

# CytoX-Violet™ Cell Proliferation/ Cytotoxicity Assay Kit

Base Catalog # OP-0005

## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

The CytoX-Violet™ Cell Proliferation/Cytotoxicity Kit is very suitable for the quick measurement of cell viability, proliferation, and cytotoxicity in both adherent and suspension cells.

## KIT CONTENTS

Components	5 x 96 assays OP-0005-1	10 x 96 assays OP-0005-2	20 x 96 assays OP-0005-3
Indicator Solution	11 ml	22 ml	50 ml

## SHIPPING & STORAGE

Upon receipt, store the kit at 4°C away from light. The kit is stable for up to 1 year from the date of shipment, when stored properly.

## GENERAL PRODUCT INFORMATION

**Usage Limitation:** The *CytoX-Violet*<sup>™</sup> Cell Proliferation/Cytotoxicity Kit is for research use only and is not intended for diagnostic or therapeutic application.

**Intellectual Property:** *CytoX-Violet*<sup>™</sup> is a trademark of EpigenTek, Inc.

## A BRIEF OVERVIEW

The *CytoX-Violet*<sup>™</sup> Cell Proliferation/Cytotoxicity Assay Kit utilizes a special dye composition to measure cell viability, proliferation, and cytotoxicity in both adherent and suspension cells by measuring the activity of the cellular dehydrogenases. The fluorometric change of the *Cyto-Violet*<sup>™</sup> indicator solution is directly proportional to cell viability or proliferation. The *CytoX-Violet*<sup>™</sup> has only one step and is proven to be very simple, accurate, and yield reproducible results.

## PROTOCOL

1. Set up assay 96-well plate containing cells in 180 µl of culture medium.
2. Add 20 µl of tested compounds or vehicle control and incubate in 37°C incubator for a desired time.
3. Remove assay plate from incubator and add 20 µl of **Indicator Solution** to each well. Incubate for 3-4 hours in 37°C incubator again.
4. Measure fluorescence at 560/590 nm.
5. Calculate cytotoxicity:

$$\text{Survival \%} = \frac{\text{fluorescent intensity (treated - blank)}}{\text{fluorescent intensity (untreated control - blank)}} \times 100\%$$

## RELATED PRODUCTS

OP-0004      *CytoX-Red*<sup>™</sup> Cell Proliferation/ Cytotoxicity Assay Kit