

**Description**

Expanded Human Peripheral Blood NK Cells are NK cells enriched and expanded from human PBMCs using the NK Expansion Kit (BPS Bioscience 78927), a K562 feeder cell-based system for about 2 weeks, and cryopreserved. Human Peripheral Blood NK Cells are >90 % pure NK cells (CD3<sup>-</sup>CD56<sup>+</sup> cells), as measured by flow cytometry analysis. They can be used in NK cytotoxicity assays or ADCC (antibody-dependent cellular cytotoxicity), after thaw and recovery in NK Cell Basal Medium for 24 hours or can be further expanded using the BPS Bioscience NK Expansion Kit.

**Background**

NK (natural killer) cells are part of the innate immune system. They function in a histocompatibility complex-independent mode and derive from the hematopoietic lineage. They are the first line of defense against cancer. Expression of marker CD56 correlates with NK cell functionality: the CD56<sup>bright</sup> subset accounts for about 5% of the population and is less cytotoxic than the CD56<sup>dim</sup> subset. Cytotoxicity can happen by the release of perforin and granzyme, while activation by KARs (killer activating receptors) leads to release of Fas Ligand, TRAIL (TNF-related apoptosis-inducing ligand) and TNF $\alpha$  (tumor necrosis factor-alpha). In a suppressive tumor microenvironment, NK cells can become inhibited and unable to fight cancer cells. Several clinical trials have focused on using *ex vivo* generated NK cells alone or in combination with other approaches. NK cells can be generated *ex vivo* from peripheral blood, umbilical cord blood, iPS cells or immortalized NK cell lines. The ability to generate a number of pure cells high enough for human dosage often requires the use of growth factors such as IL-2 (interleukin 2) or IL-15, and feeder cells. The use of NK cells or CAR (chimeric antigen receptor)-NK cells is an expanding area holding great promise in cancer therapy.

**Application**

- Use in NK cell cytotoxicity assays.
- Use in NK cell mediated ADCC assays.

**Materials Provided**

Components	Format
1 vial of frozen cells	Vial contains 5 x 10 <sup>6</sup> cells in 1 ml of CryoStor <sup>®</sup> CS10 (Stemcell Technologies #100-1061)

**Mycoplasma Testing**

The cells have been screened to confirm the absence of Mycoplasma species.

**Storage Conditions**

NK Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a -80°C freezer for long term storage. Contact technical support at [support@bpsbioscience.com](mailto:support@bpsbioscience.com) if the cells are not frozen in dry ice upon arrival.

**Materials Required but Not Supplied**

These materials are not supplied with NK cells but necessary for NK characterization and cytotoxicity assays. BPS Bioscience's reagents are validated and optimized for use and are highly recommended for the best results.

Name	Ordering Information
NK Cell Basal Medium	BPS Bioscience #82164
Growth-Arrested NK Feeder Cells	<a href="#">BPS Bioscience #78712</a>
Anti-NCAM1 Antibody, PE-Labeled	<a href="#">BPS Bioscience #101673</a>
Anti-CD3 Antibody, FITC-Labeled	<a href="#">BPS Bioscience #102008</a>
Human Interleukin-2	<a href="#">BPS Bioscience #90184</a>
500 x CFSE	BPS Bioscience #82177
1000 x 7-AAD	BPS Bioscience #82178
K562	ATCC #CCL-243
eGFP/Firefly Luciferase K562 Cell Line	BPS Bioscience #78911
eGFP/Firefly Luciferase RS4; 11 Cell Line	BPS Bioscience #78926
Clear-bottom, white 96-well tissue culture-treated plate	Corning #3610
U-bottomed 96-well plate	Corning #3799
ONE-Step™ Luciferase Assay System	<a href="#">BPS Bioscience #60690</a>
Thaw Medium 2	BPS Bioscience #60184
Luminometer	

**Recommended NK Medium:** NK Cell Basal Medium (BPS Bioscience #82164) supplemented with Interleukin-2 (BPS Bioscience #90184). NK cells can be further expanded using BPS Bioscience feeder cell-based NK Expansion Kit (BPS Bioscience #78927).

