

Ready·Sep·Go

Fast and Easy Isolation Of B Cells



 **STEMCELL™**
TECHNOLOGIES

Table of Contents

- 3 Cutting-Edge B Cell Research
- 4 Isolate B Cells In As Little As 25 Minutes
- 5 Research and Applications
- 6 EasySep™: Immunomagnetic B Cell Isolation Without Columns
- 6 RoboSep™: Fully Automated B Cell Isolation
- 7 RosetteSep™: One-Step Untouched B Cell Isolation From Whole Blood
- 8 Human B Cells: Starting with Whole Blood
- 9 Human B Cells: Starting with PBMC or Leukapheresis Samples
- 10 Human Naïve and Memory B Cells: Starting with PBMC or Leukapheresis Samples
- 11 Human CD138⁺ Plasma Cells: Starting with PBMC, Whole Blood or Bone Marrow
- 12 Mouse B Cells: Starting with Spleen or Other Tissues
- 13 Related Products
- 14 Product Listing
- 15 References

Cutting-Edge B Cell Research

Begins With Fast And Easy B Cell Isolation

B cells express cell surface immunoglobulin receptors that recognize specific antigenic epitopes, and are a pivotal component of the adaptive immune system and in particular, humoral immune responses. B cells mediate many processes necessary for immune homeostasis including antibody production, antigen presentation, cytokine secretion, T cell co-stimulation, and tumor immunity. Conversely, their dysregulation is the basis of several immune pathologies including autoimmunity, leukemia / lymphoma, and multiple myeloma.

Isolation of highly purified B cells is the first step towards successful B cell research. Isolate whole B cell populations as well as various B cell subsets directly from human whole blood, buffy coat, leukapheresis samples, or PBMC, and from mouse spleen or other tissues, with high purity and recovery using the fast and easy B cell isolation kits from STEMCELL Technologies.

With the unique cell separation platforms **EasySep™**, **RosetteSep™** and **RoboSep™**, STEMCELL Technologies is poised to support your immunological research endeavors.

Ready·Sep·Go

Isolate B Cells In As Little As 25 Minutes

Isolate whole B cell populations or specific B cell subsets in as little as 25 minutes using a B cell separation kit from our complete range.

- Whole B cell populations
- Naïve B cells
- Memory B cells
- CD138⁺ plasma cells
- B cells for CLL and multiple myeloma research
- B cells for HLA analysis
- Customized B cell isolations



Isolate B Cells In As Little As 25 Minutes

With High Purity And Preserved Viability

Fast and Easy

Magnetic or immunodensity separation without columns

High Purity and Recovery

Get high recovery, and purity of up to 99% pure B cells

B Cells are Immediately Ready for Use

Magnetic particles do not interfere with downstream applications such as flow cytometry

B Cells Retain Functionality

Isolated B cells have been used in various scientific investigations, including publications that:

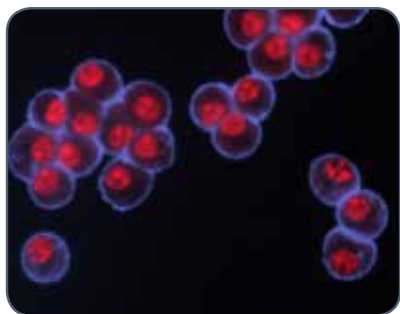
- Identify intermediate cell populations in naïve B cell development¹
- Investigate modulators of human B cell differentiation^{2,3}
- Demonstrate the role of innate immunity signaling pathways in B cell tolerance⁴
- Characterize the mechanism of plasma and memory B cell selection⁵
- Evaluate gene expression in B cell subsets⁶



Research and Applications

Isolated B Cells Are Suitable For All Downstream Analysis

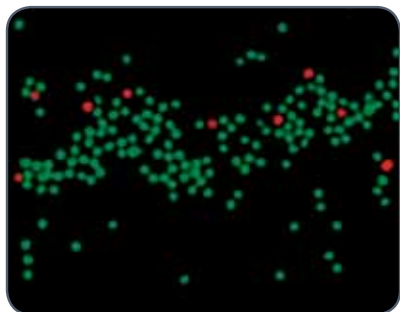
The gentle, column-free cell separation platforms **RosetteSep™**, **EasySep™** and **RoboSep™** each preserve B cell viability, ensuring that purified B cells are suitable for functional and biological studies. Both enriched and positively selected B cells are immediately ready for downstream assays. Negatively enriched B cells are untouched by EasySep™ reagents. Positively selected B cells are labeled with small **EasySep™** magnetic particles that are flow cytometry compatible. B cells can be routinely and reliably isolated at up to 99% purity, making them highly complementary to molecular diagnostic techniques used for hematologic malignancy analysis.



