

CytoX-Red™ Cell Proliferation/ Cytotoxicity Assay Kit

Base Catalog # OP-0004



PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The CytoX-Red™ Cell Proliferation/Cytotoxicity Assay Kit is very suitable for quick measurement of cell viability, proliferation, and cytotoxicity.

110 Bi County Blvd. Ste. 122, Farmingdale, NY 11735

Tel: 1-877-374-4368 ■ Fax: 1-718-484-3956 ■ E-mail: info@epigentek.com ■ Web: www.epigentek.com
© Epigentek Group Inc. All rights reserved. Products are for research use only.

Page 1

Printed 2017-09-12
OP-0004

KIT CONTENTS

Components	3 X 96 assays OP-0004-1	5 X 96 assays OP-0004-2	10 X 96 assays OP-0004-3
Indicator Solution	16 ml	26 ml	5 x 10 ml
Dissolve Solution	30 ml	2 x 25 ml	4 x 25 ml
Reservoirs	2	4	6
User Guide	1	1	1

SHIPPING & STORAGE

Store all components at room temperature for up to 1 year.

GENERAL PRODUCT INFORMATION

Usage Limitations: The CytoX-Red™ Cell Proliferation/Cytotoxicity Assay Kit is for research use only and is not intended for diagnostic or therapeutic application.

Intellectual Property: CytoX-Red™ is a trademark of Epigentek, Inc.

A BRIEF OVERVIEW

Protein synthesis rate is demonstrated to be proportional to cell proliferation and is a very sensitive cytotoxic marker. The CytoX-Red™ Cell Proliferation/Cytotoxicity Assay Kit provides a simple and convenient method for measuring cell viability, proliferation, and cytotoxicity in adherent or monolayer cells through quantifying total protein contents. The Assay using CytoX-Red™ Cell Proliferation/Cytotoxicity Assay Kit is proven to be accurate and yields reproducible results. A complete proliferation/cytotoxicity assay can be finished within 20 minutes. Thus, the CytoX-Red™ provides a considerably faster procedure than any competitor's kits. The kit has the following features:

- *The fastest procedure.* The proprietary composition of the kit allows the fixation step to be eliminated and dye incubation time to be much shorter. Thus, a complete proliferation/cytotoxicity assay with this kit can be finished within 20 minutes.
- *High sensitivity.* The cell number required for signal generation can be as low as 500 cells

PROTOCOL

1. Set up the assayed 96-well plate containing cells in 180 μ l of culture medium.
2. Add 20 μ l of the tested compounds or vehicle control and incubate in 37°C incubator for desired time.
3. Remove the growth medium from the monolayer cells and wash cells with PBS once.

4. Add 50 μ l of indicator solution to each well and incubate for 10-15 min at room temperature.
5. Discard the solution and wash the plate gently with tap water 3-4 times.
6. Add 100 μ l of dissolving solution to dissolve red staining.
7. Measure O.D at 570 nm.
8. Calculation of cell viability or cytotoxicity:

$$\% \text{ viability} = \frac{\text{O.D (treated - blank)}}{\text{O.D (control - blank)}} \times 100\%$$

RELATED PRODUCTS

OP-0005

CytoX-Violet™ Cell Viability/ Cytotoxicity Assay Kit

