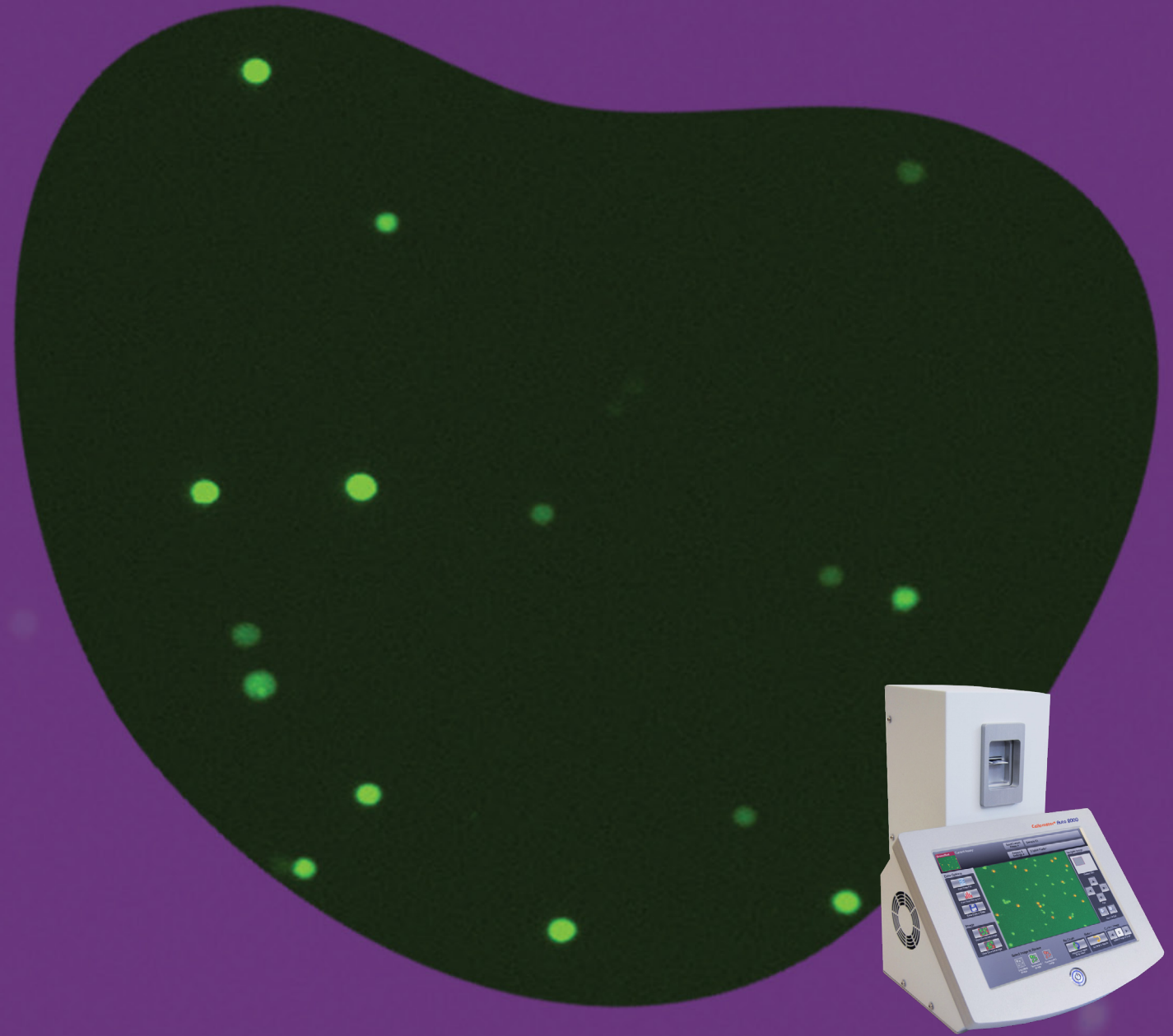


Cell viability  
counter for  
primary cell  
analysis.