B Cell Functionality

Isolated B cells are highly functional and have been used for:

- Lipid antigen presentation to NKT cells⁷
- Intracellular calcium flux assays⁸
- Hybridoma generation⁹
- Adoptive transfer¹⁰



HLA Testing

The HLA B cell isolation products are specifically optimized for isolating B cells for HLA testing purposes and have been formulated for:

- Serology-based assays
- Flow Cytometry Crossmatch (FCXM)
- Chimerism analysis



Hematologic Malignancies

Purifying malignant B cells can greatly improve the accuracy and reliability of molecular diagnostic studies. Isolate plasma cells and B cells for the study of:

- Multiple Myeloma
- Chronic Lymphocytic Leukemia

EasySep™

Immunomagnetic B Cell Isolation Without Columns

EasySep™ is a fast and easy way to get highly purified B cells from whole blood, buffy coat, fresh or previously frozen peripheral blood mononuclear cells (PBMC), leukapheresis samples, and mouse spleen or other tissue samples. Cells are crosslinked to EasySep™ magnetic particles using the Tetrameric Antibody Complex (TAC) technology, and then separated from unwanted cells with an EasySep™ magnet without the use of columns.

EasySep™ Human B Cell Enrichment Procedure

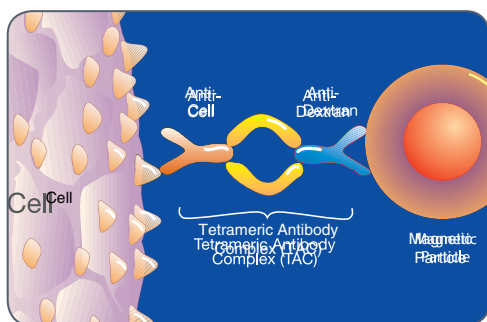
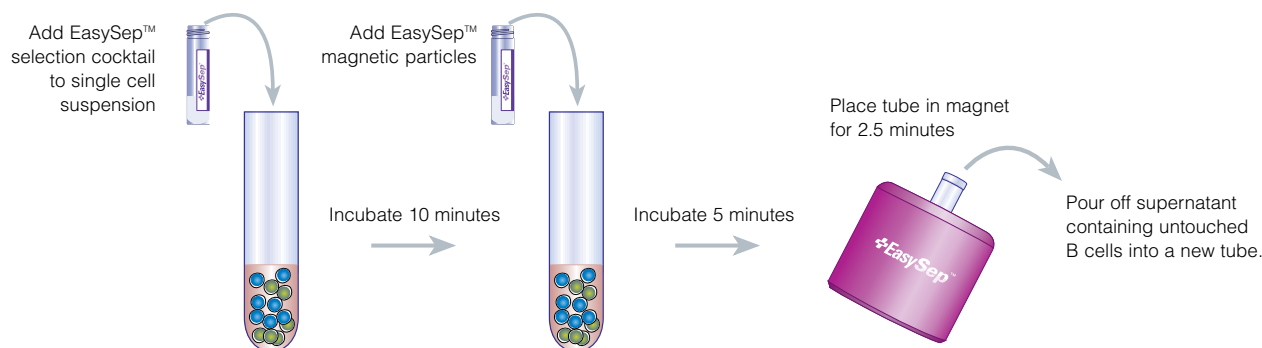


FIGURE 1. Tetrameric Antibody Complex (TAC) crosslinking a cell to a dextran coated magnetic particle.

VIDEO

EasySep™: Powerful Immunomagnetic Isolation of Virtually Any Cell Type

www.stemcell.com/EasySepVideo

SCAN ME ►



RoboSep™

Fully Automated B Cell Isolation



Streamline your B cell separations using RoboSep™, the fully automated cell separator. By performing all EasySep™ cell labeling and separation steps, RoboSep™ maintains the speed and simplicity of EasySep™ while offering walk-away automation, enabling the high-throughput, versatile isolation of highly purified cells. Minimize sample handling and eliminate cross-contamination while isolating B cells with just 5 minutes of “hands-on” technician time.

RosetteSep™

One-Step Untouched Human B Cell Isolation From Whole Blood

RosetteSep™ is a rapid immunodensity procedure for the isolation of untouched B cells directly from whole blood. It eliminates the need for a separate magnetic separation step by isolating cells during a standard density centrifugation Ficoll™ step, significantly reducing handling time and maximizing convenience. RosetteSep™ crosslinks unwanted cells to red blood cells present in the sample to form immunorosettes. When centrifuged over a density medium, the unwanted cells pellet along with the red blood cells, leaving desired cells at the interface between the plasma and the density medium.

