

EXPANSION SYSTEM FOR PRODUCING A LARGE AMOUNT (>10⁹ CELLS) AND HIGH PURITY (>90%) OF HUMAN CD3⁺CD56⁺NK CELLS FROM PBMCs AND THEIR THERAPEUTIC APPLICATION DUE TO ADCC ACTIVITY IN XENOGENIC MOUSE MODEL



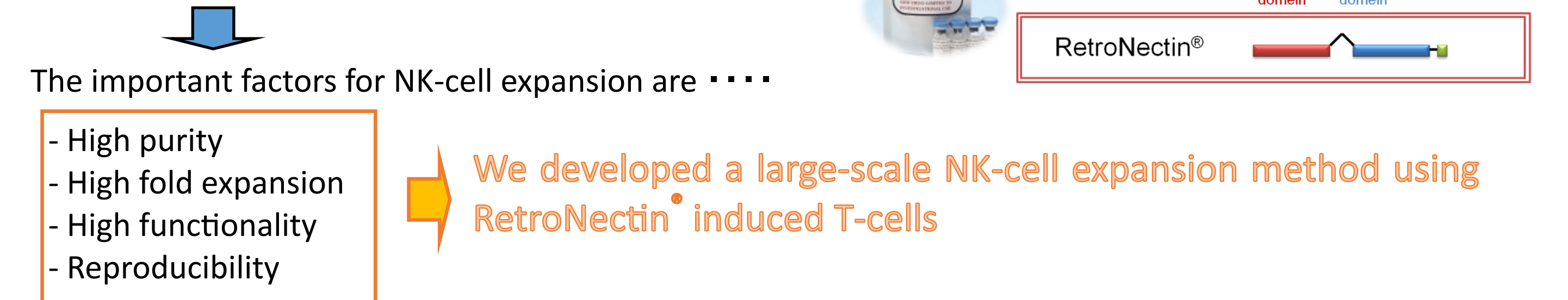
Tatsuji Enoki, Mitsuko Ideno, Akiko Kato, Junichi Mineno
CDM Center, Takara Bio Inc.

Abstract

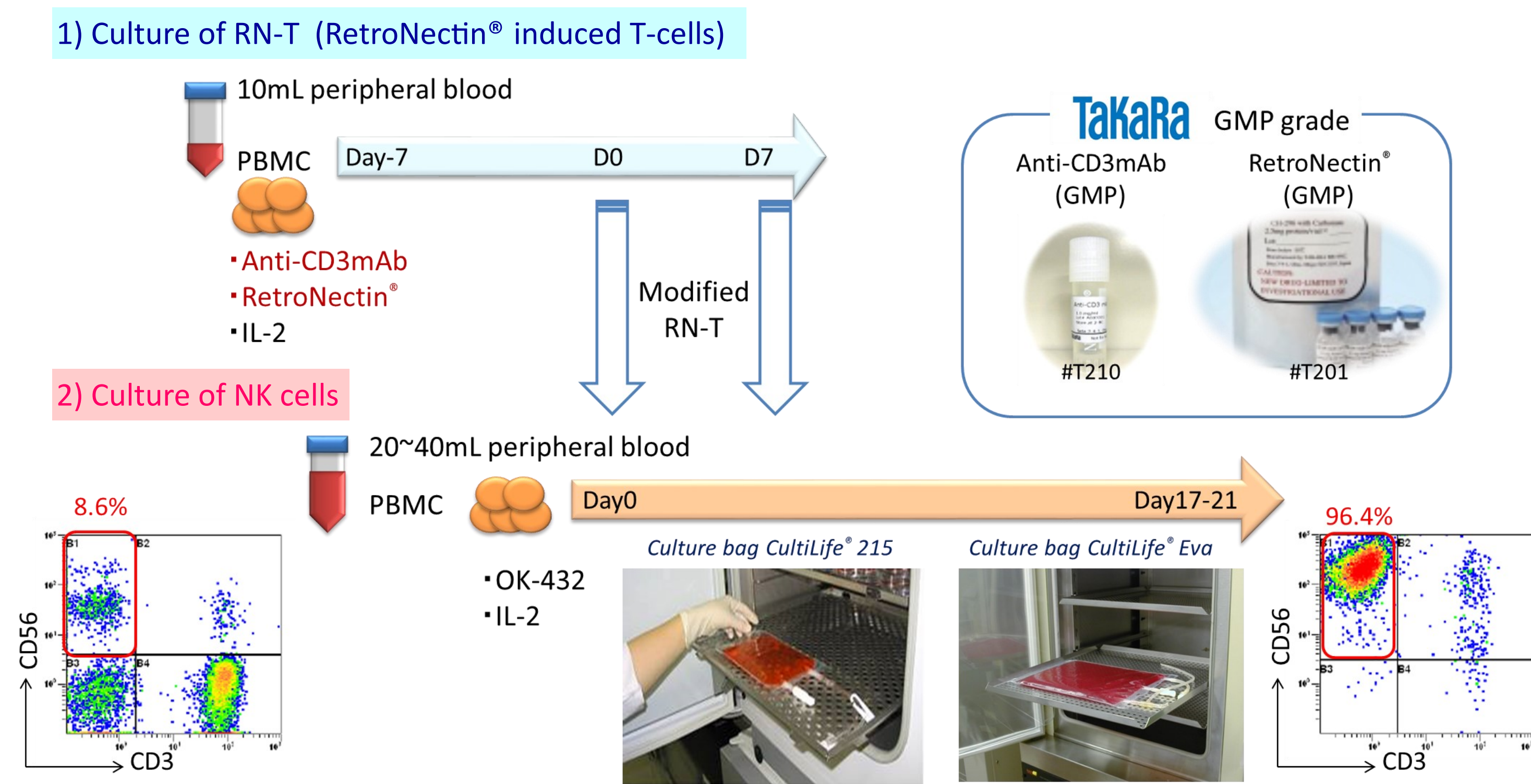
Human natural killer (NK) cells are one of attractive candidates for cell-based therapy against any cancers due to their strong cytotoxicity. However, there are few convenient and efficient method to obtain a large amount and high purity of functional NK cells from peripheral blood mononuclear cells (PBMCs) derived from a small amount of blood. Thus, we have developed a robust NK-cell expansion method using OK-432, IL-2 and RetroNectin[®] induced T (RN-T) cells as a stimulator. RN-T cells were prepared by previously established co-stimulation method using anti-CD3mAb and RetroNectin[®], and treated to suppress the growth potential (modified RN-T cells). NK cells could be expanded from PBMCs stimulated with modified RN-T cells, OK-432 and IL-2, then cultured for more than 16 days. In our large-scale culture system using gas-permeable culture bag (CultiLife[™]215 and CultiLife[™]Eva), we could obtain 10⁹ – 10¹⁰ cells containing a high proportion (>90%) of CD3-CD56+ NK cells from 50mL of peripheral blood. Furthermore, almost all cells displayed functional cell surface molecules such as NKG2D and CD16 implicated in cytotoxicity and antigen dependent cell cytotoxicity (ADCC). Thus, we investigated the antitumor effect of the expanded NK cells combined with Trastuzumab against HER2-positive human gastric cancer cell line NCI-N87 in hIL-2 Tg NOG mice (hIL-2-NOG mice; Central Institute for Experimental Animals). In this experiment, we used purified NK cells to reduce GVHD risk caused by human CD3+ cells including in the expanded cells. As a result, the combination of the NK cells and Trastuzumab dramatically enhanced the antitumor activity compared with each treatment alone. The chimerism of human NK cells in mouse peripheral blood was observed during the observation period without any GVHD symptoms and functional NK-cell surface markers such as CD16 and NKG2D also expressed in human NK cells. Furthermore, human NK cells were observed into tumor tissue even in 3 months after administration. Overall, we have established a robust NK-cell expansion system and the expanded cells showed strong antitumor activity in a xenogenic mouse model. It is considered that our expansion system could be used for chimeric antigen receptor (CAR)-NK cell processing or pluripotent stem cell derived NK-cell manufacture for future application.

