Quick PROTOCOL

Non-Radioactive Cell Proliferation Assay

Preparation of Assay Plates

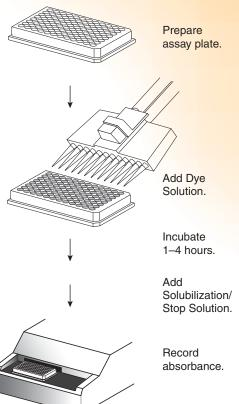
1. Prepare 96-well assay plates containing cells in 100µl culture medium, test compounds and appropriate controls.

Color Development and Recording of Data

- 1. Add 15µl of Dye Solution to each well.
- 2. Incubate the plate at 37° C for 1–4 hours in a humidified CO₂ incubator.
- 3. Add 100µl of Solubilization/Stop Solution to each well. The colored formazan product is stable at 4°C, and absorbance can be recorded in 1 hour or up to several days later.
- 4. Record the absorbance at 570nm using a 96-well plate reader. A reference wavelength between 630–750nm may be used.

Note: To use this system with different volumes, please refer to Section 5 of TB112.

See additional protocol information in Technical Bulletin #TB112, available upon request from Promega or online at **www.promega.com**





Printed in USA. Revised 12/124 Part# 9FB045 © 2000–2012 Promega Corporation. All Rights Reserved.