RosetteSep™ Human B Cell Enrichment Procedure

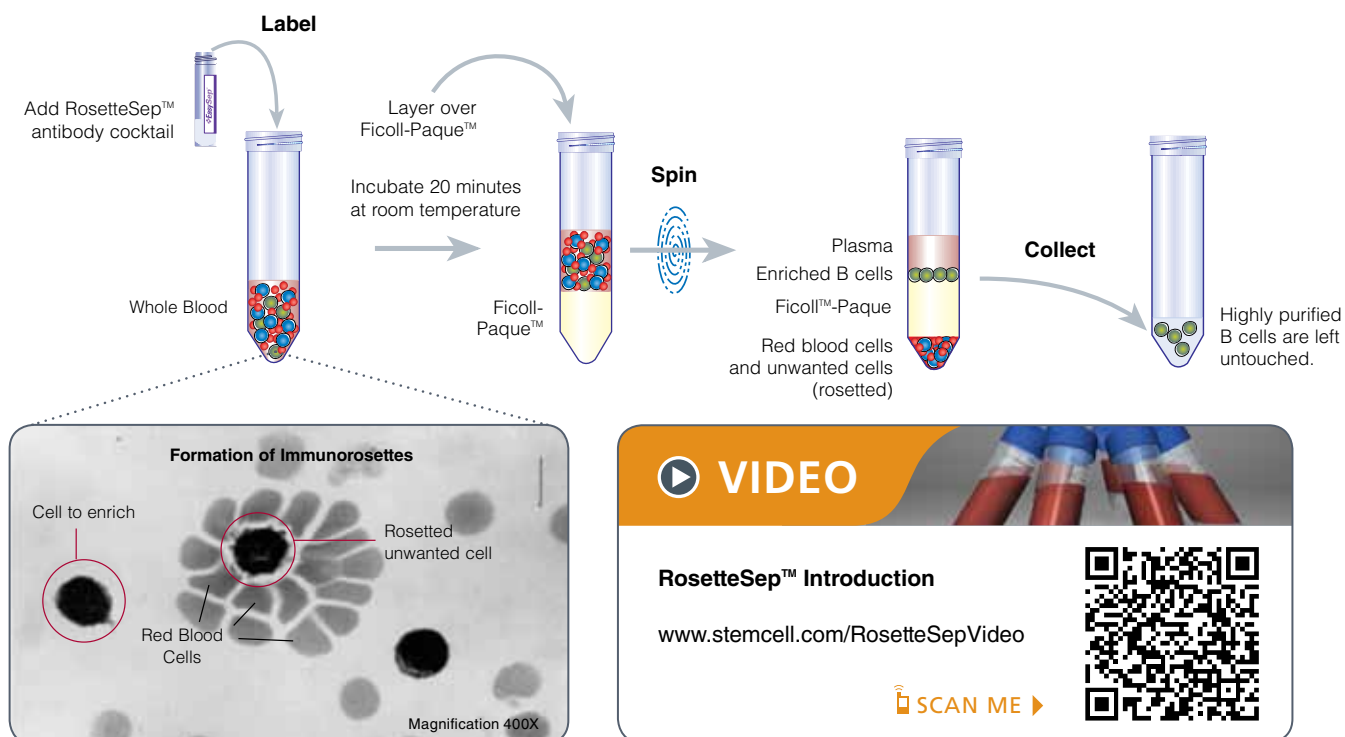


FIGURE 2. Image of a blood sample after addition of the RosetteSep™ cocktail, and prior to centrifugation over Ficoll™.

Did You Know?

Although RosetteSep™ has been optimized for use with whole blood, cells can be enriched from other sources (for example, buffy coat or leukapheresis samples.) The concentration of nucleated cells in the sample should not exceed 5×10^7 cells/mL, and red blood cells (RBCs) should be present at a ratio of at least 30 - 50 RBCs per nucleated cell.

Ficoll™ and Ficoll-Paque™ are trademarks of GE Healthcare Ltd.

Human B Cells

Starting With Whole Blood

Isolate untouched human B cells by negative selection, or positively select CD19⁺, CD20⁺, or CD19⁺/CD20⁺ cells directly from whole blood.

RosetteSep™ Human B Cell Enrichment Cocktail (Catalog #15024)

Isolate untouched human B cells directly from whole blood without magnets or special equipment.

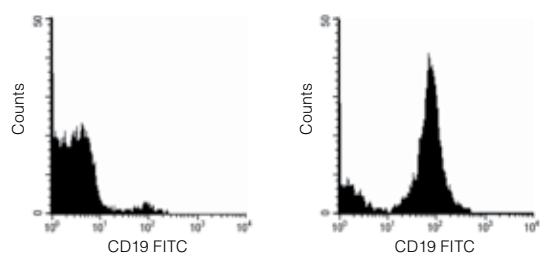
HOW IT WORKS:

The B cell population is enriched by negative selection using RosetteSep™, which depletes unwanted cells and leaves the desired B cells untouched by antibody.

Typical FACS Histogram Results with RosetteSep™ Human B Cell Enrichment Cocktail

Start: 6% CD19⁺ cells

Enriched: 81% CD19⁺ cells



Starting with fresh whole blood, the CD19⁺ cell content of the enriched fraction typically ranges from 81 - 83%.

