

# A single-step enrichment method for generating highly purified mouse NK cell populations by negative selection

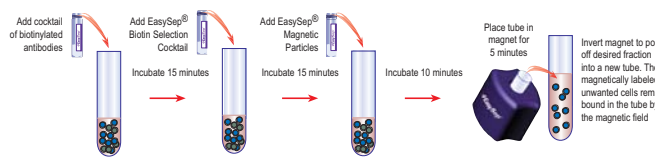
Jimmy Tran<sup>1</sup>, Sam K.P. Kung<sup>1</sup>, Ning Yuan<sup>2</sup>, Graeme Milton<sup>2</sup>, Albertus Wognum<sup>2</sup>, Allen Eaves<sup>3</sup>, Terry Thomas<sup>2</sup> and Benoit Guibault<sup>2</sup>.

<sup>1</sup>Department of Immunology, Faculty of Medicine, University of Manitoba, Winnipeg, Canada. <sup>2</sup>STEMCELL Technologies Inc., 570 W. 7th Avenue, Vancouver, BC, V5Z 1B3, Canada. <sup>3</sup>Terry Fox Laboratory 675 W. 10th Avenue, Vancouver, BC V5Z 1L3, Canada

## Summary

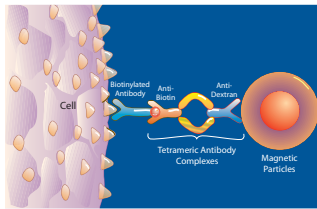
Natural Killer (NK) cells play a critical role in innate immunity and tumor immunology. They are particularly difficult to isolate because they are rare and express several markers found on other cell types. Using column-free immunomagnetic cell separation technology (EasySep<sup>®</sup>), we sought to develop a single-step negative selection method for NK cell isolation, where negative selection enables the isolation of cells with minimal impact on cell function. Unwanted cells were cross-linked to magnetic particles using a highly optimized cocktail of 10 biotinylated antibodies. This enabled us to achieve NK cell purities of  $92.3 \pm 2.6\%$  ( $n=6$ ; defined as %CD49b<sup>+</sup>CD3<sup>neg</sup>). The isolated cells retained the expression of several markers associated with an NK phenotype, including NKG2D and LY49C/I/F/H. The purified cells could be readily expanded when cultured in the presence of IL-2 (> 8-fold expansion after 5 days). <sup>51</sup>Cr release assays demonstrated the cytotoxic capacity of the isolated cells, and NK-specific killing was confirmed using both NK-sensitive and NK-resistant cell lines. In conclusion, we present here a novel and unique enrichment method for the isolation of pure NK cells in a single step.

## Methods



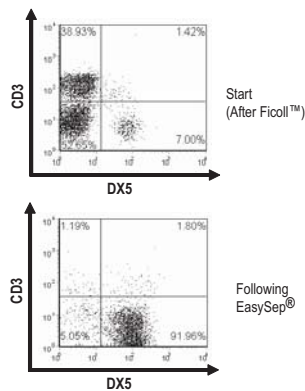
**Figure 1. EasySep<sup>®</sup> procedure for negative selection of mouse cells**

Mouse cells are targeted for immunomagnetic depletion by first labeling the cells with a cocktail of anti-cell biotinylated antibodies, while cells of interest are not targeted for depletion. Labeled (unwanted) cells are then crosslinked to EasySep<sup>®</sup> magnetic particles. The tube containing the labeled cell suspension is placed directly in the specially designed EasySep<sup>®</sup> Magnet. This light, handheld magnet is gently inverted to remove the cells that are not bound to the magnetic particles and the desired cells are poured off into a new tube.



**Figure 2. Schematic drawing of EasySep<sup>®</sup> magnetic labelling of mouse cells**

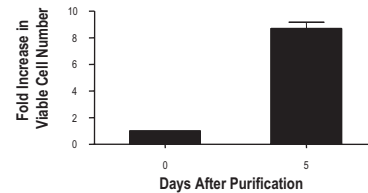
Tetrameric Antibody Complexes are comprised of two mouse IgG<sub>1</sub> monoclonal antibodies held in tetrameric array by two rat anti-mouse IgG<sub>1</sub> antibody molecules. One mouse antibody recognizes biotin while the other recognizes dextran on the EasySep<sup>®</sup> magnetic particles.



