

# Cell Viability Cytotoxicity (CVC) Reagent

Catalog Number CVC-000X

(For Research Use Only)

# Introduction

Signosis' Cell Viability/Cytotoxicity (CVC) Reagent is a colorimetric WST-8 based assay that provides a simple and accurate method to measure cell proliferation. This is a tetrazolium reduction assays based on the cleavage of the tetrazolium salt WST-8 to water-soluble, orange formazan dye by cellular mitochondrial dehydrogenases. Intensity of dye is a direct proportion of increased activity of the mitochondrial dehydrogenases and therefore number of viable cells which can be quantified by measuring the absorbance at 450 nm.

### Materials provided with the kit

1 ml (CVC-0001) or 5 ml (CVC-0005) of Cell Viability Cytotoxicity (CVC) Reagent.
o Store at 4°C.

### Material required but not provided

- 96 well plate
- Plate reader (450 nm filter)
- CO<sub>2</sub> incubator
- Multi-channel pipettes

# **Assay Procedure**

- 1. The day before performing the assay, trypsinize the cells and seed each well of a 96 well plate with around 1-5 x  $10^4$  cells in 100µl medium.
- 2. Incubate the plate in a humidified incubator at 37°C with 5% CO<sub>2</sub> overnight.
- 3. If not testing substances, skip to step 4. Otherwise, add 10  $\mu$ l of each substance at various concentrations to each well and incubate as appropriate for your substances.
- Carefully add 10 μl of CVC reagent to each well. Note: Do not introduce bubbles to the wells, this will affect the readings.
- 5. Incubate the plate in a humidified incubator at  $37^{\circ}$ C with 5% CO<sub>2</sub> for 1-3 hours.
- 6. Measure the absorbance at 450nm in a microplate reader.