

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.

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.

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
User Guide	1	1	1

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*[™] Cell Proliferation/Cytotoxicity Kit is for research use only and is not intended for diagnostic or therapeutic application.

Intellectual Property: *CytoX-Violet*[™] is a trademark of Epigentek, Inc.

A BRIEF OVERVIEW

The *CytoX-Violet*[™] 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*[™] indicator solution is directly proportional to cell viability or proliferation. The *CytoX-Violet*[™] 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 μ l of culture medium.
2. Add 20 μ l 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 μ l 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*[™] Cell Proliferation/ Cytotoxicity Assay Kit