 3.2 The fluorescence reading in each well indicates the cell number in that well.
- 3.3 Calculate the percentage of cell viability for samples and controls based on the following formula:

$$\% \text{ Cell viability} = 100 \times (F_{\text{sample}} - F_o) / (F_{\text{ctrl}} - F_o)$$

$F_{\text{sample}}$  is the fluorescence reading in the presence of the test compound.

$F_{\text{ctrl}}$  is the fluorescence reading in the absence of the test compound (vehicle control).

$F_o$  is the averaged background (non-cell control) fluorescence intensity.



**Figure 1.** CHO-K1 cell number response was measured with Cell Meter™ Fluorimetric Cell Cytotoxicity Assay Kit. CHO-K1 cells at 0 to 10,000 cells/well/100  $\mu$ L were seeded overnight in a Costar black wall/clear bottom 96-well plate. The cells were incubated with 20  $\mu$ L/well of Assay Solution (Component A) for 3 hours at 37  $^{\circ}$ C. The fluorescence intensity was measured at Ex/Em = 540/590 nm with NOVOstar instrument (BMG Labtech). The fluorescence intensity was linear ( $R^2 = 0.998$ ) to the cell number as indicated. The detection limit was 60 cells/well ( $n=6$ ). The insert shows the enlargement of the lower end of the cell number response.

## References

1. R. M. DeBAUN and G. de STEVENS, Arch.Biochem. 31, 300-308 (1951).
2. Nachlas, M.M. et al. Anal. Biochem. 1, 317-26 (1960).
3. D. W. De Jong and W. G. Woodlief, Biochim.Biophys.Acta 484, 249-259 (1977).
4. T. Mosmann, J.Immunol.Methods 65, 55-63 (1983).
5. C. Korzeniewski and D. M. Callewaert, J.Immunol.Methods 64, 313-320 (1983).
6. Decker, T. and Lohmann-Matthes, M.L. J. Immunol. Meth. 115, 61-9 (1988).
7. C. Legrand et al., J.Biotechnol. 25, 231-243 (1992).
8. G. Konjevic, V. Jurisic, I. Spuzic, J.Immunol.Methods 200, 199-201 (1997).
9. Singer, C.A. et al. J. Neurosci. 19, 2455-63 (1999).

**Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.**