revvity



Cellometer® Auto 2000 cell viability counter



# Optimized analysis of primary cells

## Features of the Cellometer Auto 2000

**Dual fluorescence and brightfield imaging:** Staining of both live and dead cells in heterogeneous samples

**Integrated system:** Simple, space-saving design

**User-friendly touchscreen and assay selection:** Enhanced inter-operator reproducibility, minimal training, auto-save option

**Fast results:** Obtain cell images, counts, size measurements, and viability calculations in 30 seconds

**Small sample size:** Only 20  $\mu$ l of sample

**Broad dynamic range:** Measurable concentration range of  $1 \times 10^5$  to  $1 \times 10^7$  cells/mL using Revvity's patent-pending de-clustering function

**Many compatible dyes:** Trypan blue, AO, PI, EB, 7AAD, AO/PI, AO/EB, Calcein AM, CFDA, Calcein AM/PI, CFDA/PI



| Cellometer Auto 2000 cell viability counter





# Imagry for greater results

## Advantages of Cellometer Auto 2000 cell viability counter

### Cell imaging

- Verify cell morphology and counted live/dead cells
- Export cell images for presentations and publications

### Pattern recognition software

- Accurately count cells in clumps
- Count irregular-shaped cells
- Eliminate debris from cell counts
- Differentiate cells based on size

### Automated data management

- Pre-set assays and automated reports
- Archive sample images and auto-save results

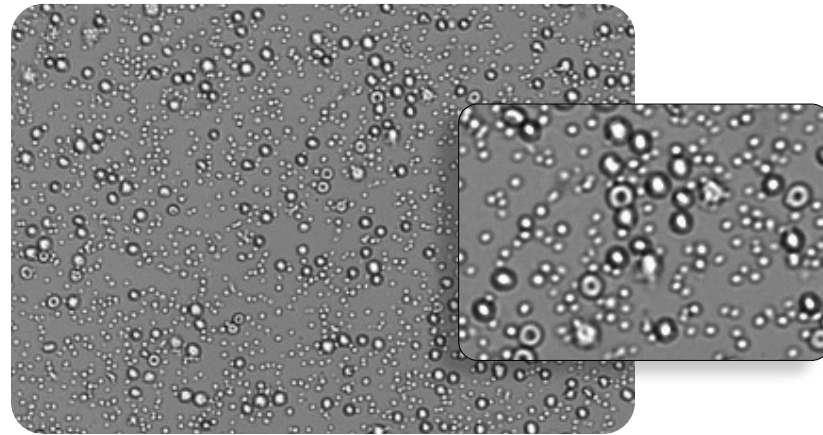
### Maintenance-free system

- Disposable counting chambers – no wash steps
- No required instrument maintenance

“I like the Cellometer Auto 2000 because it eliminates manual counting and our counts are consistent between users. We count cells from primary samples and I like that the RBCs are not counted when we use the AO/PI stain”

*- Moffitt Cancer Center*

# Primary cell analysis



## **PBMC analysis in the presence of red blood cells**

Measure PBMCs from whole blood without lysing. Obtain baseline PBMC concentration and viability prior to biomarker studies.

## **Nucleated cell concentration and viability**

Evaluate cord blood and bone marrow samples

## **GFP transfection efficiency and viability**

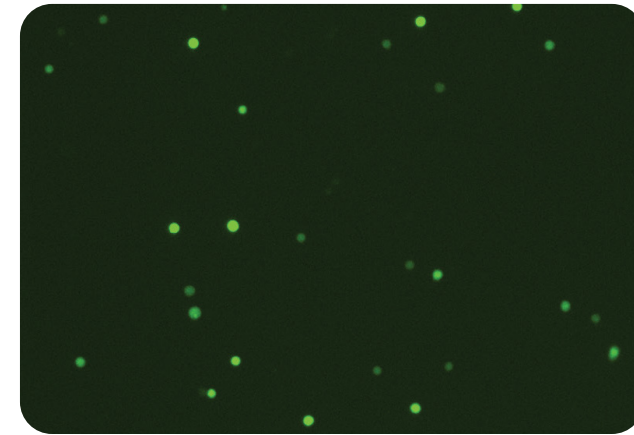
Quickly and easily monitor DNA, RNA, and siRNA transfection

