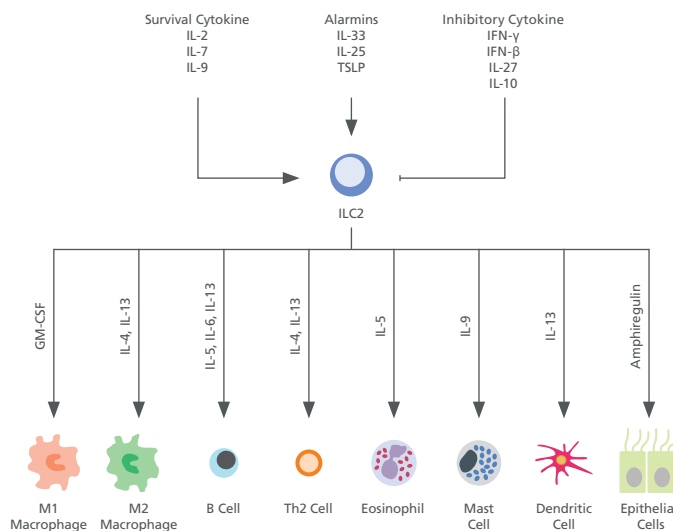


# Gating Strategy for the Identification of ILC2s

## Background

Studies starting in 2010, led several research groups to identify a small population of immune cells that have characteristics of lymphoid cells but lack lineage markers (Lin-) and re-arranged antigen specific cell surface receptors.<sup>1,2</sup> As these cells were identified they were referred to by different names such as natural helper cells,<sup>3,4</sup> nuocytes,<sup>5</sup> innate lymphoid cells<sup>6</sup> and innate helper 2 cells.<sup>7</sup> In 2013 a uniform nomenclature was proposed to classify these cells as group 2 innate lymphoid cells (ILC2s).<sup>8</sup> ILC2s are part of the larger ILC family, that according to this new nomenclature, is divided into three major groups: group 1 ILCs (ILC1s), group 2 ILCs (ILC2s) and group 3 ILCs (ILC3s).<sup>8</sup> The groupings are based on transcription factor expression, surface receptors and cytokine-secreting profiles that mirror the profiles of T helper subsets, Th1, Th2 and Th17 respectively. Accordingly, ILC2s produce Th2 associated cytokines including interleukin (IL)-4, IL-5, IL-9 and IL-13 and are involved in type 2 immune response (Figure 1). By surface staining, human ILC2s are negative for lineage markers and stain positive for CD45, CD294, CD127 and CD161. Mouse ILC2s are negative for lineage markers and positive for CD45, ICOS and ST2 (at varying levels). Since their initial identification, we now know that ILC2s are important effector cells implicated in innate immunity and are required for anti-helminth immunity, allergic inflammation and tissue repair.<sup>1</sup> More information on the development, biology and function of the distinct ILC subsets, including ILC2s, can be found in several publications.<sup>1,2,9,10</sup>



**Figure 1. Group 2 Innate Lymphoid Cells (ILC2s)**

ILC2s respond to a diverse range of stimuli including survival cytokines, such as IL-2, IL-7 and IL-9, and alarmins such as IL-33, IL-25 and TSLP. ILC2s then produce type-2 cytokines, including IL-4, IL-5, IL-9 and IL-13 and thereby promote activation of macrophages, B cells, CD4<sup>+</sup> T cells, basophils, eosinophils, mast cells, dendritic cells and epithelial cells. GM-CSF, granulocyte macrophage colony-stimulating factor; IFN<sub>γ</sub>, interferon-γ; IL, interleukin; TSLP, thymic stromal lymphopoietin.

## The Challenge

ILC2s are widely distributed throughout the body. They are enriched at barrier surfaces such as the skin, intestine and lung, but are also present in adipose tissue and peripheral blood.<sup>1</sup> Even though ILC2s are found in different tissues they comprise less than 1% of the total leukocyte population in human and mice.<sup>11,12</sup> The low frequency and lack of specific surface markers for ILC2s and ILCs in general, creates challenges when it comes to identification, isolation and further characterization of these cells. Currently, ILC2s are isolated using multicolor flow cytometric cell sorting but this method is time-consuming, expensive and can result in low cell recovery. To address this challenge, STEMCELL Technologies has developed products to enrich for ILC2 populations from human and mouse tissue using the RosetteSep™ and EasySep™ technology (see page 4). Pre-enrichment of ILC2s will reduce sorting time and improve purity and recovery. The enriched ILC2s can then be further prepared for analysis or sorting using flow cytometry.