The issues of NK-cell expansion for therapeutic application

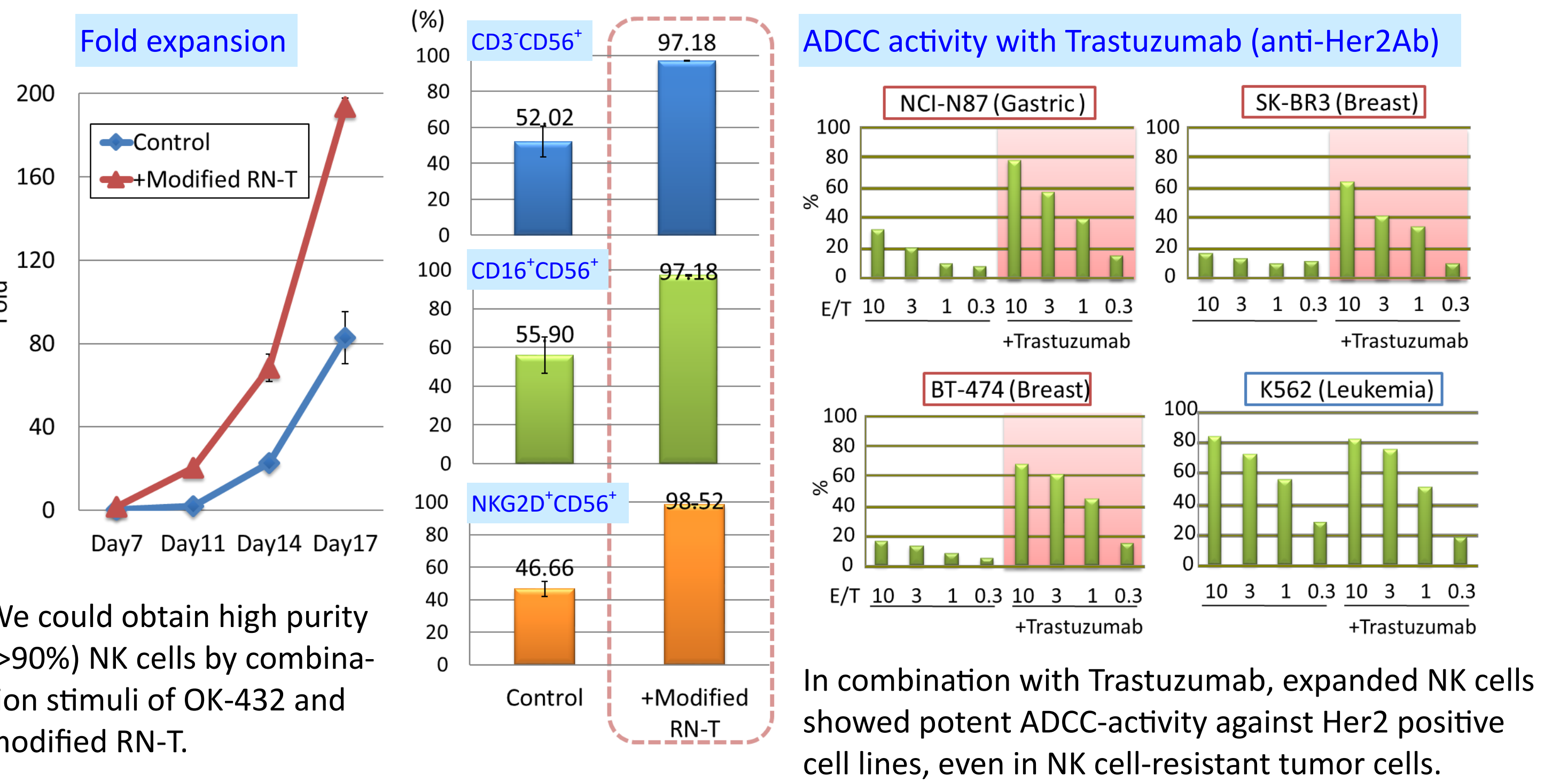
- ✓ Low purity of NK cells after expansion
- ✓ Low proliferation of NK cells