## **Analysis of clumpy and irregular-shaped cells**

Revvity's exclusive pattern-recognition software enables accurate analysis of >98% of mammalian cell types

## **Cell line analysis**

Automatically capture fluorescent cell images, concentration, Trypan blue or PI viability, and mean diameter in 30 seconds!



## **Primary cell analysis**

Accurate concentration and % viability for primary cells (PBMCs, stem cells, splenocytes, neural cells, and more)

## **Analysis of cells from heterogeneous samples**

- Whole Blood
- Peripheral Blood
- Cord Blood
- Bone Marrow



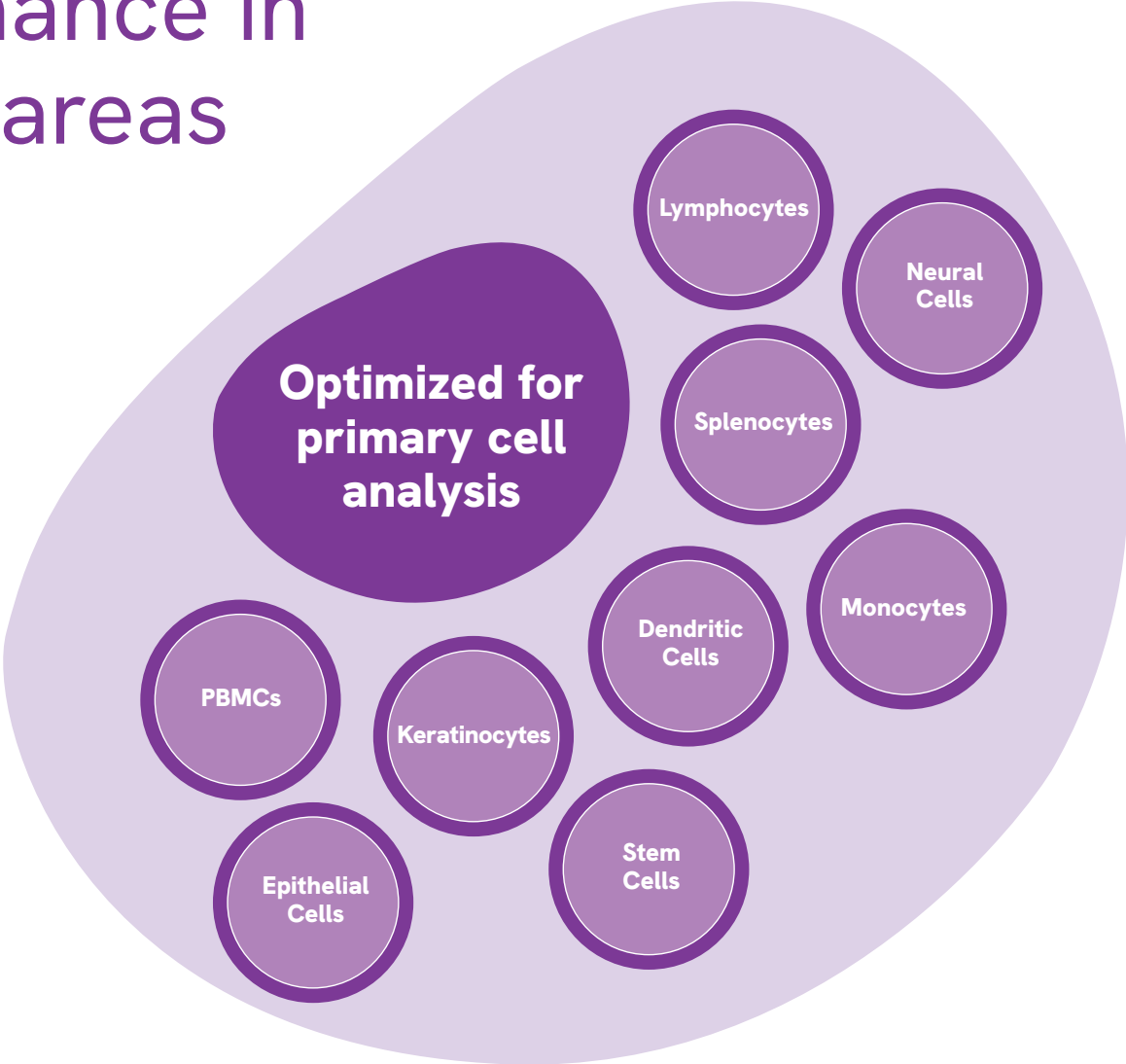


# Proven performance in many research areas

- **Clinical Immunology:** PBMCs
- **Regenerative Medicine:** stem cells
- **Transplantation:** nucleated cells