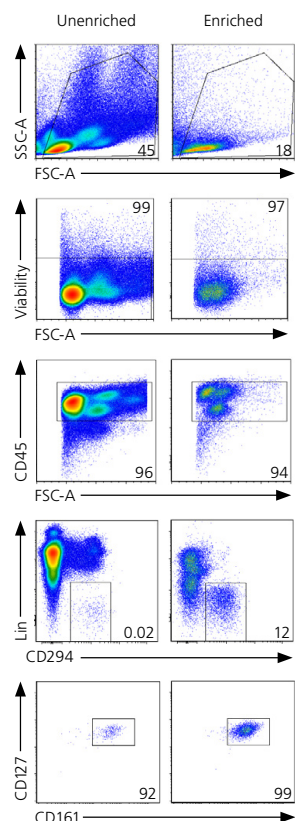
This Technical Bulletin Shows the Gating Strategy for Sorting or Analyzing Pre-enriched Human and Mouse ILC2s by Flow Cytometry.

## Gating Strategy for the Identification of Human and Mouse ILC2s

### Cell Pre-Enrichment and Antibody Labeling

Traditionally, to sort or analyze ILC2s by flow cytometry, one would begin by collecting the tissue of interest and prepare a single cell suspension using a suitable protocol. The entire sample is then labeled with the appropriate fluorochrome-conjugated antibodies and a viability dye. Recommended antibodies for labeling human and mouse cells are shown in Table 1 and Table 2, respectively. Note that both human and mouse ILC2s are defined as lineage-negative (Lin-) cells. In order to gate the Lin- cell population more easily, ensure that all antibodies recognizing lineage markers have the same fluorochrome (e.g. FITC). The user will then proceed to a flow cytometer and continue with the analysis or sorting of rare ILC2s.

Since ILC2s are present at a low frequency, we recommend performing a pre-enrichment step using EasySep™ or RosetteSep™ ILC2 enrichment kits (see page 4) before labeling cells with fluorochrome-conjugated antibodies. By removing the bulk of unwanted cells and pre-enriching for ILC2s, the user is able to process their sample more quickly while improving the recovery of ILC2s on a flow cytometer.



**Figure 2. Gating Strategy for ILC2s From Human Blood Samples.**

Enriched cells were obtained using RosetteSep™ Human ILC2 Enrichment Kit (Catalog #15382) as shown in page 4. For the unenriched sample, cells were isolated from human whole peripheral blood using density gradient centrifugation. The frequency of ILC2s (Lin-CD45+ CD294+ CD127+ CD161+) in the enriched fraction typically ranges from 0.44 - 53% and is donor dependent. In this example, the percentage of ILC2s in the unenriched and enriched fractions corresponds to 0.02 and 12%, respectively. FACS plots were generated using FCS Express 5 Software (De Novo Software).

### Identification of Human ILC2s from Blood

A representative gating strategy to identify human ILC2s from unenriched and enriched human whole blood samples is shown in Figure 2. The same gating strategy can be applied when analyzing cells from leukapheresis samples (not shown). Following the gating strategy shown on Figure 2, start on the SSC-A versus FSC-A plots and create a gate to include all leukocytes based on cell granularity and cell size. Alternatively, create a smaller gate around only the lymphocytes (not shown). When sorting cells, remove doublets and gate on single cells (singlets) using either the FSC-W versus FSC-H or SSC-W versus SSC-H plots (not shown). Then, select for viable cells and then for CD45+ cells. Next, create a gate on the Lin- CRTH2+ (CD294) population. Within this subset the ILC2 population is positive for CD127 and CD161.

