

## Introduction

Major advances have been made in harnessing natural killer (NK) cells in cancer immunotherapy in recent years. Regulated by their germ-line encoded activating and inhibitory receptors, NK cells can recognize and eliminate tumor cells rapidly without prior sensitization. Clinical evidence has shown that donor-derived NK cells have low risk in inducing graft-versus-host-disease as well as a reduced risk of life-threatening cytokine storms during

therapy. Both properties make NK cells ideal for allogeneic application. To enhance the anti-tumor specificity and efficiency, NK cells can be further modified by chimeric antigen receptors (CARs). Promising clinical outcome of CAR-engineered NK cells has been recently reported for patients with relapsed or refractory B cell malignancies.

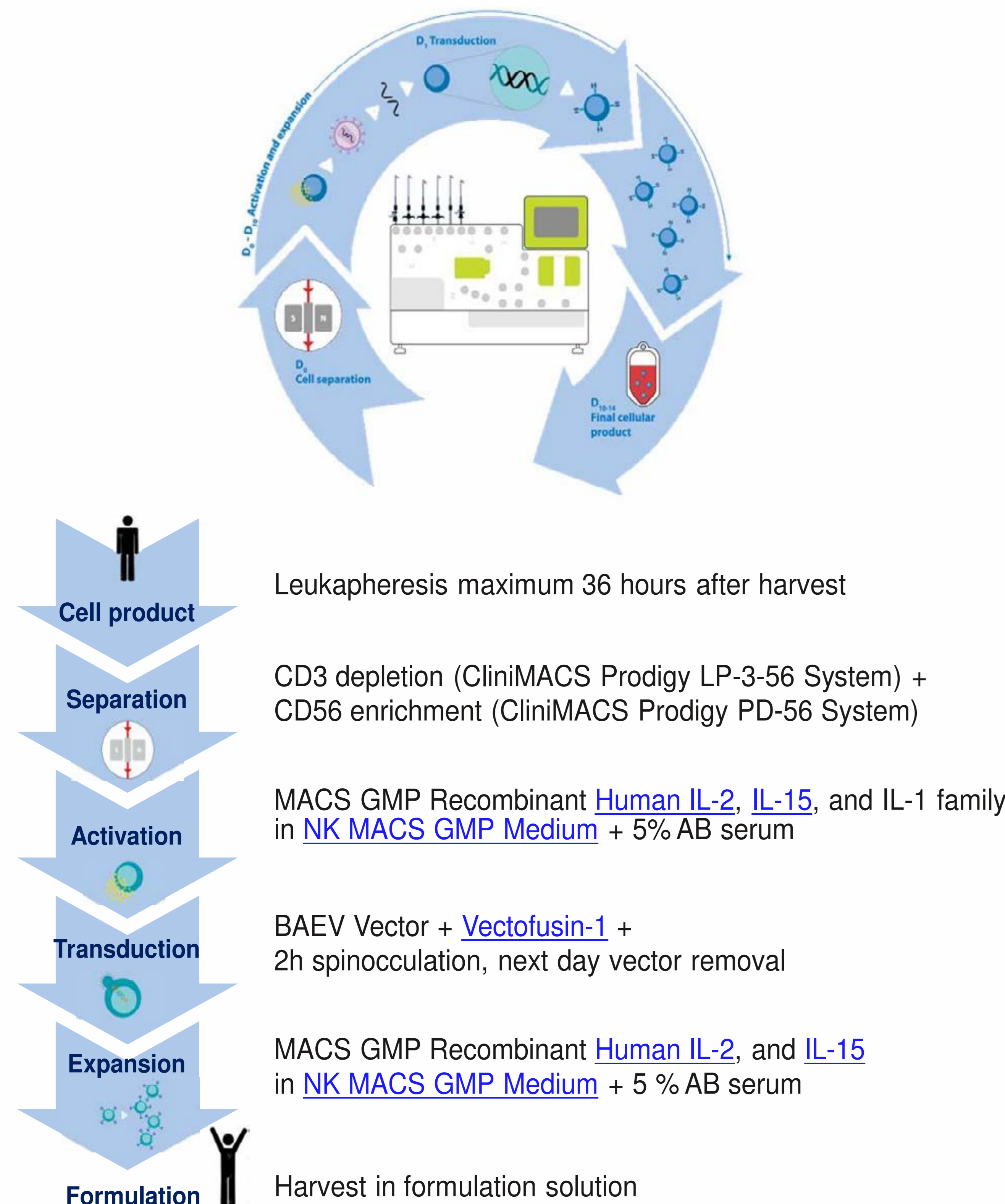
## Methods

### 1 Leukapheresis derived NK cell culture using CliniMACS Prodigy®

NK cells can be derived from leukapheresis using the CliniMACS Prodigy® System. The first step in this CliniMACS Prodigy® Natural Killer cell Engineering (NKE) System takes place in CliniMACS Prodigy® TS 320 using CliniMACS Prodigy® LP-3-56 System to deplete T cells. After successfully depleting T cells, resulting in less than 0.1 %

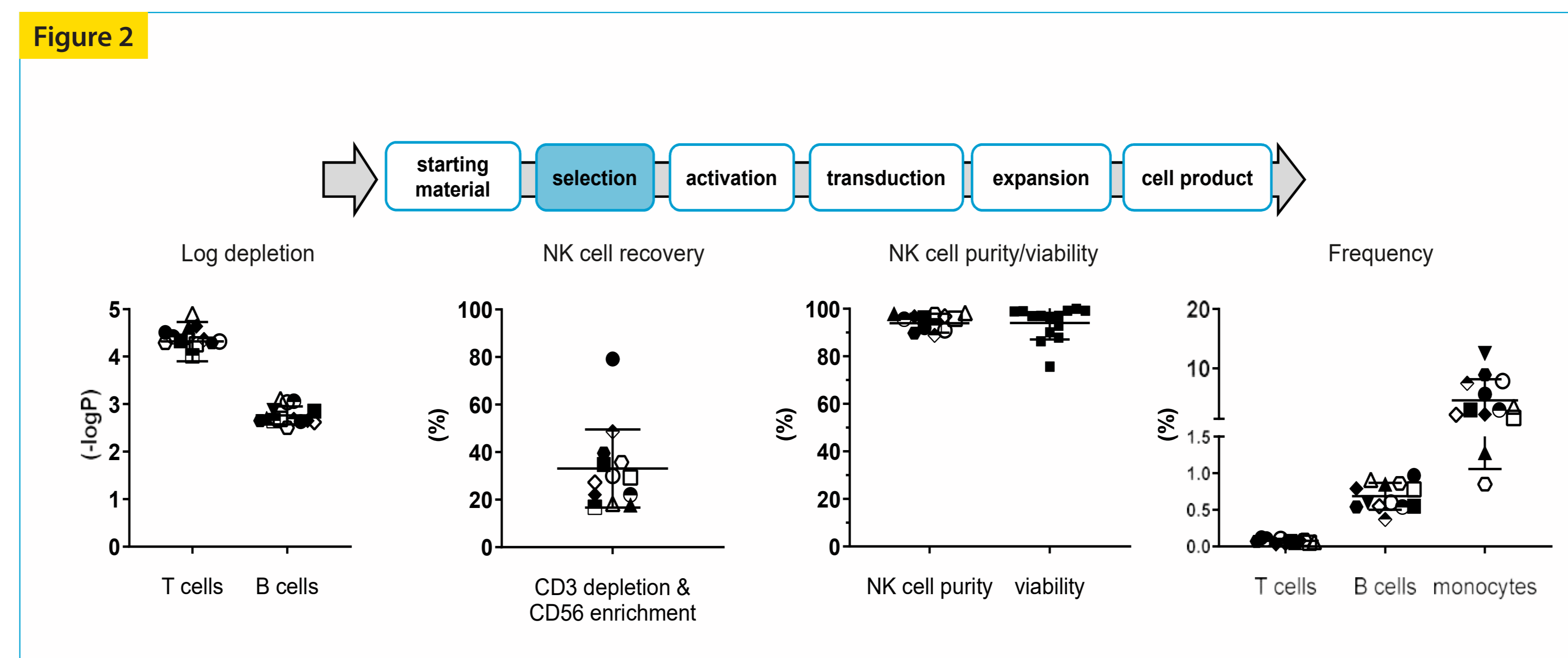
CD3<sup>+</sup> cells, the CD56<sup>+</sup> NK cells are enriched using CliniMACS Prodigy® TS 520 within the CliniMACS Prodigy® PD-56 Engineering System. Subsequent to NK cell cytokine based activation the NK cell transduction is performed at culture day 2. The CAR<sup>+</sup>/NK cells are cultivated and harvested at culture day 14 (fig. 1).