- **Vaccine Development:** splenocytes
- **Oncology:** cell lines
- **Basic Research:** primary cells / cell lines

“My colleague and I purchased a Cellometer Auto 2000 cell counter and we are using it now. It has facilitated our work greatly. We routinely process PBMCs from both fresh whole blood and from frozen stock. The Cellometer has made it much easier to get cell numbers and viability percentages for use in downstream applications such as IVS and Elispot.

- Human Longevity, Inc



# Dual fluorescence

Primary cell viability in heterogeneous samples live/dead cell concentration using AO/PI

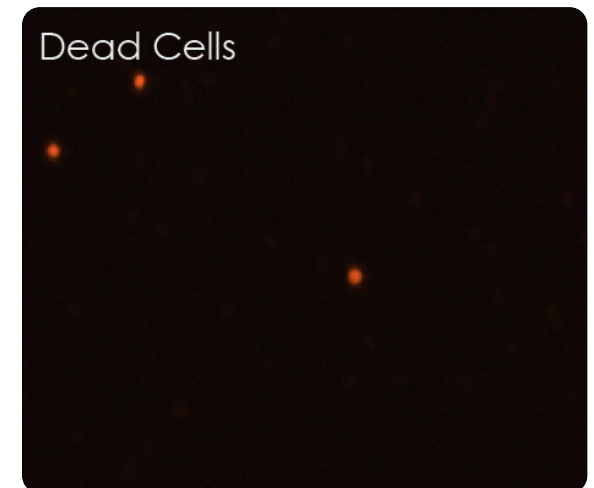
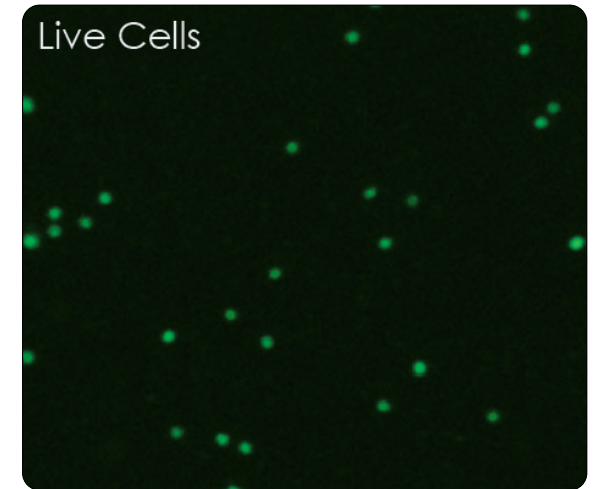
## Why isn't Trypan blue recommended for viability analysis of primary cells?