This cocktail has been used to isolate B cells for:

- Intracellular calcium flux assays⁸
- Immunofluorescence and FISH analysis¹¹
- Chromatin immunoprecipitation analysis¹¹

EasySep™ Human Whole Blood CD19 Positive Selection Kit (Catalog #18084)

Isolate CD19⁺ human B cells directly from whole blood.

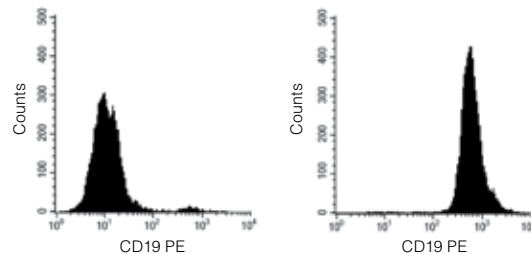
HOW IT WORKS:

CD19⁺ B cells are isolated by positive selection using EasySep™.

Typical FACS Histogram Results with EasySep™ Human Whole Blood CD19 Positive Selection Kit.

Start: 2% CD19⁺ cells

Selected: 99% CD19⁺ cells



Starting with fresh whole blood, the CD19⁺ cell content of the selected fraction typically ranges from 94 - 99%.

This kit has been used to isolate B cells for:

- Hybridoma generation⁹
- Therapeutic monoclonal antibody production⁹
- Hybridoma optimization using morphogenics⁹

Also Available

EasySep™ Human Whole Blood CD20 Positive Selection Kit (Catalog #18685)

Isolate CD20⁺ B cells directly from whole blood. Starting with fresh whole blood, the CD20⁺ cell content of the selected fraction typically ranges from 96 - 99%.

B Cell Kits for HLA Analysis

See page 14.

Human B Cells

Starting With PBMC Or Leukapheresis Samples

Isolate untouched human B cells (with or without CD43 depletion) by negative selection directly from PBMC or leukapheresis samples.

EasySep™ Human B Cell Enrichment Kit (Catalog #19054)

Isolate untouched B cells from PBMC or leukapheresis samples.

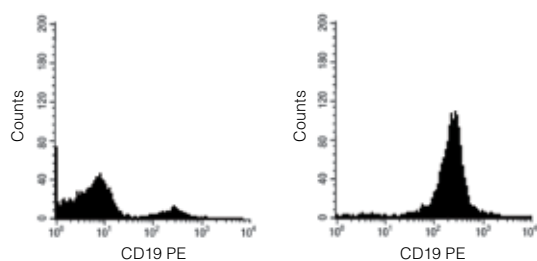
HOW IT WORKS:

The B cell population is enriched by negative selection using EasySep™, which depletes unwanted cells and leaves desired B cells untouched by antibody.

Typical FACS Histogram Results with EasySep™ Human B Cell Enrichment Kit

Start: 8% CD19⁺ cells

Enriched: 99% CD19⁺ cells



Starting with fresh or previously frozen PBMC, the CD19⁺ cell content of the enriched fraction typically ranges from 95 - 99%.

This kit has been used to isolate B cells for:

- Lipid antigen presentation to NKT cells⁷
- DNase protection assays¹²
- Gene expression profiling by microarray analysis¹³

EasySep™ Human B Cell Enrichment Kit without CD43 Depletion (Catalog #19154)

CLL
Research
Compatible

Isolate untouched human B cells from PBMC samples that include B cells that express CD43 (e.g. CLL samples).

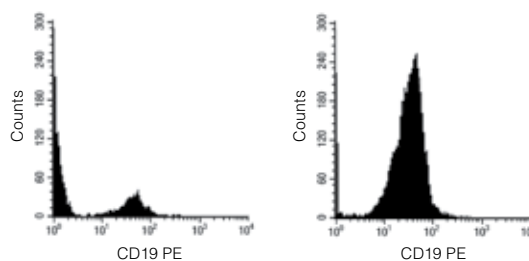
HOW IT WORKS:

The B cell population, including CD43⁺ B cells, is enriched by negative selection using EasySep™, which depletes unwanted cells and leaves desired B cells untouched by antibody.

Typical FACS Histogram Results with EasySep™ Human B Cell Enrichment Kit (without CD43 Depletion)

Start: 11% CD19⁺ cells

Enriched: 96% CD19⁺ cells



Starting with fresh or previously frozen PBMC, the CD19⁺ cell content of the enriched fraction typically ranges from 87 - 98%.

This kit has been used to isolate B cells for:

- Quantitative rt-PCR analysis in Chronic Lymphocytic Leukemia (CLL)¹⁴
- Analyzing the specificity of cytotoxic drugs for CLL¹⁵
- Gene expression profiling in CLL¹⁶

Also Available

EasySep™ Human CD19 Positive Selection Kit (Catalog #18054)

Isolate CD19⁺ B cells from PBMC. Starting with fresh or previously frozen PBMC, the CD19⁺ cell content of the selected fraction typically ranges from 97 - 99%.