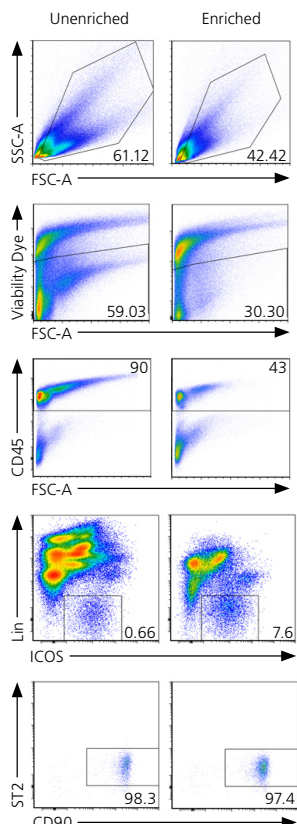
**Table 1. Antibodies for Flow Cytometric Analysis and Cell Sorting of Human ILC2s**

Labeling Antibody	Clone	Fluorochrome	Catalog #
Anti-Human CD45	HI30	PE	60018
Anti-Human CD127 (IL-7Rα)	A019D5	PE-Cy7	N/A
Anti-Human CD161 (KLRB1)	HP-3G10	PerCP-Cy5.5	N/A
Anti-Human CD294 (CRTH2)	BM16	APC	N/A
For Lineage-Specific Antigen Labeling			
Anti-Human CD1a	HI149	FITC	N/A
Anti-Human CD3	UCHT1	FITC	N/A
Anti-Human CD4	RPA-T4	FITC	N/A
Anti-Human CD11c	3.9	FITC	N/A
Anti-Human CD14	M5E2	FITC	60004
Anti-Human CD16	3G8	FITC	60041
Anti-Human CD19	HIB19	FITC	60005
Anti-Human CD34	581	FITC	60013
Anti-Human CD94	DX22	FITC	N/A
Anti-Human CD123	6H6	FITC	60110
Anti-human CD303H	201A	FITC	N/A
Anti-Human FcεR1α	AER-37	FITC	N/A
Anti-Human TCR-α/β	IP26	FITC	N/A
Anti-Human TCR-γ/δ	B1	FITC	N/A

## Identification of Mouse ILC2s from Lung

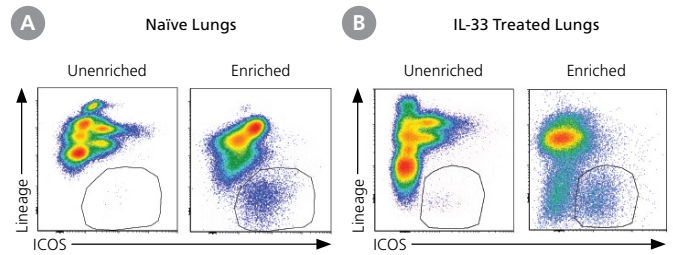
A representative gating strategy to identify mouse ILC2s from unenriched and enriched single-cell suspension from lungs is shown in Figure 3. The same gating strategy may be applied for cells from other tissues. However, the starting frequency and cell types will vary between different tissues. Therefore, ILC2s may have a somewhat different profile and the plots may look slightly different from what is shown in Figure 3. Start on the SSC-A versus FSC-A plots and create a gate to include all leukocytes. Alternatively, create a smaller gate around the lymphocytes only (not shown). When sorting cells, remove doublets and gate on single cells (singlets) using either the FSC-W versus FSC-H or SSC-W versus SSC-H plots. Then, select for viable cells and CD45<sup>+</sup> cells. Next, create a gate on the Lin<sup>-</sup> ICOS<sup>+</sup> population. Within this subset the ILC2 population is positive for CD90 and ST2. Mouse ILC2s from the lungs are described as Lin<sup>-</sup>, CD45<sup>+</sup>, ICOS<sup>+</sup>, ST2<sup>+</sup>, CD90<sup>+</sup>. ILC2s are also positive for transcription factor, GATA-3. Other markers commonly used to identify mouse ILC2s include KLRG1, CD127 and c-Kit.<sup>3,4</sup>

Intranasal administration of IL-33 in mice has been shown to increase the frequency of ILC2s found in the lung.<sup>12</sup> To get familiar with ILC2 enrichment and sorting, IL-33 treated lungs can be used as positive control (Figure 4).



**Figure 3. Gating Strategy for ILC2s From Mouse Lungs.**

Single cell suspension from C57Bl/6J mouse lung tissue were enriched using EasySep™ Mouse ILC2 Enrichment Kit (Catalog #19842) as shown in page 4, or left unenriched. The ILC2 content (Lin<sup>-</sup>CD45<sup>+</sup> ICOS<sup>+</sup> ST2<sup>+</sup>) in the enriched fraction typically ranges from 2.2 - 7.1%. In this example, the percentage of ILC2s in the unenriched and enriched fractions correspond to 0.8% and 16.65% of CD45<sup>+</sup> cells, respectively. FACS plots were generated using FCS Express 5 Software (De Novo Software).