Trypan blue dye enters and stains all cells with a compromised membrane, including both nucleated and non-nucleated cells, such as red blood cells. For the most accurate calculation of nucleated cell viability, fluorescent nuclear staining dyes are required.

**Dual-fluorescence viability**, using acridine orange (AO) and propidium iodide (PI), is the recommended method for accurate viability analysis of primary cells, such as PBMCs, splenocytes, and stem cells, in samples containing debris and unwanted non-nucleated cell types including red blood cells.

Acridine orange (AO) and propidium iodide (PI) are nuclear staining (nucleic acid binding) dyes. AO is permeable to both live and dead cells and stains all nucleated cells to generate green fluorescence. PI enters dead cells with compromised membranes and stains all dead nucleated cells to generate red fluorescence.

Because mature mammalian red blood cells do not contain nuclei, only live and dead mononuclear cells produce a fluorescent signal. There is no need to lyse red blood cells, saving time and eliminating an extra sample preparation step.





# Results that count

The results indicate the accuracy of the Cellometer Auto 2000 instrument in assessing the viability of Jurkat cells using PI for cell viability. Four measurements were performed for each sample. The viability average was calculated and plotted. The results show the reliability and accuracy of the Cellometer Auto 2000 in measuring cell concentration and viability of mammalian cells.

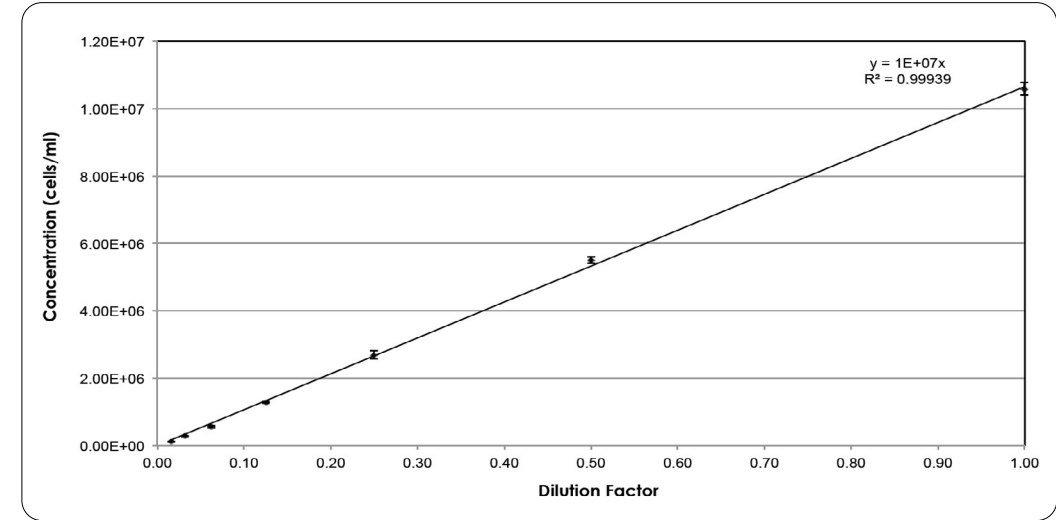


Figure 1: Table of results for cell concentration. Data shown depicts the dynamic range for cell concentration measurements on Cellometer Auto 2000.

The concentration can be measured from  $1 \times 10^5$  -  $1 \times 10^7$  cells / mL without further dilution. The %CV at each concentration was below 10%. This data set was taken on a concentration series of primary mouse splenocytes.

Table 1: Results for cell viability using PI only.