B Cell Kits for HLA Analysis

See page 14.

Human Naïve and Memory B Cells

Starting with PBMC or Leukapheresis Samples

Isolate specific human B cell subsets from PBMC or leukapheresis samples. Isolate untouched naïve B cells by negative selection or positively select CD27⁺ memory B cells from a pre-enriched B cell population.

EasySep™ Human Naïve B Cell Enrichment Kit (Catalog #19254)



Isolate untouched human naïve B cells from PBMC or leukapheresis samples.

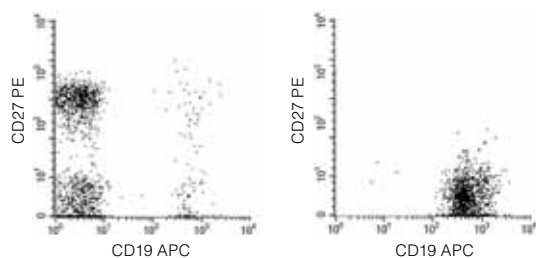
HOW IT WORKS:

The naïve B cell population is enriched by negative selection using EasySep™ which depletes unwanted cells and leaves desired cells untouched by antibody.

Typical FACS Profile Results with EasySep™ Human Naïve B Cell Enrichment Kit

Start:
4% CD19⁺ CD27⁻ cells

Enriched:
97% CD19⁺ CD27⁻ cells



Starting with fresh or previously frozen PBMC, the CD19⁺CD27⁻ cell content of the enriched fraction typically ranges from 92 - 98%.

EasySep™ Human Memory B Cell Isolation Kit (Catalog #18164)



Isolate human memory B cells from PBMC or leukapheresis samples.

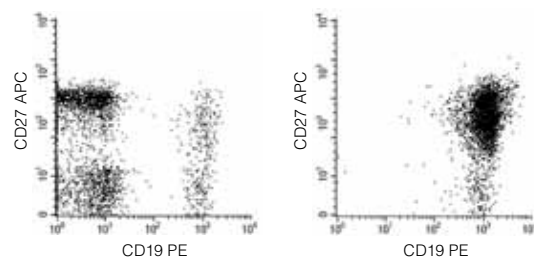
HOW IT WORKS:

The B cell population is first enriched by negative selection. CD27⁺ memory B cells are then isolated using EasySep™ CD27 positive selection.

Typical FACS Profile Results with EasySep™ Human Memory B Cell Isolation Kit

Start:
6% CD19⁺ CD27⁺ cells

Selected:
92% CD19⁺ CD27⁺ cells



Starting with fresh or previously frozen PBMC, the CD19⁺CD27⁺ cell content of the enriched fraction typically ranges from 85 - 95%.

Did You Know?

We also provide comprehensive protocols on how to isolate highly purified naïve and memory B cells from the same sample. Contact techsupport@stemcell.com for a copy.

Human CD138⁺ Plasma Cells

Starting with PBMC, Whole Blood or Bone Marrow

Isolate CD138⁺ (normal and malignant) plasma cells directly from bone marrow, PBMC, or whole blood.

EasySep™ Human CD138 Positive Selection Kit (Catalog #18357)



For Multiple
Myeloma
Research

Isolate CD138⁺ plasma cells from PBMC or bone marrow.

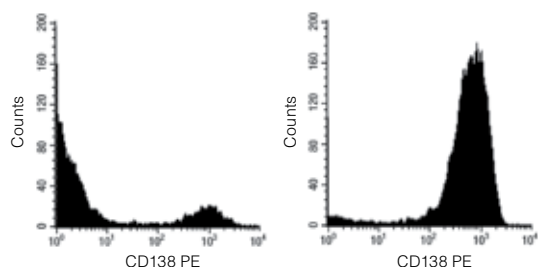
HOW IT WORKS:

CD138⁺ (Syndecan-1) cells are isolated by positive selection using EasySep™. CD138 is expressed on normal or malignant plasma cells, but not on mature B cells. Isolated cells are highly suited for multiple myeloma research.

Typical FACS Histogram Results with EasySep™ Human CD138 Positive Selection Kit

Start: 10% CD138⁺ cells

Selected: 91% CD138⁺ cells



Starting with fresh or previously frozen PBMC, the CD138⁺/Syndecan-1 cell content of the selected fraction typically ranges from 85 - 95%. Please note that purity is highly dependent on the starting sample.

EasySep™ Human Whole Blood CD138 Positive Selection Kit (Catalog #18387)



For Multiple
Myeloma
Research

Isolate CD138⁺ plasma cells from whole blood.

