

TC20 Automated Cell Counter

BIO-RAD

With its innovative auto-focus technology and sophisticated cell counting algorithm, the TC20 Automated Cell Counter eliminates subjectivity while delivering reliable results in 30 seconds. Whether you work with cell lines or primary cells, the TC20 Counter gives you accurate, highly reproducible cell counts. Depending on the complexity of your samples, you can let the cell counting algorithm do all the work or select a population of interest by adjusting cell size gates.

Avoid the tedium of manually counting cells and move on to other important tasks. It all adds up to a better workflow for your workday.

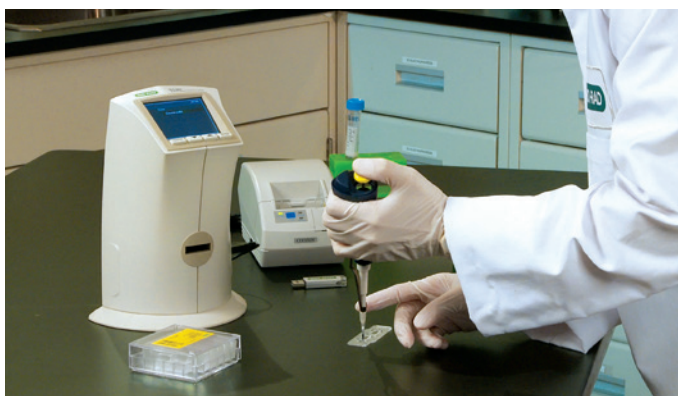
TC20 Automated Cell Counter





Key features and benefits include:

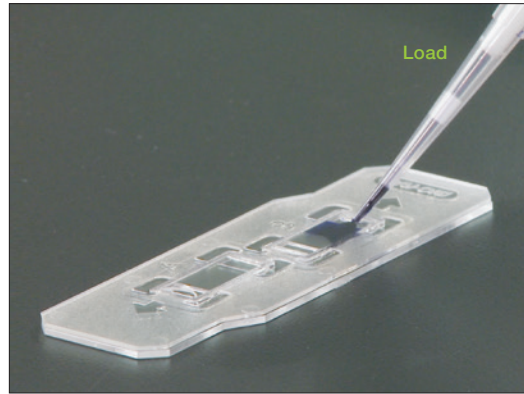
- **Compatible with a broad range of cell sizes and types** — counts cell lines, primary cells, and stem cells
- **Innovative auto-focus technology** — removes the variation associated with manual focusing and leads to precise cell counts in 30 seconds
- **Cell size gates** — user selects a population of interest in complex samples, such as primary cells, or lets the cell counting algorithm do all the work
- **Cell viability** — analyzes cells accurately using multifocal plane analysis
- **Easy archiving and analysis** — store up to 100 counts in onboard memory to view any time, or export counts to your PC for analysis in the optional TC20 Data Analyzer Software



The TC20 Cell Counter is an extremely compact, stand-alone instrument that does not require a computer to operate. With its small footprint, the TC20 Cell Counter conveniently fits any laboratory setting with limited benchspace. Its fast setup and intuitive operation let you quickly and easily start counting cells.

Reliably Accurate

Accelerate your research with automated cell counting. The TC20 Cell Counter performs automated counting of cells in one simple step, immediately initiating a count when you insert the slide. Its built-in auto-focus and sophisticated image analysis algorithm produce accurate, consistent cell counts in 30 seconds.



Load the sample onto a slide.



Insert the slide into the TC20 Cell Counter.

Accurate and Reproducible Results

The TC20 Automated Cell Counter uses microscopy with auto-focus that analyzes multiple focal planes to identify the best plane. Without requiring any user input, the sophisticated cell counting algorithm uses the image acquired from the best focal plane to identify cells and exclude debris, thereby calculating the total cell count. The auto-focus leads to highly reproducible cell counts with reduced user-to-user variability compared to a hemocytometer and cell counters with manual focus.

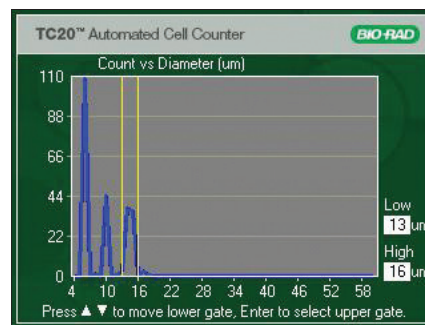
In less than 30 seconds, the TC20 Cell Counter provides accuracy comparable to results obtained with a hemocytometer. It can count cells with a 6–50 μm cell diameter and within a broad concentration range of 5×10^4 – 1×10^7 cells/ml, which eliminates the need to dilute cells, thus reducing the errors associated with sample dilutions prior to counting. In addition, the counting algorithm counts individual cells within clusters of up to five cells, providing accurate counts without the need to extensively declump cells prior to loading.