### Figure 1 CliniMACS Prodigy Natural Killer cell Engineering System



## Results

### 2 NK isolation using CliniMACS Prodigy® NKE System

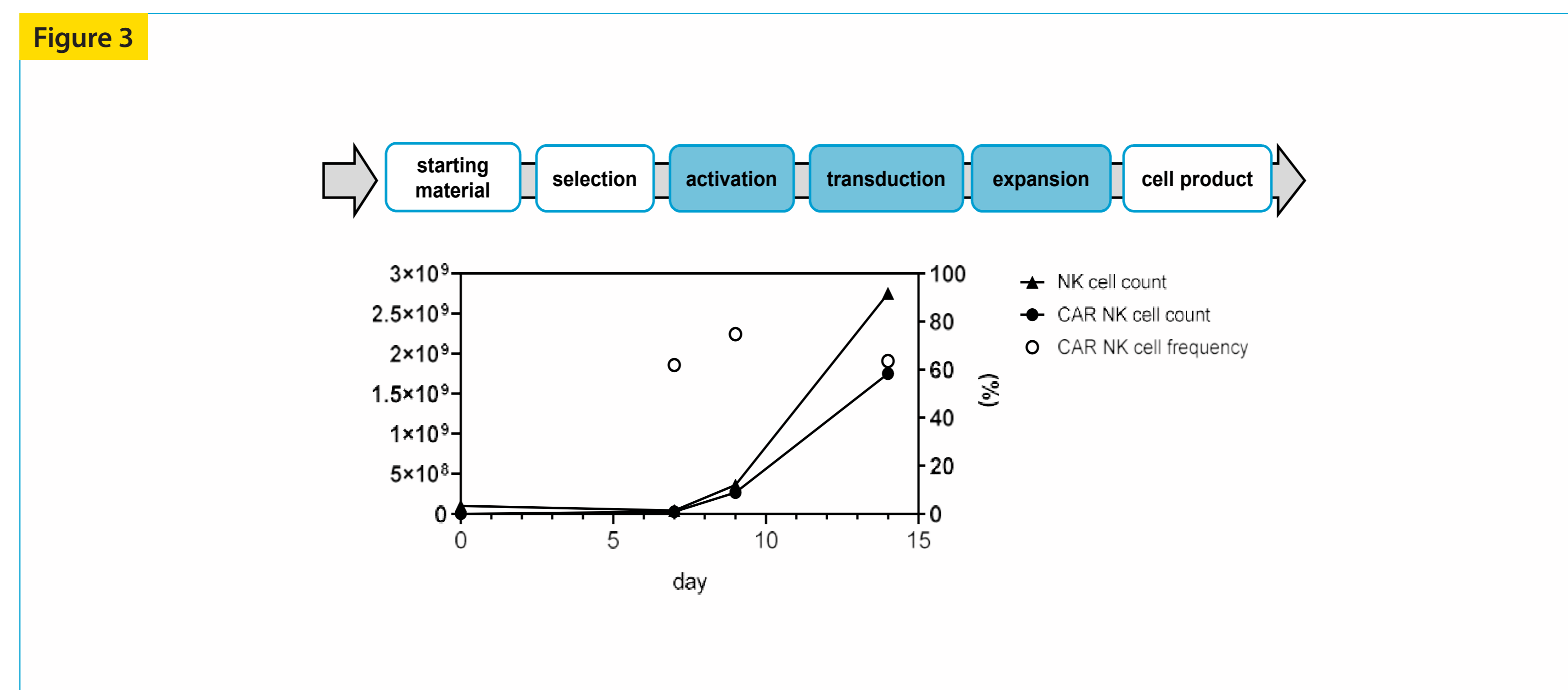


The CliniMACS Prodigy® LP-3-56 System enables a fully automated, two-step NK cell separation in a single tubing set (CliniMACS Prodigy® TS 320). The process is designed in distinct blocks, enabling the use of CD3 depletion as a stand-alone process, as it is used in the CliniMACS Prodigy® NKE System.

The CliniMACS Prodigy® PD-56 Engineering System enables a fully automated NK cell separation with subsequent culture. Starting with CD3 pre-depleted material

