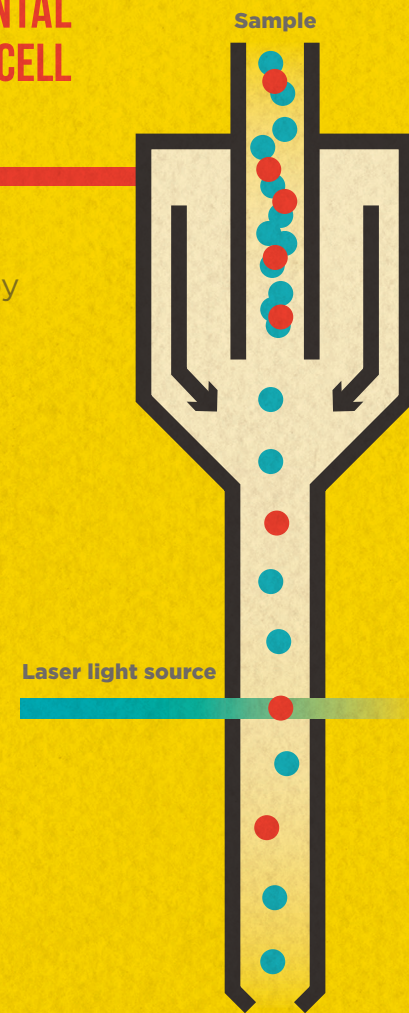
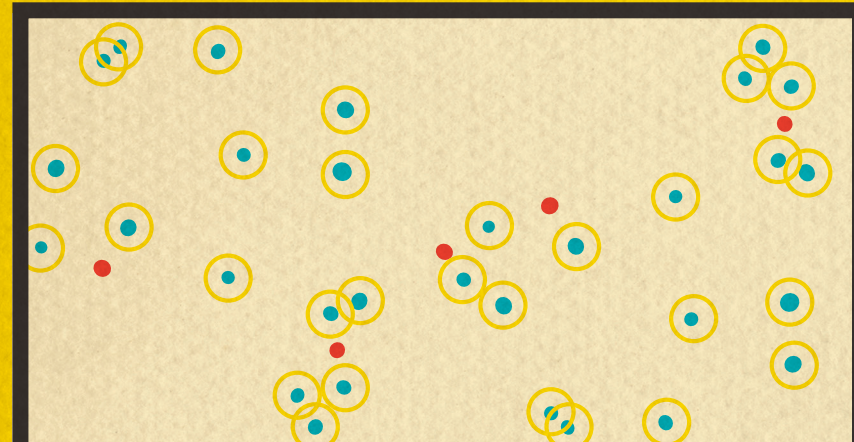


ACCURATE CELL COUNTING IS CRITICAL TO ENSURE EXPERIMENTAL SUCCESS AND REPRODUCIBILITY FOR CELL-BASED ASSAYS AND CELL CULTURE APPLICATIONS.

IMAGE-BASED COUNTERS

Image-based systems use bright field or fluorescent microscopy to capture an image of the cells. Some systems operate on a flow-based imaging methodology where cells are drawn into a capillary and the cells are imaged and counted as they pass a detector. Cell viability can be calculated using dye exclusion methods, such as Trypan Blue. Software analyzes the images based on system specific parameters such as cell diameter, brightness, and circularity to determine the number of cells and cell viability.



# MASTERING CELL COUNTING

COULTER COUNTERS



Coulter counters measure changes in electrical resistance to determine the number, volume and size of the cells in the sample. Some Coulter counters offer the ability to distinguish live cells from dead cells and cell debris.

FLOW CYTOMETERS

Flow cytometers are not dedicated cell counters and report on relative values, such as the percent of cells in a given sample that have specific properties. The volume of sample counted needs to be determined to calculate the absolute cell count. To accomplish this, samples need to be spiked with fluorescent counting beads as a control.

CALCULATING THE NUMBER OF CELLS IS REQUIRED FOR MONITORING CELL PROLIFERATION & VIABILITY, OPTIMIZING CELL CULTURE CONDITIONS, AND PREPARING CELL-BASED ASSAYS.

CELL COUNTING PROCEDURES NEED TO BE:

- Accurate
- Reproducible
- Cell type independent, i.e., take into consideration cell property variables such as cell clumping, cell shape, and cell populations of varying size

HEMOCYTOMETER

DIMENSIONS	AREA	VOLUME AT 0.1 MM DEPTH
1 x 1 mm	1 mm <sup>2</sup>	0.0001 mL
0.25 x 0.25 mm	0.0625 mm <sup>2</sup>	0.00000625 mL
0.20 x 0.20 mm	0.04 mm <sup>2</sup>	0.000004 mL
0.05 x 0.05 mm	0.0025 mm <sup>2</sup>	0.00000025 mL

1. Apply the cell suspension so that the sample fills the chamber.  
*Tip:* Dilute the cell suspension if necessary. Cells should be uniformly distributed without clumping or overlap on the grid.
2. Using a microscope, count the number of cells in 4 squares.  
*Tip:* The lower the concentration of cells, the more squares should be counted to reduce statistical errors.  
*Tip:* Only count cells that are within the square or on the bottom or left lines; do not count cells that touch the top or right lines of the square.
3. Calculate the density of cells in the suspension.

CELL ANALYSIS

**CELL BRIGHTNESS**

Cells & Clutter Counted

Only Cells Counted

**CELL SHARPNESS**

Sharp & Fuzzy Cells Counted

Fuzzy Cells Not Counted

**CELL SPOT AREA**

2% Area - Viable

10% Area - Dead

**DECLUSTER DEGREE**

None

Low

Medium

**MINIMUM CIRCULARITY**

Value of 1 Fewer Cells Counted

Value of .05 More Cells Counted

**SPOT BRIGHTNESS**

Default Value 75 - Classified Viable

Default Value 90 - Classified Dead

## AUTOMATED CELL COUNTING



VS.

## MANUAL CELL COUNTING



AUTOMATED CELL COUNTERS OFFER AN EFFICIENT AND RELIABLE WAY TO QUICKLY COUNT AND ANALYZE A CELL POPULATION.

MANUAL CELL COUNTING OFFERS AN ACCESSIBLE WAY TO DETERMINE THE CONCENTRATION OF CELLS IN A LIQUID SAMPLE, REQUIRING JUST A LIGHT MICROSCOPE AND HEMOCYTOMETER.

$$\text{Cell density (cells/mL)} = \frac{(\text{Average number of cells counted per square}) \times (\text{Dilution factor})}{\text{Volume of square (mL)}}$$

Cell viability can be calculated by treating the cell suspension with cellular dyes, such as Trypan Blue. Live cells do not take up Trypan Blue, dead cells will stain blue.

$$\text{Cell viability (\%)} = \frac{\text{Total viable cells (unstained)}}{\text{Total cells (stained + unstained cells)}} \times 100$$