The TC20 Counter uses a patented* counting slide designed to ensure an even distribution of cells throughout the counting chamber, regardless of the user's pipetting style, leading to accurate and consistent cell counts.

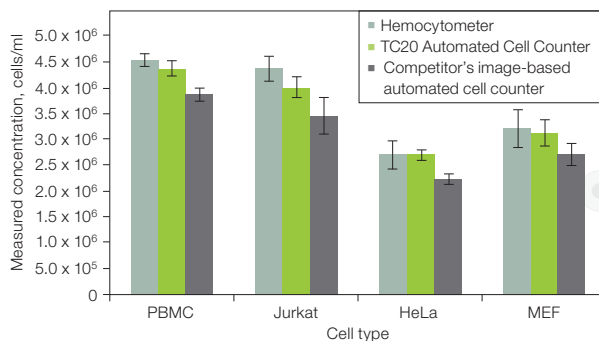
The accurate, consistent cell counts from the TC20 Counter lead to more reproducible results with downstream processes and analysis, saving money and time by enabling successful experiments the first time.

Cell Size Gating

For complex samples composed of multiple cell populations, such as primary cells, users can adjust the cell size gates to define the population of interest that will be counted. When counting multiple sample replicates, the TC20 Cell Counter can save the positions of gates and apply them to subsequent counts.

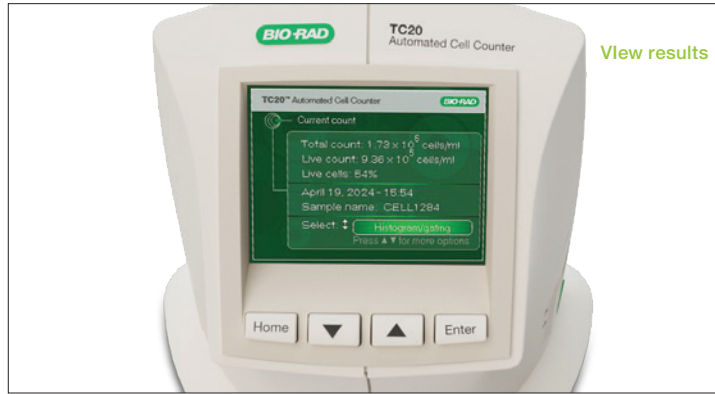
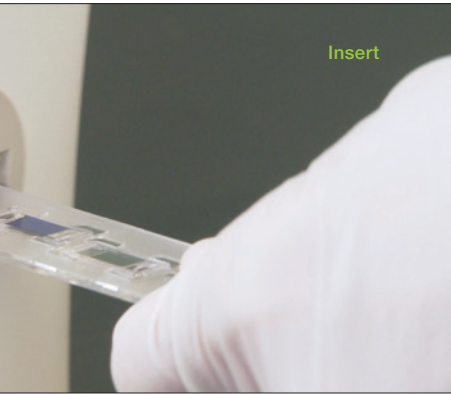


When user-defined gates are enabled you can select a population of interest by adjusting the position of the cell size gates.



The TC20 Cell Counter demonstrates accurate cell counts across a range of cell sizes. Small (PBMC, Jurkat), medium (HeLa), and large (MEF) cells were counted with a hemocytometer, a TC20 Automated Cell Counter, and a competitor's image-based automated cell counter. The TC20 Counter and hemocytometer cell counts showed no statistically significant differences. Precision is indicated by the standard deviations; error bars represent average standard deviations. Cell counts on the TC20 Counter were performed on one instrument with four sample replicates.

* U.S. patent 8,852,524 B2.

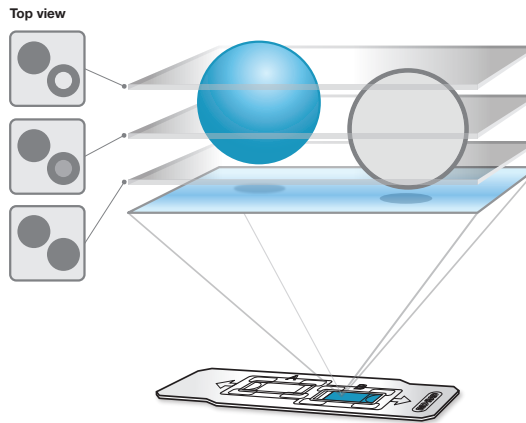


TC20 Automated Cell Counter; counting automatically begins.