#### Cell Thawing and Culture Protocol:

- The expansion fold obtained will vary, depending on the source of NK cells and donor.
  - It is recommended that cell cultures are monitored daily, and they are split and fed in order to keep optimal cell density (< 2 million/ml).
  - Flow cytometry analysis for typical NK markers, such as CD3 and CD56, can be performed to monitor NK purity and determine NK fold expansion.
  - Expanded Human Peripheral Blood NK Cells can be expanded further for 3~5 weeks, or other desired expansion period, and then used in downstream applications. Generally, Expanded Human Peripheral Blood NK Cells can reach >90% purity (CD3<sup>-</sup>CD56<sup>+</sup> cells) and reach >20-fold expansion after further expansion for 1 week.
1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. As soon as the cells are thawed (it may be slightly faster or slower than 60 seconds), quickly transfer the entire contents of the vial to a tube containing 10 ml of pre-warmed NK Cell Basal Medium.  
**Note: Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.**
  2. Immediately spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in 20 ml of pre-warmed NK Cell Basal Medium **supplemented with 20 ng/ml Interleukin-2.**
  3. Transfer cells into one T75 flask and grow in a 5% CO<sub>2</sub> incubator at 37°C.

4. After 24 hours in a 5% CO<sub>2</sub> incubator at 37°C, NK cells can be directly used in NK cytotoxicity assays or ADCC assays.

*Note: For further expansion follow instructions described below.*

5. Add Growth-Arrested NK Feeder Cells to NK cells at a cell-to-cell ratio of 1:1, and grow the cells in a 5% CO<sub>2</sub> incubator at 37°C.
6. Determine cell density every 2-3 days. When the cell density reaches >2 million/ml, dilute cell to 0.25-0.5 x 10<sup>6</sup> cells/mL with NK Cell Basal Medium supplemented with 20 ng/ml IL-2.
7. Incubate the cells at 37°C and 5% CO<sub>2</sub> in a humidified incubator.
8. Refresh medium of NK cells every 2-3 days and refresh feeder cells by providing NK cells with a 1:1 ratio of feeder cells:NK cells weekly.

#### Validation Data

- The following experiments are an example of co-culture assay used to evaluate the cytotoxicity of NK cells using eGFP-Firefly Luciferase K562 as target cells.
- K562, a human erythromyeloblastoid leukemia cell line, is a NK target due to the lack of HLA expression on the cell surface. RS4;11, a lymphoblast cell line that expresses HLA-C alleles, that bind the most expressed KIRs (killer-cell immunoglobulin-like receptors), are NK resistant. RS4;11 cells were used as negative control in the NK cytotoxicity assays.

#### Luciferase activity-based NK cytotoxicity assay using eGFP-Luciferase K562 Cell Line as target cells.

- The assay should include a “Minimum Viability Control” (or MIN), “Maximum Viability Control” (or MAX) and “Test” conditions.
  - Samples and controls should be run in triplicate.
1. Harvest eGFP-Luciferase K562 cells and seed at 5,000 cells/well in 50 µl of Thaw Medium 2 in a 96-well white, clear bottom tissue culture plate.
  2. Prepare Thaw Medium 2 with 2% SDS (50 µl/well of “Minimum Viability Control” wells).
  3. Resuspend expanded NK cells in Thaw Medium 2 at the appropriate concentrations to reach the desired Effector:Target (E:T) ratios (50 µl/well).
  4. Add 50 µl of NK cells to the target cells by carefully pipetting NK cells to the “Test” wells containing eGFP-Luciferase K562 cells. The total volume of each well is now 100 µl.
  5. Add 50 µl of Thaw Medium 2 to the “Maximum Viability Control” wells.
  6. Add 50 µl of Thaw Medium 2 with 2% SDS to the “Minimum Viability Control” wells.
  7. Incubate the plate at 37°C with 5% CO<sub>2</sub> for 4 hours.

8. Add 100 µl of ONE-Step™ Luciferase reagent.
9. Incubate for 15-30 minutes.
10. Measure luminescence signal in a microplate reader capable of reading luminescence.