Sample	N Value	Average live cell concentration	% Viability	CV of concentration	CV of viability
A	4	4.20E+06	91.1	10%	2%
B	4	1.06E+06	22.7	7%	1%
C	4	3.27E+06	57.5	7%	7%



# Features

**Analysis of multiple species of stem cells**

**Easily edit and import assays**

**User-friendly touch screen**

**Total nucleated cell count & viability**

**One-Step cell concentration and viability**

**Images for data verification**

**Cell size histograms**

**GFP transfection efficiency and % viability**

**Primary splenocyte concentration & viability**

**Accurate PBMC counts in the presence of red blood cells**

**Cell Diameter Histogram (Percentage of Total Count)**

Size Microns (um)	Percent (%)
7.5	1
8.0	2
8.5	3
9.0	4
9.5	5
10.0	6
10.5	8
11.0	10
11.5	12
12.0	15
12.5	18
13.0	20
13.5	22
14.0	23
14.5	24
15.0	22
15.5	18
16.0	15
16.5	12
17.0	10
17.5	8
18.0	6
18.5	5
19.0	4
19.5	3
20.0	2
20.5	1
21.0	1
21.5	1
22.0	1
22.5	1
23.0	1
23.5	1
24.0	1
24.5	1

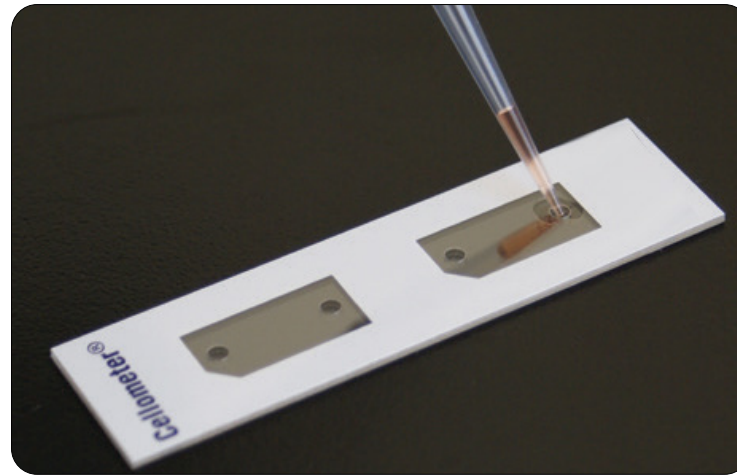
For research use only. Not for use in diagnostic procedures.





