

Genesis Cell Isolation System with Celselect Slides™

**Capture more.
Recover more.
Accomplish more.**

Add new depth to rare cell studies with the efficiency, ease, and versatility of Celselect Slides.

The process of capturing cells — particularly those that are rare, treated, or part of a mixed population — is essential for research into their structure, function, and diversity.

The Genesis Cell Isolation System, using Celselect Slides, offers gentle, efficient size-based cell capture without the need for sample preprocessing, making it an ideal solution for capturing rare and circulating tumor cells from liquid biopsies and various other liquid samples.



Celselect Slides Enrichment Kit 2.0



Kit for cell capture and recovery for downstream applications

Celselect Slides Enumeration Indirect Stain Kit 2.0



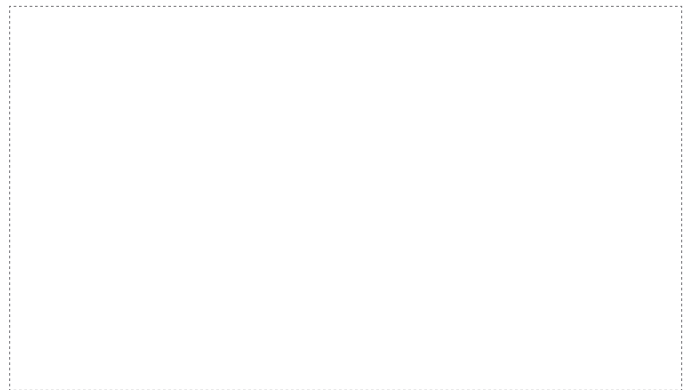
Kit for cell capture, staining via immunolabeling with secondary antibody, and on-slide imaging

Celselect Slides Enumeration Direct Stain Kit 2.0

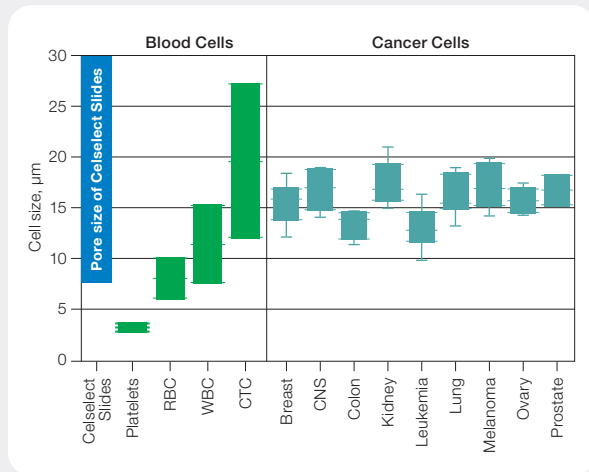


Kit for cell capture, staining via immunolabeling with primary antibody, and on-slide imaging

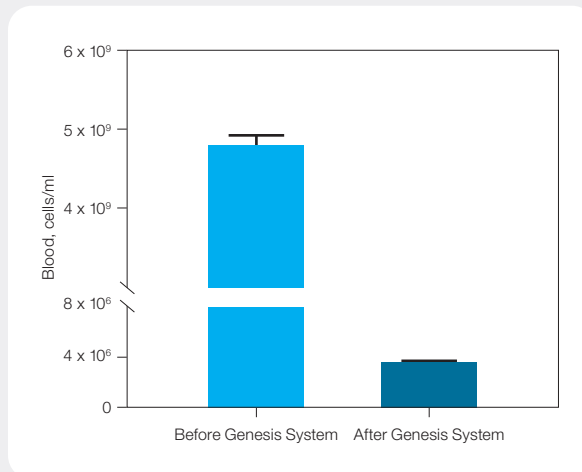
- Efficient capture of cells (8–30 μm)
- Smaller components (<8 μm) flow through slide
- Sample-type flexibility
- On-slide imaging ready
- Automated processing of up to 10 ml of sample



Rare Cell Capture Simplified

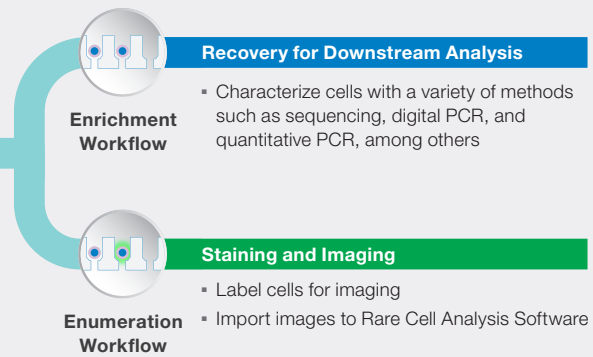
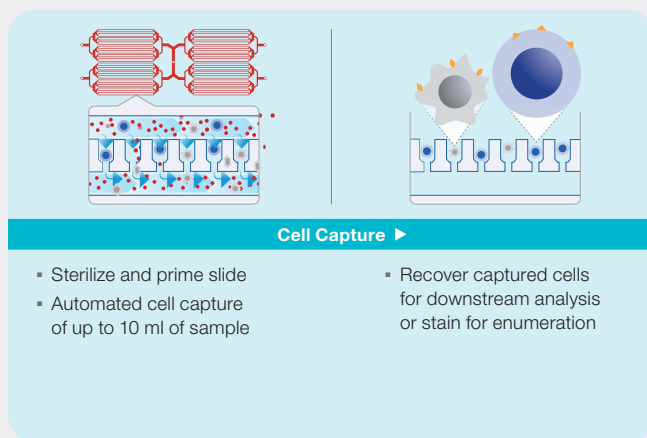


Pore size of Celselect Slides is suitable for diverse cell capture. CNS, central nervous system; CTC, circulating tumor cells; RBC, red blood cells; WBC, white blood cells.



Efficient blood cell removal: 99.9% elimination of red blood cells, white blood cells, platelets, and other components.

One System, Two Validated Workflows



Ordering Information

Catalog #	Description
12019822	Genesis Cell Isolation System , for automated capture of cells by size
17009455	Celselect Slides Enrichment Kit 2.0 , for cell capture and recovery for downstream applications
17009802	Celselect Slides Enumeration Indirect Stain Kit 2.0 , for cell capture and staining via labeled secondary antibody
17009841	Celselect Slides Enumeration Direct Stain Kit 2.0 , for cell capture and staining via labeled primary antibody



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

BIO-RAD and CELSELECT SLIDES 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.



**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