**Figure 3. Isolation of NK cells using the EasySep<sup>®</sup> Mouse NK Cell Enrichment Kit**

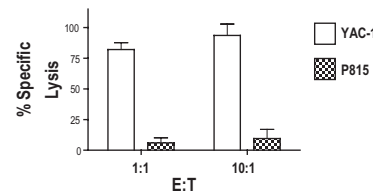
NK cells were isolated from splenocytes of either C57BL/6 or BALB/c mice. Red blood cells were initially separated from splenocytes by Ficoll-Paque™. NK cells were then purified using the EasySep<sup>®</sup> Mouse NK Cell Enrichment Kit (STEMCELL Technologies). NK purity was assessed by surface staining with DX5-PE and CD3-PECy5 and detection using flow cytometry.

## Results



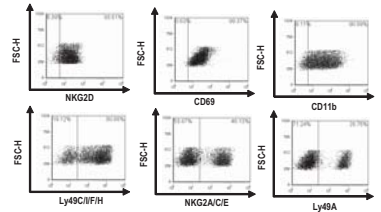
**Figure 4. Isolated NK cells proliferate and expand in number when cultured in IL-2 supplemented medium**

Isolated NK cells were cultured in complete media (RPMI 1640 supplemented with 10% fetal bovine serum containing penicillin, streptomycin, glutamine and β-mercaptoethanol) supplemented with 1000 IU/mL IL-2 for 5 days. NK cell number was determined by viability counting with Trypan Blue. Values are expressed as a ratio of the number of cells obtained after culture relative to the number plated cells on day 0, and are averages of 3 experiments with error bars indicating standard deviation from the mean.



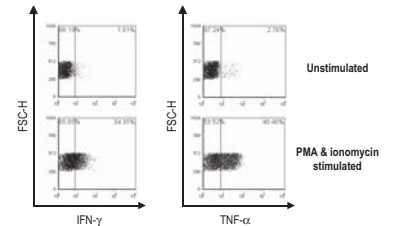
**Figure 5. Culture-expanded NK cells display NK cell-mediated cytotoxicity and target specificity**

Isolated NK cells were cultured in complete media supplemented with 1000 IU/mL IL-2 for 5 days. Activated NK cells were then used as effectors in a 4-hour <sup>51</sup>Cr release assay using either the NK-sensitive YAC-1 line or NK-resistant P815 line.



**Figure 6. Isolated NK cells retain their polyclonal nature following expansion**

Isolated NK cells were cultured in media supplemented with 1000 IU/mL IL-2 for 5 days. NK cell surface receptor expression profiling was performed by staining with fluorochrome conjugated anti-NKG2D, CD69, CD11b, Ly49C/I/F/H, NKG2A/C/E, and Ly49A, and detection using flow cytometry analysis. Similar results were obtained in three separate experiments.



**Figure 7. Expanded NK cells produce IFN-γ and TNF-α when stimulated with PMA and ionomycin**

Purified NK cells were cultured in media supplemented with 1000 IU/mL IL-2 for 5 days. Activated NK cells were stimulated in ionomycin and phorbol 12-myristate 13-acetate (PMA) overnight to induce IFN-γ and TNF-α production. The cells were intracellularly stained with IFN-γ-AlexaFluor and TNF-α-PE and analyzed by flow cytometry.

## Conclusions

- Mouse NK cells can be highly purified in a single step using the EasySep<sup>®</sup> Mouse NK Cell Enrichment Kit
- The purified cells can be expanded in culture while retaining the phenotypic and functional properties of mouse NK cells.

Supported by the Manitoba Health Research Council and The Dean of Medicine Strategic Research Fund. Ficoll™ is a trademark of GE Healthcare Ltd.



WWW.STEMCELL.COM

IN NORTH AMERICA TOLL-FREE 1. 800.667.0322 TOLL-FREE 1. 800.567.3899 F. 1. 604.877.0713 E. 1. 604.877.0704 I.INFO@STEMCELL.COM  
IN EUROPE T. +33 (0)4 76 04 75 30 F. +33 (0)4 76 18 99 63 E.INFO@STEMCELLFRANCE.COM  
IN AUSTRALIA TOLL-FREE 1. 800.060.350 T. 07.5474.5042 E.INFO.AUS@STEMCELL.COM