derived from CliniMACS Prodigy® LP-3-56 System, the CD56<sup>+</sup> enrichment, using CliniMACS Prodigy® PD-56 Engineering System, results in a highly pure (93 %) and viable (93 %) NK cell product (fig. 2 NK cell purity/viability). The recovery of NK cells in mean is about 33 % (fig. 2 NK cell recovery). An average log depletion of T cells of 4.3 and B cells of 2.7 (fig. 2 Log depletion) results in a low frequency of contaminating T and B cells as well as monocytes (fig. 2 Frequency).

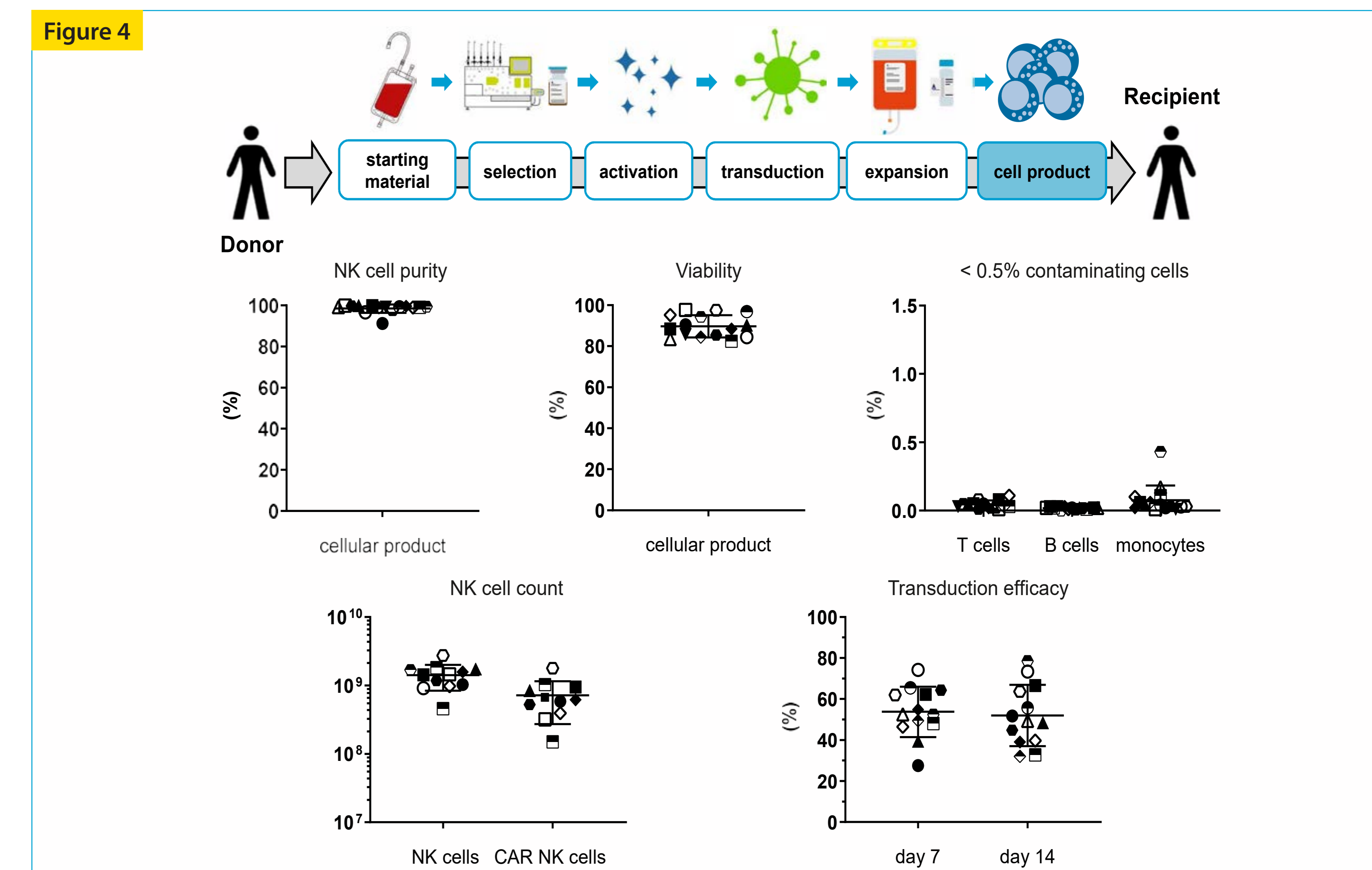
### 3 NK cell culture and CAR generation in CliniMACS Prodigy® NKE System



