



## Insect Cell Analyses Using the Beckman Coulter® Vi-CELL™

### VI-CELL XR™

Part of the Cell Lab family



#### Insect Cell Analyses

Insect cells are used to produce therapeutic proteins. The specific gene coding the protein of interest is introduced into the cell by a DNA virus of the family Baculoviridae. Protein production via this mechanism, virus infection of the insect cell, is referred to as the Baculovirus Vector System. The most common insect cell used is the SF9 cell or *Spodoptera frugiperda* species.

#### Equipment Used

Beckman Coulter  
Vi-CELL XR  
Validated Vi-CELL XR reagent pack

#### Instrument Settings

When using the Vi-CELL most cells can be analyzed using the default instrument settings. In select cases, certain cell types may require modification of the default analyses parameters. The Vi-CELL software allows for easy user selection of the optimum parameters thus ensuring accurate results. The Vi-CELL settings for SF9 insect cells are the following:

<b>Min. Size (<math>\mu\text{m}</math>) = 9</b>	<b>Cell Brightness = 85</b>
<b>Max. Size (<math>\mu\text{m}</math>) = 50</b>	<b>Cell Sharpness = 100</b>
<b>Number of Images = 50</b>	<b>Viable Cell Spot Brightness = 75</b>
<b>Aspirate Cycle = 1</b>	<b>Min. Circularity = 0</b>
<b>Trypan Blue Mixing Cycle = 2</b>	<b>Decluster Degree = Low</b>

#### Results

The results reported by the Vi-CELL are shown in the Results Section of Figure 1. Percent viability and concentration were 95.1% and  $2.01 \times 10^6/\text{ml}$ . The Vi-CELL also reports viable cell concentration, mean cell diameter, and cell circularity. Graphs may be expanded via a mouse click. Real time cell images may be archived for future re-analyses.

#### Conclusion

The Beckman Coulter Vi-CELL automates the standard manual method, providing accurate results and removing the subjective nature inherent in the manual method.

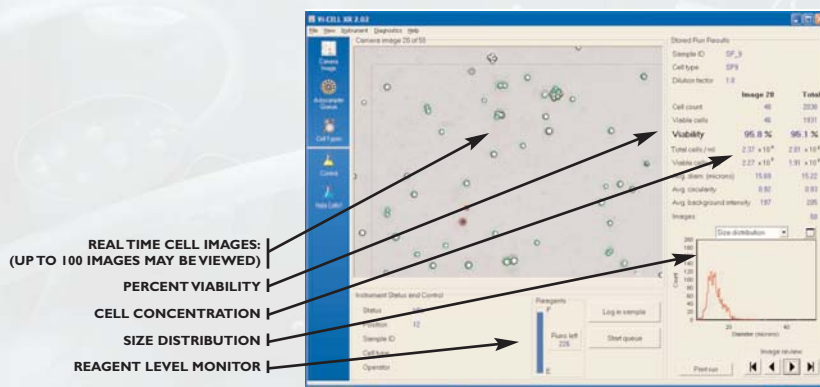


Figure 1. Viable cells are circled in green and non-viable in red.

## THE VI-CELL

The Vi-CELL automates the widely accepted Trypan Blue Dye Exclusion method. The Vi-CELL combines the state of the art in imaging technology, proprietary algorithm and fluidics management. At the heart of the Vi-CELL is the customized liquid handling system. This system, which allows sample aspiration, reagent handling and subsequent instrument cleaning, is fully automated. Once the cellular suspension has been aspirated and mixed with the trypan blue dye, it is pumped to the flow cell for imaging. The Vi-CELL can analyze up to 100 images for a given analysis increasing total volume from 15 to 30 times over the manual method with result in a less than 2.5 minutes.

## MANUAL TRYPAN BLUE DYE EXCLUSION METHOD

As mentioned, the standard method for measuring cell viability is the Trypan Blue Dye Exclusion method. Trypan blue stain (0.4%) is mixed with an equal volume of cells. Viable cells, given their intact membranes, exclude the trypan blue stain; non-viable cells, membrane permeable, stain dark blue. The manual method, however, requires a technician, using a hemacytometer and microscope, to enumerate both stained and unstained cells and manually calculate the percent viability. In addition to being labor intensive, this technique has substantial accuracy error due to its subjective nature.