Obtain a total cell count (without trypan blue) or total and live cell counts (with trypan blue) in 30 seconds.

Cell Viability

The TC20 Cell Counter can count samples with or without trypan blue. It auto-detects the presence of trypan blue in the sample to assess cell viability via trypan blue exclusion. Along with the total cell count, the TC20 Counter assesses cell viability and provides a live cell count and percentage of live cells. The conventional method of analyzing viability using a single focal plane can lead to inaccurate conclusions because light scattering and the alignment of cells at different heights in a counting chamber — live cells may appear to be dead and vice versa. To determine if cells are viable, the TC20 Counter analyzes each cell using images acquired from multiple focal planes during the focusing step.

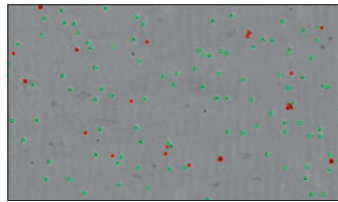


Effect of multifocal plane analysis on live/dead count accuracy for HeLa cells.

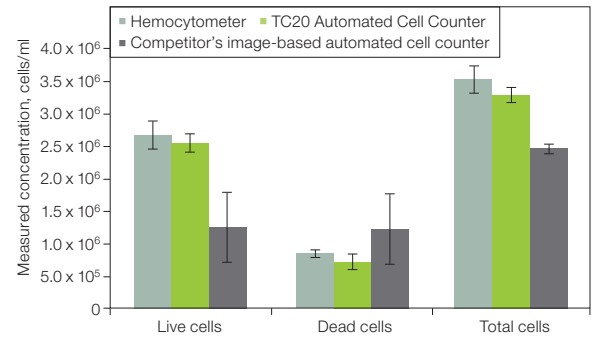
Viability Assessment	Live Cells, %	Dead Cells, %
Single best focal plane*	35	65
Multiple focal planes	65	35

* On single focal plane, 30% of live cells were misidentified as dead cells.

Multifocal plane analysis.



Assessment of cell viability of Pan T cells via trypan blue exclusion. Annotated image is from a cell count exported from the TC20 Counter and viewed as a JPEG image on a computer. Green circles, live cells; red circles, dead cells.



The TC20 Cell Counter demonstrates accurate counts of viable cells. Pan T cells mixed with trypan blue (1:1) were counted with a hemocytometer, a TC20 Automated Cell Counter, and a competitor's image-based automated cell counter. The TC20 Counter and hemocytometer cell counts showed no statistically significant differences. Precision is indicated by the standard deviations; error bars represent average standard deviations. Cell counts on the TC20 Counter were performed on four different instruments with five sample replicates.



Typical Experiment Workflow

Bio-Rad offers powerful building blocks for your cellular research, providing the flexibility and reliability you need to accelerate discovery. The TC20 Automated Cell Counter takes the guesswork out of cell counting, giving you confidence in your results. Use it in combination with other analysis tools from Bio-Rad to streamline your experiments and get more reproducible results.



Grow cells



Count cells



Perform siRNA

Analysis Options

After viewing cell count results, an image of the counted cells can be viewed on the TC20 Counter.

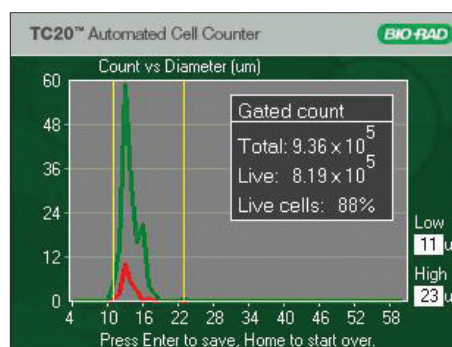
A JPEG file of the image containing a metatag of the count data is available for immediate export to a USB flash drive or, to save every image, automatic data export can be enabled. Use TC20 Data Analyzer Software to review single images, perform multi-file comparative analyses, and generate reports. Go to bio-rad.com/TC20DataAnalyzer to download the software.