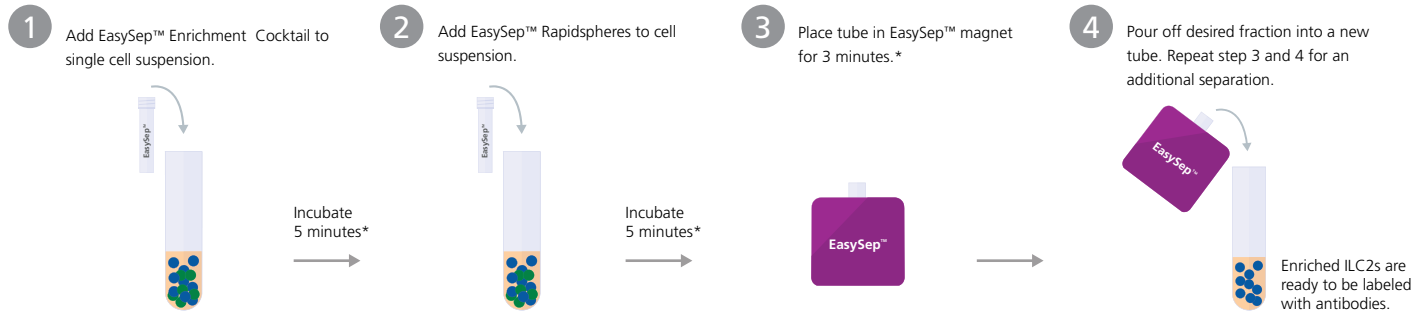
**Figure 4. ILC2s From Lungs From Naive and IL-33 Treated Mice.**

ILC2s were isolated from the lungs of naive or IL-33 treated C57Bl/6J mice as previously described.<sup>13</sup> Multiple lungs were pooled (unenriched sample) and a portion of the pooled sample was then enriched using EasySep™ Mouse ILC2 Enrichment Kit (Catalog #19842).

**Table 2. Antibodies for Flow Cytometric Analysis and Cell Sorting of Mouse ILC2s**

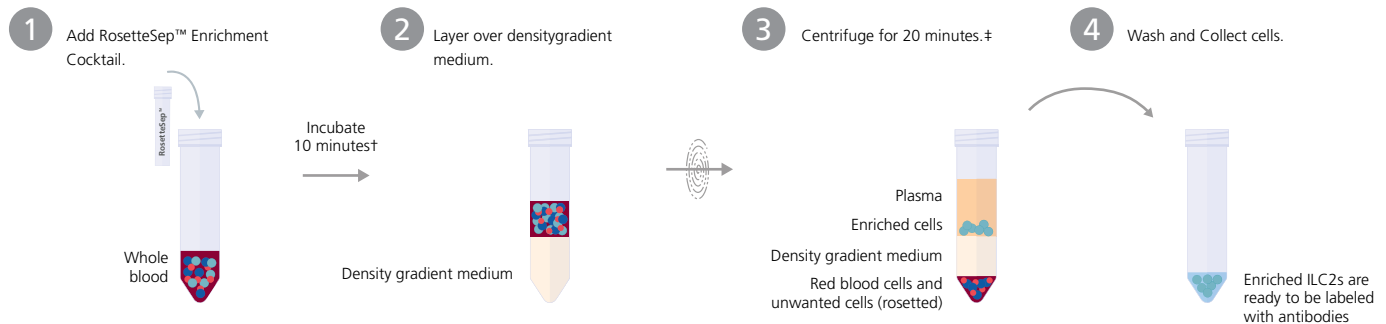
Labeling Antibody	Clone	Fluorochrome	Catalog #
Anti-Mouse CD45	30-F11	Brilliant Violet 421	60030
Anti-Mouse CD90.2 (Thy-1.2)	53-2.1	APC	60115
Anti-Mouse CD278 (ICOS)	C398.4A	PE	N/A
Anti-Mouse ST2	RMST2-2	PerCP-eFluor® 710	N/A
For Lineage-Specific Antigen Labeling			
Anti-Mouse CD3e	145-2C11	FITC	N/A
Anti-Mouse CD4	H129.19	FITC	N/A
Anti-Mouse CD11b	M1/70	FITC	60001
Anti-Mouse CD11c	N418	FITC	60002
Anti-Mouse CD19	1D3	FITC	60112
Anti-Mouse Ly-6G	RB6-8C5	FITC	60028
Anti-Mouse NK1.1 (CD161)	PK136	FITC	60103
Anti-Mouse TER119	TER-119	FITC	60033
Anti-Mouse TCR-β	H57-597	FITC	N/A
Anti-Mouse TCR-γ/δ	GL3	FITC	60104