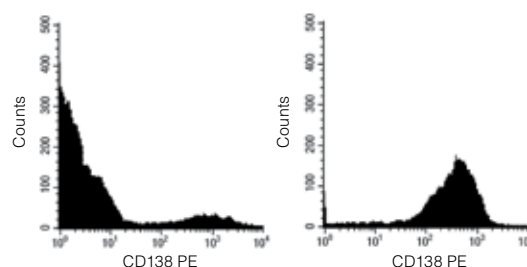
HOW IT WORKS:

CD138⁺ (Syndecan-1) cells are isolated by positive selection using EasySep™. CD138 is expressed on normal or malignant plasma cells, but not on mature B cells. Isolated cells are highly suited for multiple myeloma research.

Typical FACS Histogram Results with EasySep™ Human Whole Blood CD138 Positive Selection Kit

Start: 6% CD138⁺ cells*

Selected: 94% CD138⁺ cells



Starting with fresh whole blood spiked with a multiple myeloma cell line, the CD138⁺/Syndecan-1 cell content of the selected fraction typically ranges from 89 - 98%.

*Residual red blood cells were removed by lysis prior to flow cytometry.

These kits have been used to isolate plasma cells for research of multiple myeloma¹⁷, primary systemic amyloidosis¹⁸ and plasma cell leukemia¹⁹.

Also Available

RosetteSep™ Human Multiple Myeloma Cell Enrichment Cocktail (Catalog #15129)

Isolate untouched multiple myeloma cells (B cells and plasma cells) from fresh human bone marrow aspirates by negative selection, without the use of magnets or special equipment.

Mouse B Cells

Starting With Spleen Or Other Tissues

Isolate untouched mouse B cells by negative selection, or positively select CD19⁺ B cells directly from single cell suspensions of spleen or other tissues.

EasySep™ Mouse B Cell Enrichment Kit (Catalog #19754)

Isolate untouched mouse B cells from single cell suspensions of spleen or other tissues.

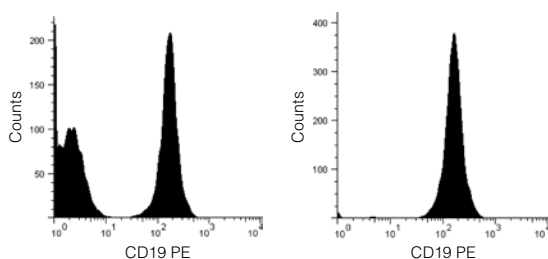
HOW IT WORKS:

The B cell population is enriched by negative selection using EasySep™ which depletes unwanted cells and leaves desired cells untouched by antibody.

Typical FACS Histogram Results with EasySep™ Mouse B Cell Enrichment Kit

Start: 50% CD19⁺ cells

Enriched: 96% CD19⁺ cells



This kit has been used to isolate primary B cells for:

- Protein extraction and immunoblotting^{20,21}
- Analysis of class-switching recombination^{22,23}
- Retroviral transduction²²

EasySep™ Mouse CD19 Positive Selection Kit (Catalog #18754)

Isolate CD19⁺ mouse B cells from single cell suspensions of spleen or other tissues.

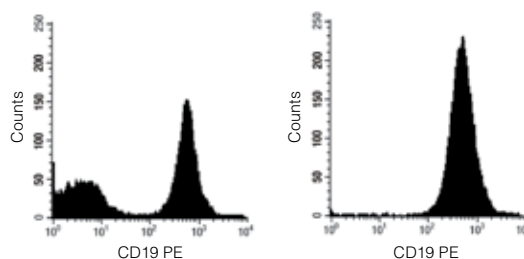
HOW IT WORKS:

CD19⁺ B cells are isolated by positive selection using EasySep™.

Typical FACS Histogram Results with EasySep™ Mouse CD19 Positive Selection

Start: 54% CD19⁺ cells

Selected: 99% CD19⁺ cells



This kit has been used to isolate primary B cells for:

- Co-culture with tumor cells²⁴
- Quantitative RT-PCR²⁵
- Adoptive Transfer¹⁰

Related Products

Equipment

PRODUCTS	DESCRIPTION	CATALOG #
EasySep™ Magnet	Immunomagnetic Column-Free Magnet (for 5 mL tubes)	18000
"The Big Easy" EasySep™ Magnet	Immunomagnetic Column-Free Magnet (for 14 mL tubes)	18001
"Easy50" EasySep™ Magnet	Immunomagnetic Column-Free Magnet (for 50 mL tubes)	18002
EasySep™ Multistand	Holds up to 4 EasySep™ or 4 "The Big Easy" EasySep™ Magnets for separating up to 4 samples at once	18010
	18010 with 4 EasySep™ Magnets	18004
	18010 with 4 "The Big Easy" EasySep™ Magnets	18100
"EasyPlate" EasySep™ Magnet	Immunomagnetic Column-Free Magnet (holds a 96-well plate)	18102
RoboSep™	The Fully Automated Cell Separator	20000