- Prepare**
  - Automated liquid handling
  - Automated lysing
  - General purpose centrifugation
  - High performance centrifugation
  - Ultracentrifugation
- Identify**
  - Automated fluorescence microscopy
  - Cell counting
  - Cell markers
  - Cell viability analysis
  - Flow cytometry
  - Monoclonal antibodies
- Probe**
  - Automated liquid handling
  - Flow cytometry
  - Microarray technology
  - Monoclonal antibodies
  - Signal transduction assays
- Sort**
  - Cell sorters
  - Micro-piezo electric tips
  - Reagents (various)
- Evaluate**
  - Monoclonal antibodies
  - Multi-mode plate reading
  - Genomics solutions
  - Proteomics solutions
  - Software informatics
- Diagnose**
  - Automated liquid handling
  - Flow cytometry
  - Immunoassays
  - Monoclonal antibodies
  - Software algorithms

## VI-CELL TECHNICAL SPECIFICATIONS

### INSTRUMENT FUNCTION:

Concentration Range:  
 $5 \times 10^4$  to  $1 \times 10^7$  cells / mL  
 \*Counting Accuracy:  $\pm 6\%$

### OPERATING SYSTEM:

Windows® 98  
 Windows® 2000  
 Windows® XP

### INSTRUMENT TYPE:

Video imaging through a quartz flow cell

### POWER REQUIREMENTS:

Power 50 watts  
 (65 Watts Max.)  
 Voltages 100V, 120V,  
 220V or 240V 50/60 Hz

### TEMPERATURE:

10° to 40° C (50° to 104° F)

### WEIGHT:

11.3kg (25lb)

### UNIT DIMENSIONS:

44.5cm (17.5") height  
 38cm (15") width  
 41cm (16") depth

## VI-CELL SERIES

	PN	AUTO SAMPLE	SIZE RANGE (µm)	SAMPLE VOLUME (mL)	ANALYSIS TIME (Min)	VIABILITY RANGE	IMAGING TECHNOLOGY
VI-CELL XR	383556	Yes	2-70	0.5	<2.5	0-100	Auto-focus routine Firewire Camera 1394 X 1040 CCD array
VI-CELL AS	6605769	Yes	5-70	1.0	<3.5	0-100	Manual focus routine Image frame grabber 640 X 480 CCD array
VI-CELL S	383080	No	5-70	1.0	<3.5	0-100	Manual focus routine Image frame grabber 640 X 480 CCD array
VI-CELL XR QUAD PACK	383722						
VI-CELL AS, S QUAD PACK	383198						
VI-CELL CONCENTRATION CONTROL	175478						
VI-CELL FOCUS CONTROL	175474						



*Developing innovative solutions in  
Systems Biology.*

**Innovate** **Automate**  
SIMPLIFY

Beckman Coulter, Inc. • 4300 N. Harbor Boulevard, Box 3100 • Fullerton, California 92834-3100  
 Sales & Service: 1-800-742-2345 • Telex: 678413 • Fax: 1-800-643-4366 • www.beckmancoulter.com

### Worldwide Biomedical Research Division Offices:

Australia (61) 2 9844-6000 Canada (905) 819-1234 Caribbean and South America 1-305-380-4709 China (86) 10 6515 6028  
 Eastern Europe, Middle East, North Africa (41) 22 994 07 07 France 01 49 90 90 00 Germany 49 21 513335 Hong Kong (852) 2814 7431/2814 0481  
 Italy 02-953921 Japan 03-5404-8359 Mexico and Central America (52) 55-560 57770 Netherlands 0297-230630 Singapore (65) 6339 3633  
 South Africa, Sub-Saharan Africa (27) 11-805-2014/5 Spain 91 3836080 Sweden 08-564 85 900 Switzerland 0800 850 810  
 Taiwan (886) 2 2378 3456 Turkey 90 216 309 1900 U.K. 01494 441181 U.S.A. 1-800-742-2345