Results from 100 previous counts are stored in the TC20 Cell Counter. Previous count results can be exported via the USB port and opened in a Microsoft Excel spreadsheet.

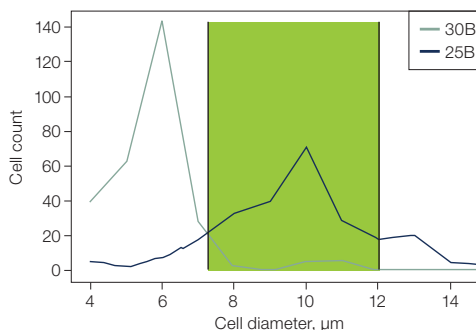
You can assign a name to each sample. To save time when working with multiple sample replicates, enable the automatic name serialization option to populate numerically serialized names.

The optional thermal label printer can be connected to the TC20 Counter to print results onto labels that can be placed into a laboratory notebook for traceability of the count results.

Users can access the onboard dilution calculator to determine volume adjustments required to achieve the cell concentration needed for the next experiment.



Histogram of live and dead cell counts (sample with trypan blue) in the Current count screen.



Single image file analysis using TC20 Data Analyzer Software. Distribution of a total cell population is shown.





A transfection



Count cells

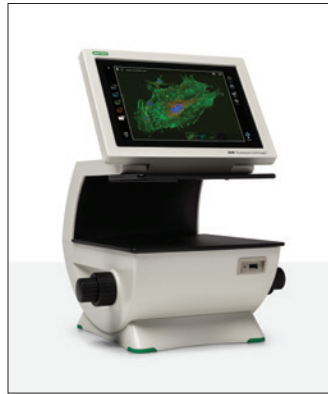


Image cells



Use qPCR to quantitate effect of silencing

Counting Slides

The TC20 Automated Cell Counter uses a patented* counting slide designed to ensure even distribution of cells throughout the counting chamber, regardless of the user's pipetting style, leading to accurate and consistent cell counts.

Disposable counting slides eliminate setup, cleaning, and maintenance steps, and also minimize exposure to biohazardous samples. The dual-chamber slide can provide counts for two separate samples or dilutions. Each chamber requires only 10 μ l, saving precious cells.



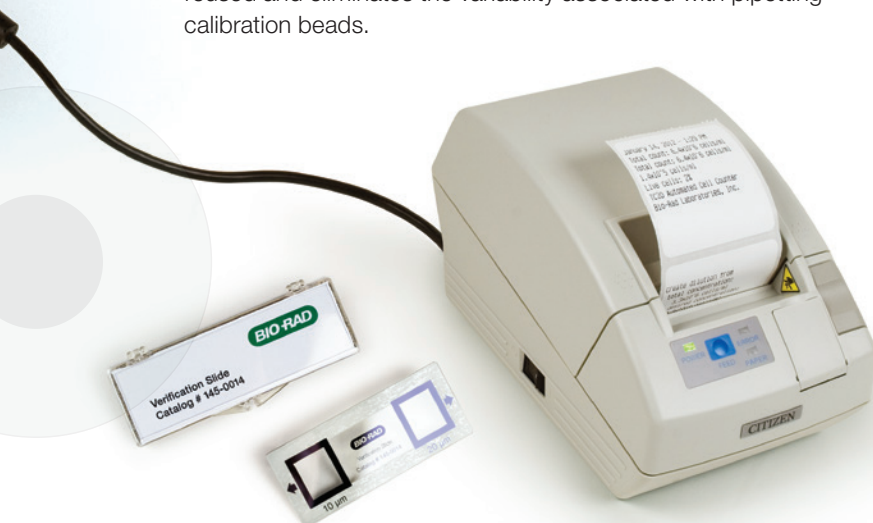
Verification

The ready-to-use verification slide qualifies the functionality of the TC20 Cell Counter. It also serves as a positive control for cell counting experiments. The verification slide can be reused and eliminates the variability associated with pipetting calibration beads.

Counting Bio-Plex™ Magnetic or Polystyrene Beads

The Bio-Plex Suspension Array System is a powerful tool for measuring analyte concentration. While a variety of validated assays is currently available, many Bio-Plex users are developing custom assays for which it is essential to accurately and consistently determine bead concentration to ensure that uniform quantities of antibody are used in the bead-antibody conjugation step. The TC20 Automated Cell Counter is an excellent tool for bead counting and provides speed, accuracy, and ease of use.

* U.S. patent 8,852,524 B2.