NK cells derived from CD3 depletion and CD56 enrichment are activated at day 0 with MACS GMP Recombinant Human IL-2, MACS GMP Recombinant Human IL-15 and a 3<sup>rd</sup> cytokine from IL-1 family. The transduction process using baboon-enveloped lentiviral vector takes place at day 2 of culture in the presence of MACS GMP Vectofusin-1<sup>®</sup> and a subsequent spinoculation for 2h. A cell wash at day 3 removes excess lentiviral vector as well

as the 3<sup>rd</sup> cytokine. The CAR<sup>+</sup>/NK cell culture is continued until day 14 using MACS GMP Recombinant Human IL-2 and IL-15 in NK MACS GMP Medium with 5 % AB serum. Figure 3 exemplifies a CAR<sup>+</sup>/NK growth curve resulting in 2.8x10<sup>9</sup> NK cells with a transduction efficiency of 64 % (1.8x10<sup>8</sup> CAR-NK cells) at day 14 without using feeder cells for expansion.

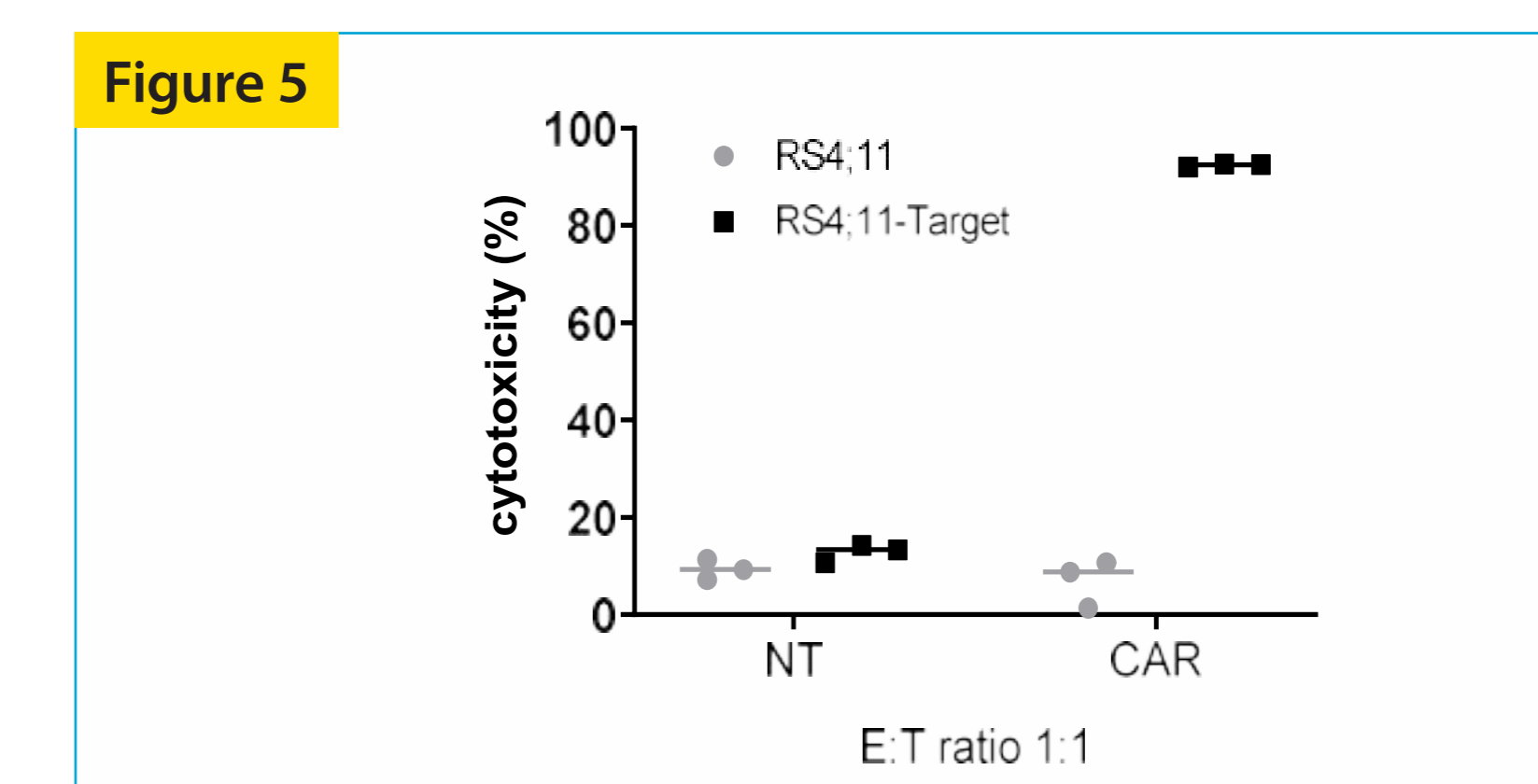
### 4 Quality assessment of final CAR-NK cell product



The transduction of NK cells in a GMP-compliant system resulted in a highly pure (>99 %) and viable (90 %) NK cell product with low CD3<sup>+</sup> cell content (mean 0.09 %). On average the expansion resulted in 1.19x10<sup>9</sup> NK cells

including 5.7x10<sup>8</sup> CAR-NK cells. The transduction efficacy during culture is high (mean 50 %) and stable compared between day 7 and 14.

### 5 Functional assessment of final CAR-NK cell product



To test the in-vitro killing capability, the generated CAR<sup>+</sup>/NK cells were challenged using a resistant tumor cell line. After 24 h and an E:T ratio of 1:1 no killing of the tumor cell line using not-transduced (NT) NK cells was observed, the tumor model therefore works well. The transduced CAR-NK cells show no cytotoxicity in the absence of the target on the tumor cells, but show CAR-mediated NK cytotoxicity in the presence of the target. Apart from the big safety aspect of the CAR-NK cells they show a high killing capability only in the presence of the target.

## Conclusion

- Development of novel process for automated NK purification, transduction and cultivation in a closed GMP compatible system.
- High level of automation enables standardized, consistent and operator-independent genetic engineering of NK cells.
- NK cell transduction using baboon-enveloped LV

- results in high and stable transduction efficacy.
- After 14 days of culture, the generated CAR-NK cells show high purity and viability as well as in-vitro functionality.

To get more information contact Martha Elia Luevano Salinas: marthal@miltenyibiotec.de