## EasySep™ Protocol



\*Times shown are typical for EasySep™ Mouse ILC2 Enrichment Kit with the Purple EasySep™ Magnet. Times will vary depending on the specific kit and the magnet used.

## RosetteSep™ Protocol



†Times shown are typical for RosetteSep™ Human ILC2 Enrichment Kit.

‡Use SepMate™ to reduce centrifugation time to 10 minutes with brake on.

## Product Listing

Product Name	Source	Catalog #
RosetteSep™ Human ILC2 Enrichment Kit	Whole Blood	15382
EasySep™ Human ILC2 Enrichment Kit	Leukapheresis Sample (Leukopak)	17972
EasySep™ Mouse ILC2 Enrichment Kit	Mouse Single-Cell Suspension	19842

### Advantages of ILC2 Enrichment Before Sorting

**EASY.** Enrich cells with EasySep™ or RosetteSep™ without the need for columns.

**FAST.** Reduce the time spent sorting cells.

**EFFICIENT.** Improve the purity and yield of sorted ILC2s.

## References

- Artis D & Spits H. (2015) The biology of innate lymphoid cells. *Nature* 517(7534): 293–301.
- Spits H & Cupedo T. (2012) Innate lymphoid cells: emerging insights in development, lineage relationships, and function. *Annu Rev Immunol* 30(1): 647–75.
- Moro K et al. (2010) Innate production of T(H)2 cytokines by adipose tissue-associated c-Kit(+)/Sca-1(+) lymphoid cells. *Nature* 463(7280): 540–4.
- Halim TYF et al. (2012) Lung natural helper cells are a critical source of Th2 cell-type cytokines in protease allergen-induced airway inflammation. *Immunity* 36(3): 451–63.
- Neill DR et al. (2010) Nuocytes represent a new innate effector leukocyte that mediates type-2 immunity. *Nature* 464(7293): 1367–70.
- Buonocore S et al. (2010) Innate lymphoid cells drive interleukin-23-dependent innate intestinal pathology. *Nature* 464(7293): 1371–5.
- Price AE et al. (2010) Systemically dispersed innate IL-13-expressing cells in type 2 immunity. *Proc Natl Acad Sci U S A* 107(25): 11489–94.
- Spits H et al. (2013) Innate lymphoid cells--a proposal for uniform nomenclature. *Nat Rev Immunol* 13(2): 145–9.
- Walker JA et al. (2013) Innate lymphoid cells--how did we miss them? *Nat Rev Immunol* 13(2): 75–87.
- Sonnenberg GF & Artis D. (2012) Innate lymphoid cell interactions with microbiota: implications for intestinal health and disease. *Immunity* 37(4): 601–10.
- Mjösberg JM et al. (2011) Human IL-25- and IL-33-responsive type 2 innate lymphoid cells are defined by expression of CRTH2 and CD161. *Nat Immunol* 12(11): 1055–62.
- Halim TYF & Takei F. (2014) Isolation and characterization of mouse innate lymphoid cells. *Curr Protoc Immunol* 106(5): 3.25.1-13.
- Martinez-Gonzalez I et al. (2016) Allergen-Experienced Group 2 Innate Lymphoid Cells Acquire Memory-like Properties and Enhance Allergic Lung Inflammation. *Immunity* 45(1): 198–208.

Copyright © 2020 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, Scientists Helping Scientists, RosetteSep and SepMate are trademarks of STEMCELL Technologies Canada Inc. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES UNLESS OTHERWISE STATED. FOR ADDITIONAL INFORMATION ON QUALITY AT STEMCELL, REFER TO WWW.STEMCELL.COM/COMPLIANCE.