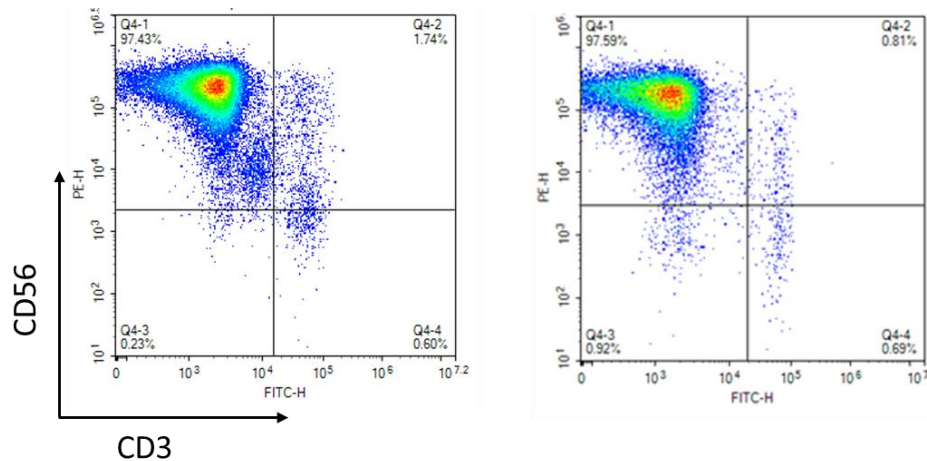
**Data Analysis:** For each target, 3 replicates of the internal references for the 0% viability background (MIN) and the 100% viability maximal signal (MAX) were run.

Percent viability = [(mean luminescence of the experimental sample - mean luminescence of MIN)/(mean luminescence of the MAX - mean luminescence of MIN)] x 100.

Percent Specific Lysis is calculated as follows:

$$\% \text{ specific lysis} = [1 - (\text{experimental value} - \text{MIN value}) / (\text{MAX value} - \text{MIN value})] \times 100.$$

### Data and Figures



*Figure 1. Flow cytometry analysis of Expanded Human Peripheral Blood NK Cells.*

Expanded Human Peripheral Blood NK cells were thawed and expanded *ex vivo* for one week (left) and two weeks (right), respectively. Cells were collected and stained with Anti-CD3 Antibody FITC-Labeled and Anti-NCAM-1 Antibody, PE-Labeled and analyzed by flow cytometry.

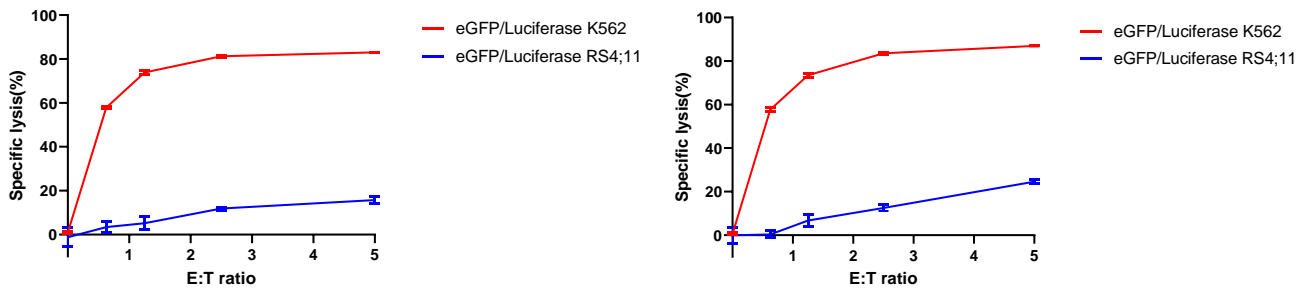


Figure 2. Luciferase-Based cytotoxicity profile of Expanded Human Peripheral Blood NK Cells.

Expanded Human Peripheral Blood NK cells were co-cultured with eGFP/Luciferase K562 target cells and eGFP/Luciferase RS4;11 control cells at indicated ratios of Effector: Target (E:T) cells at 37°C for 4 hours. ONE-Step™ Luciferase Assay was used to detect NK cell cytotoxicity on the luciferase-expressing target cell. Killing curves were obtained based on different Effector: Target (E:T) ratios.

Data shown is representative. For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

## References

Du N., et al., 2021 *Cancers (Basel)* 13 (16): 4129.

## License Disclosure

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

## Troubleshooting Guide

Visit [bpsbioscience.com/cell-line-faq](https://bpsbioscience.com/cell-line-faq) for detailed troubleshooting instructions. For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

## Related Products

Products	Catalog #	Size
NCAM1/CD56 Positive Cell Isolation Kit	78808	1 x 10 <sup>8</sup> cells/1 x 10 <sup>9</sup> cells
NKp46 CHO Cell Line (High, Medium or Low Expression)	78916	2 vials
NKG2D, Avi-Tag, Fc fusion Recombinant	100252	100 µg
NKG2D, Avi-Tag, Fc fusion, Biotin-labeled Recombinant	100313	25 µg/50 µg
NKp46 Lentivirus	78717	500 x 2
Anti-NCAM1 (CD56) IgG Antibody, Biotin-labeled	101112	100 µg