# How it works

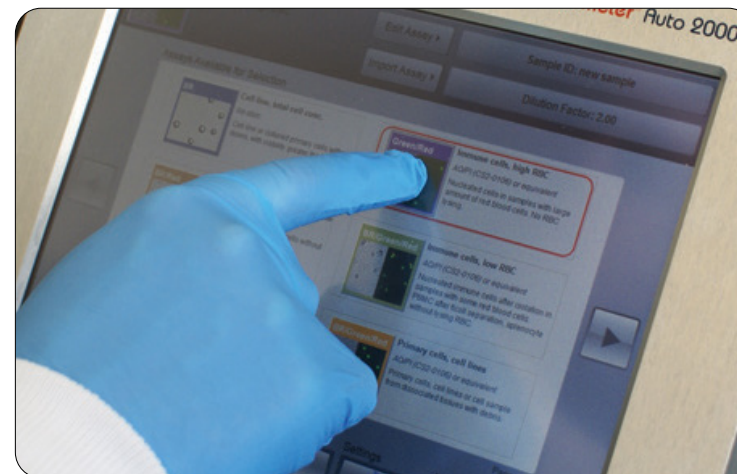
**Step 1: Pipette 20µl**



**Step 2: Insert counting chamber**



**Step 3: Select assay and click count**

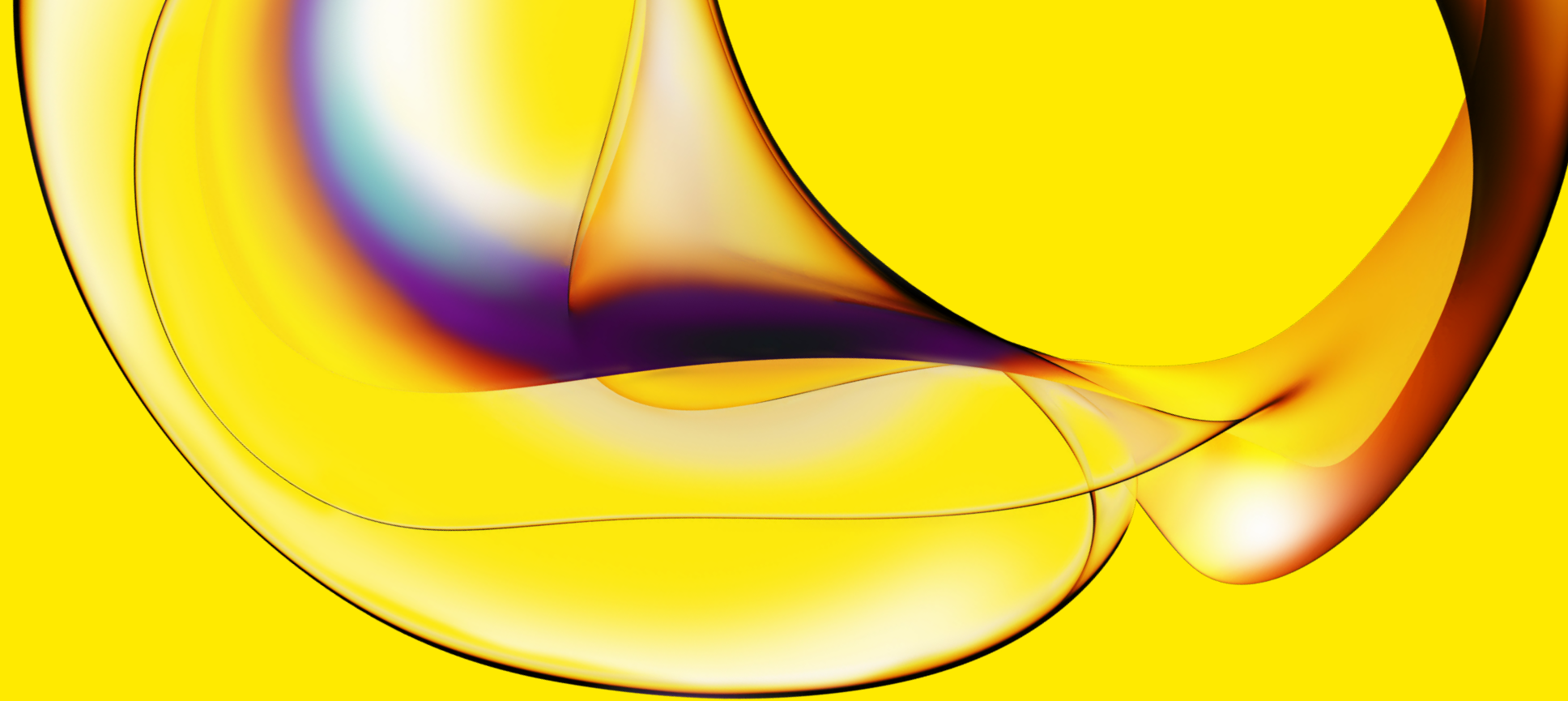


**Step 4: Get results**

**Assay: Immune cells, high RBC**

Sample ID: Blood\_AOPI\_4-2  
Dilution Factor: 2.00

Count	Concentration
=====	=====
Total: 340 cells	1.18x10 <sup>6</sup> cells/mL
Live: 324 cells	1.12x10 <sup>6</sup> cells/mL
Dead: 16 cells	5.53x10 <sup>4</sup> cells/mL



[www.revivity.com](http://www.revivity.com)

revivity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revivity.com](http://www.revivity.com)

For a complete listing of our global offices, visit [www.revivity.com](http://www.revivity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.

1167057