Antibodies

PRODUCTS	DESCRIPTION	CATALOG #
CD19 Antibody, Clone 4G7, PE-Conjugated	Mouse Monoclonal Antibody to Human CD19 - PE Conjugated	10509
CD19 Antibody, Clone 4G7, FITC-Conjugated	Mouse Monoclonal Antibody to Human CD19 - FITC Conjugated	10409
CD20 Antibody, Clone L27, PE-Conjugated	Mouse Monoclonal Antibody to Human CD20 - PE Conjugated	10510
CD20 Antibody, Clone L27, FITC-Conjugated	Mouse Monoclonal Antibody to Human CD20 - FITC Conjugated	10410

Product Listing

Complete Product Listing

SPECIES	STARTING SAMPLE	B CELL PHENOTYPE	PRODUCT	CATALOG #
Human	Whole blood	All B cells	RosetteSep™ Human B Cell Enrichment Cocktail	15024
		Multiple myeloma cells	RosetteSep™ Human Multiple Myeloma Cell Enrichment Cocktail	15129
		CD19 ⁺	EasySep™/RoboSep™ Human Whole Blood CD19 Positive Selection Kit	18084/18084RF
		CD20 ⁺	EasySep™/RoboSep™ Human Whole Blood CD20 Positive Selection Kit	18685/18685RF
		CD138 ⁺ (plasma cells)	EasySep™/RoboSep™ Human Whole Blood CD138 Positive Selection Kit	18387/18387RF
	PBMC	All B cells	EasySep™/RoboSep™ Human B Cell Enrichment Kit	19054/19054RF
		All B cells without CD43 depletion	EasySep™/RoboSep™ Human B Cell Enrichment Kit without CD43 Depletion	19154/19154RF
		Naïve B cells	EasySep™/RoboSep™ Human Naïve B Cell Enrichment Kit	19254/19254RF
		Memory B cells	EasySep™/RoboSep™ Human Memory B Cell Isolation Kit	18164/18164RF
		CD19 ⁺	EasySep™/RoboSep™ Human CD19 Positive Selection Kit	18054/18054RF
		CD138 ⁺ (plasma cells)	EasySep™/RoboSep™ Human CD138 Positive Selection Kit	18357/18357RF
	Custom	EasySep™/RoboSep™ Human Custom Enrichment Kit		19309/19309RF
		EasySep™/RoboSep™ Human Custom Positive Selection Kit		18309/18309RF
Mouse	Spleen or other tissue	All B cells	EasySep™/RoboSep™ Mouse B Cell Enrichment Kit	19754/19754RF
		CD19 ⁺	EasySep™/RoboSep™ Mouse CD19 Positive Selection Kit	18754/18754RF
	Custom	EasySep™/RoboSep™ Mouse Custom Enrichment Kit		19709/19709RF
		EasySep™/RoboSep™ Mouse Custom Positive Selection Kit		18709/18709RF

B Cell Kits for HLA Analysis

STARTING SAMPLE	B CELL PHENOTYPE	PRODUCT	CATALOG #
Whole blood	All B cells	RosetteSep™ HLA B Cell Enrichment Cocktail	15064HLA
		EasySep™/RoboSep™ HLA B Cell Enrichment: Complete Processing Kit for Whole Blood	19954HLA/19954HLARF
	CD19 ⁺ /CD20 ⁺	EasySep™/RoboSep™ HLA Whole Blood B Cell Positive Selection Kit	18184HLA/18184HLARF
	CD19 ⁺	EasySep™/RoboSep™ Human Whole Blood CD19 Positive Selection Kit	18084/18084RF
	CD20 ⁺	EasySep™/RoboSep™ Human Whole Blood CD20 Positive Selection Kit	18685/18685RF
PBMC	All B cells	EasySep™/RoboSep™ HLA B Cell Enrichment Kit	19054HLA/19054HLARF
	CD19 ⁺ /CD20 ⁺	EasySep™/RoboSep™ HLA B Cell Positive Selection Kit	18454HLA/18454HLARF
	CD19 ⁺	EasySep™/RoboSep™ Human CD19 Positive Selection Kit	18054/18054RF

