

## Cell Counting Kit 8 (CCK-8)

Catalog number: 22772, 22773  
Unit size: 1000 Tests, 10000 Tests

| Component      | Storage                                    | Amount (Cat No. 22772) | Amount (Cat No. 22773)   |
|----------------|--------------------------------------------|------------------------|--------------------------|
| CCK-8 Solution | Freeze (< -15 °C), Minimize light exposure | 1 bottle (10 mL)       | 2 bottles (50 mL/bottle) |

### OVERVIEW

Cell Counting Kit 8 (CCK-8) is a colorimetric assay used to quantitate cell proliferation, cell viability and cytotoxicity. CCK-8 utilizes the enzymatic activity of live cells to reduce a colorless compound into an orange dye, formazan. The amount of formazan produced, as determined by absorbance ( $\lambda_{460}$  nm), is directly proportional to the number of live cells in the sample. Metabolically active cells are known to produce dehydrogenase enzymes. Under neutral pH conditions and with the aid of an electron acceptor known as 1-methoxy phenazine methosulfate, NADPH-dependent cellular enzymes operate across the cell membrane to mediate the reduction of the colorless, cell-impermeant WST-8 tetrazolium salt. This yields an orange-colored formazan dye which is soluble in cell culture media or aqueous buffers, exhibiting strong absorption at 460 nm. This absorbance is directly proportional to the number of live cells present in the samples, serving as a reliable indicator of cell viability.

### AT A GLANCE

1. Prepare cells in a 96-well plate (100  $\mu$ L/well)
2. Add 10  $\mu$ L of CCK-8 solution to each well
3. Incubate at 37°C for 1 - 4 hours
4. Monitor absorbance at OD = 460 nm

**Important note:** The CCK-8 solution is stable for up to one year if stored at 4°C and protected from light. Store it at <-20°C for longer storage.

### KEY PARAMETERS

#### Absorbance microplate reader

|                   |              |
|-------------------|--------------|
| Absorbance        | 460 nm       |
| Recommended plate | Clear bottom |

### CELL PREPARATION

For guidelines on cell sample preparation, please visit: <https://www.aatbio.com/resources/guides/cell-sample-preparation.html>

### SAMPLE EXPERIMENTAL PROTOCOL

#### Cell Proliferation and Cytotoxicity Assay:

1. Plate 5000 to 10,000 cells/well in a tissue culture microplate with clear bottom.
2. Add test compounds into the cells and incubate for a desired period of time (such as 24, 48, or 96 hours) in a 37°C, 5% CO<sub>2</sub> incubator. For blank wells (medium without the cells), add the same amount of test compounds. The suggested total volume is 100  $\mu$ L for a 96-well plate or 50  $\mu$ L for a 384-well plate.  
**Note:** Each cell line should be evaluated on an individual basis to determine the optimal cell density for proliferation or cytotoxicity induction. For proliferation assays, use fewer cells; for cytotoxicity assays, use more cells to start with.
3. Add 10  $\mu$ L/well (96-well plate) or 5  $\mu$ L/well (384-well plate) of CCK-8 Solution to each well.
4. Incubate the plate at 37°C for 1 - 4 hours, protect from light.  
**Note:** The incubation time could be from 30 minutes to overnight

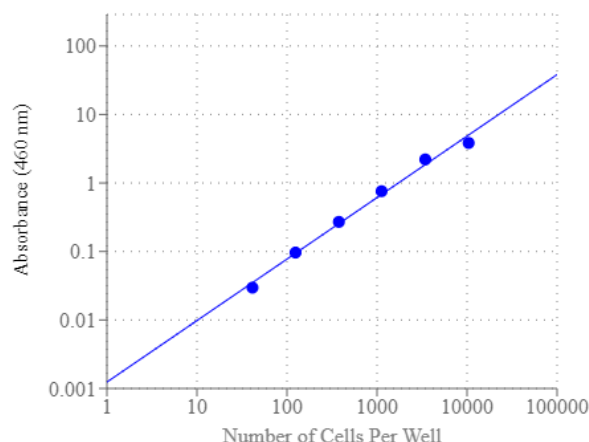
depending on the individual cell type and cell concentration used. Optimize the incubation time for each experiment.

5. Monitor the absorbance increase with an absorbance plate reader at OD = 460 nm.

#### Cell Counting Assay:

1. Prepare cell culture in a tissue culture microplate with a clear bottom. The suggested total volume is 100  $\mu$ L for a 96-well plate or 50  $\mu$ L for a 384-well plate.  
**Note:** We used serially diluted HeLa and Jurkat cell suspension in a clear bottom 96-well plate for the assay.
2. Add 10  $\mu$ L/well (96-well plate) or 5  $\mu$ L/well (384-well plate) of CCK-8 solution to each well.
3. Incubate the plate at 37°C for 1 - 4 hours, protect from light.  
**Note:** The incubation time could be from 30 minutes to overnight depending on the individual cell type and cell concentration used. Optimize the incubation time for each experiment.
4. Monitor the absorbance increase with an absorbance plate reader at OD = 460 nm.

### EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Cell number was determined with Cell Counting Kit 8 (CCK-8). HeLa cells at 40 to 10,000 cells/well/100  $\mu$ L were added in a clear bottom 96-well plate. The absorbance was measured at 460 nm using a SpectraMax reader (Molecular Devices).

### DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.