Cell Lines Counted on the TC20 Automated Cell Counter*,**

Cell Name	Morphology	Organism	Source Organ/Disease	Growth Property
CCD-1137Sk	Fibroblast	Human	Skin (foreskin)	Adherent
CHO	Epithelial	Chinese hamster	Ovary	Adherent
COS-7	Fibroblast	African green monkey	Kidney	Adherent
HeLa	Epithelial	Human	Cervix carcinoma	Adherent
Jurkat	Lymphoblast	Human	Acute T-cell leukemia	Suspension
K562	Lymphoblast	Human	Bone marrow, chronic myelogenous leukemia	Suspension
MCF-7	Epithelial	Human	Mammary gland, adenocarcinoma	Adherent
MDCK	Epithelial	Dog	Kidney	Adherent
MEF	Fibroblast	Mouse	Embryo	Adherent
mESC	Stem cell	Mouse	Embryo, primary extraction	Adherent
NIH 3T3	Fibroblast	Mouse	Embryo	Adherent
Pan T	Lymphocyte precursor	Human	Blood	Suspension
PBMC	Mixture of different blood cells	Human	Blood, primary extraction	Suspension
PP034	Lymphoid	Human	B cell	Suspension

* The TC20 Counter can count cells with a round shape after trypsinization. ** Go to bio-rad.com/TC20 for a current list of cell lines.

Other Samples Counted on the TC20 Automated Cell Counter*

Sample Name	Organism
<i>Amoeba</i> species	Unicellular protozoa
<i>Chlamydomonas</i>	Unicellular flagellates
<i>Entamoeba histolytica</i>	Unicellular protozoan
<i>Myxomycetes</i>	Slime mold spores
<i>Saccharomyces cerevisiae</i>	Unicellular fungus

* Total cell count only.

Specifications

Parameter	Specification
Counting time	30 sec
Cell concentration range	5 x 10 ⁴ –1 x 10 ⁷ cells/ml
Cell diameter range	6–50 µm
Sample volume	10 µl
Data storage	100 counts
Data export	Via USB drive
Dimensions (W x D x H)	19 x 15 x 25.4 cm (7.5 x 6 x 10 in.)
Weight	2.2 kg (4.8 lb) without the external power supply
Use	For Research Use Only. Not for use in diagnostic procedures.

Ordering Information

Catalog # Description

TC20 Automated Cell Counter

1450102	TC20 Automated Cell Counter , 120–240 V, includes instrument, power supply, USB flash drive, USB cable, 30 dual-chamber counting slides (60 counts), 1.5 ml trypan blue
1450103	TC20 Automated Cell Counter with Thermal Label Printer , 120–240 V, includes instrument, power supply, USB flash drive, USB cable, thermal label printer, 1 roll of 185 labels, 30 dual-chamber counting slides (60 counts), 1.5 ml trypan blue

Kits and Reagents*

1450003	Cell Counting Kit , includes 30 dual-chamber counting slides (60 counts), 1.5 ml trypan blue
1450014	System Test Kit , includes verification slide, instructions
1450021	Trypan Blue , 5 x 1.5 ml 0.4% trypan blue (750 counts), sterile filtered
1450022	Trypan Blue , 10 x 1.5 ml 0.4% trypan blue (1,500 counts), sterile filtered

Accessories*

1450005	Thermal Label Printer , 120–240 V, includes thermal label printer, USB cable, 1 roll of 500 labels
1450007	Thermal Printer Labels , 1 roll of 500 labels, for thermal label printer
1450011	Cell Counting Slides , 30 dual-chamber counting slides (60 counts)
1450015	Cell Counting Slides , 150 dual-chamber counting slides (300 counts)
1450016	Cell Counting Slides , 300 dual-chamber counting slides (600 counts)
1450017	Cell Counting Slides , 600 dual-chamber counting slides (1,200 counts)
1450018	Cell Counting Slides , 900 dual-chamber counting slides (1,800 counts)
1450019	Cell Counting Slides , 1,200 dual-chamber counting slides (2,400 counts)
1450020	Cell Counting Slides , 2,400 dual-chamber counting slides (4,800 counts)

* TC20 kits, reagents, and accessories are compatible with the TC10 Automated Cell Counter.

Go to bio-rad.com/TC20DataAnalyzer to download the software.

Visit bio-rad.com/TC20more for more information.

BIO-RAD and BIO-PLEX are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

For Research Use Only. Not for use in diagnostic procedures.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