References

1. Lee J, Kuchen S, Fischer R et al. Identification and characterization of a human CD5⁺ pre-naïve B cell population. *J. Immunol.* 182:4116-4126, 2009
2. Chen S, Sims G, Chen X et al. Modulatory effects of 1,25-Dihydroxyvitamin D3 on human B cell differentiation. *J. Immunol.* 179:1634-1647, 2007
3. Darce J, Arendt B, Chang SK et al. Divergent Effects of BAFF on human memory B cell differentiation into Ig-secreting cells. *J. Immunol.* 178:5612-5622, 2007
4. Isnardi I, Ng Y, Srdanovic I et al. IRAK-4- and MyD88-dependent pathways are essential for the removal of developing autoreactive B cells in humans. *Immunity* 29:746-757, 2008
5. Good-Jacobson K, Szumilas C, Chen L et al. PD-1 regulates germinal center B cell survival and the formation and affinity of long-lived plasma cells. *Nat. Immunol.* 11:535-543, 2010
6. Girschick H, Grammer A, Nanki T et al. RAG1 and RAG2 expression by B cell subsets from human tonsil and peripheral blood. *J. Immunol.* 166:377-386, 2001
7. Dumont N, Aubin E, Proul D et al. Increased secretion of hyperimmune antibodies following lipopolysaccharide stimulation of CD40-activated human B cells in vitro. *Immunology* 126:588-595, 2009
8. Isnardi I, Ng Y, Menard L et al. Complement receptor 2/CD21⁺ human naïve B cells contain mostly autoreactive unresponsive clones. *Blood* 115:5026-5036, 2010
9. Li J, Sai T, Berger M et al. Human antibodies for immunotherapy development generated via a human B cell hybridoma technology. *PNAS* 103:3557-3562, 2006
10. Gligley J, Fox B, and Bzik D. Cell-mediated immunity to toxoplasma gondii develops primarily by local Th1 host immune responses in the absence of parasite replication. *J. Immunol.* 182:1069-1078, 2009
11. Brugat T, Nguyen-Khac F, Grelrier A et al. Telomere dysfunction-induced foci arise with the onset of telomeric deletions and complex chromosomal aberrations in resistant chronic lymphocytic leukemia cells. *Blood* 116:239-249, 2010
12. Mihalcik S, Huddleston P, Wu X et al. The structure of the TNFRSF13C promoter enables differential expression of BAFF-R during B cell ontogeny and terminal differentiation. *J. Immunol.* 185:1045-1054, 2010
13. Longo N, Lugar P, Yavuz S et al. Analysis of somatic hypermutation in X-linked hyper-IgM syndrome shows specific deficiencies in mutational targeting. *Blood* 113:3706-3715, 2009
14. Asslaber D, Piñón J, Seyfried I et al. microRNA-34a expression correlates with MDM2 SNP309 polymorphism and treatment-free survival in chronic lymphocytic leukemia. *Blood* 115:4191-4197, 2010
15. Isham C, Tibodeau J, Jin W et al. Chaetocin: a promising new antimyeloma agent with in vitro and in vivo activity mediated via imposition of oxidative stress. *Blood* 109:2579-2588, 2007
16. Morabito F, Cutrona G, Gentile M et al. Definition of progression risk based on combinations of cellular and molecular markers in patients with Binet stage A chronic lymphocytic leukaemia. *Brit.J.of Haem.* 146:44-53, 2009
17. Clendening J, Pandya A, Li Z et al. Exploiting the mevalonate pathway to distinguish statin-sensitive multiple myeloma. *Blood* 115:4787-4797, 2010
18. Arendt B, Ramirez-Alvarado M, Sikkink L et al. Biologic and genetic characterization of the novel amyloidogenic lambda light chain-secreting human cell lines, ALMC-1 and ALMC-2. *Blood* 112:1931-1941, 2008
19. Ng P, Helguera G, Daniels T et al. Molecular events contributing to cell death in malignant human hematopoietic cells elicited by an IgG3-avidin fusion protein targeting the transferrin receptor. *Blood* 108:2745-2754, 2006
20. Rai D, Kim S, McKeller M et al. Targeting of SMAD5 links microRNA-155 to the TGF- β pathway and lymphomagenesis. *PNAS* 107:3111-3116, 2010
21. Kim S, Rai D, McKeller M et al. Rational combined targeting of phosphodiesterase 4B and SYK in DLBCL. *Blood* 113:6153-6160, 2009
22. Park S, Zan H, Pal Z et al. HoxC4 binds to the promoter of the cytidine deaminase AID gene to induce AID expression, class switch DNA recombination and somatic hypermutation. *Nat. Immunol.* 10:540-550, 2009
23. Wu X, Darce J, Chang SK et al. Alternative splicing regulates activation-induced cytidine deaminase (AID): implications for suppression of AID mutagenic activity in normal and malignant B cells. *Blood* 112:4675-4682, 2008
24. Inoue S, Leitner W, Golding B et al. Inhibitory effects of B cells on antitumor immunity. *Cancer Res.* 66:7741-7747, 2006
25. Makui H, Soares R, Jiang W. et al. Contribution of Hfe expression in macrophages to the regulation of hepatic hepcidin levels and iron loading. *Blood* 106:2189-2195, 2005

THE CELL EXPERTS™ | WWW.STEMCELL.COM

TOLL-FREE PHONE 1 800 667 0322 • PHONE 1 604 877 0713

TOLL-FREE FAX 1 800 567 2899 • FAX 1 604 877 0704

ORDERS@STEMCELL.COM • INFO@STEMCELL.COM

FOR FULL CONTACT DETAILS WORLDWIDE VISIT OUR WEBSITE

FOR RESEARCH USE ONLY. NOT FOR THERAPEUTIC OR DIAGNOSTIC USE. CATALOG #28786 VERSION 1.0.1 JULY 2012