Procedure for NK-cell expansion



NK-cell expansion using modified RN-T

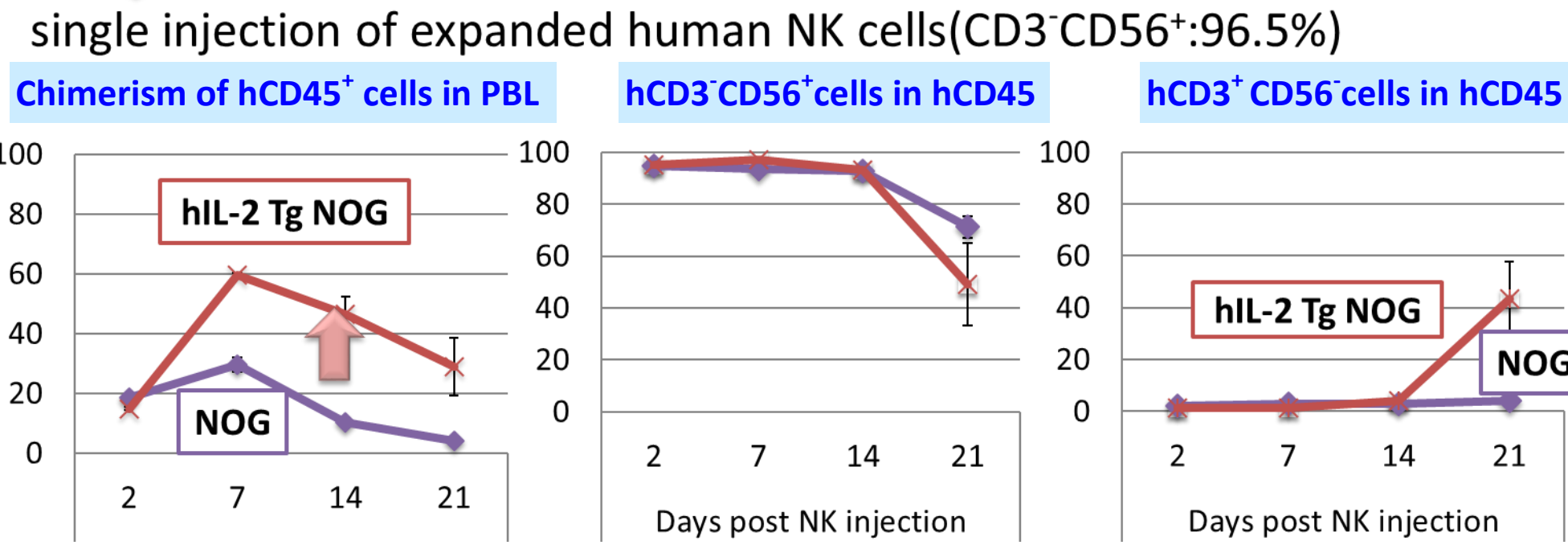


We could obtain high purity (>90%) NK cells by combination stimuli of OK-432 and modified RN-T.

hIL-2 Tg NOG mice model and study design for evaluating in vivo ADCC activity

hIL-2 Tg NOG mice model

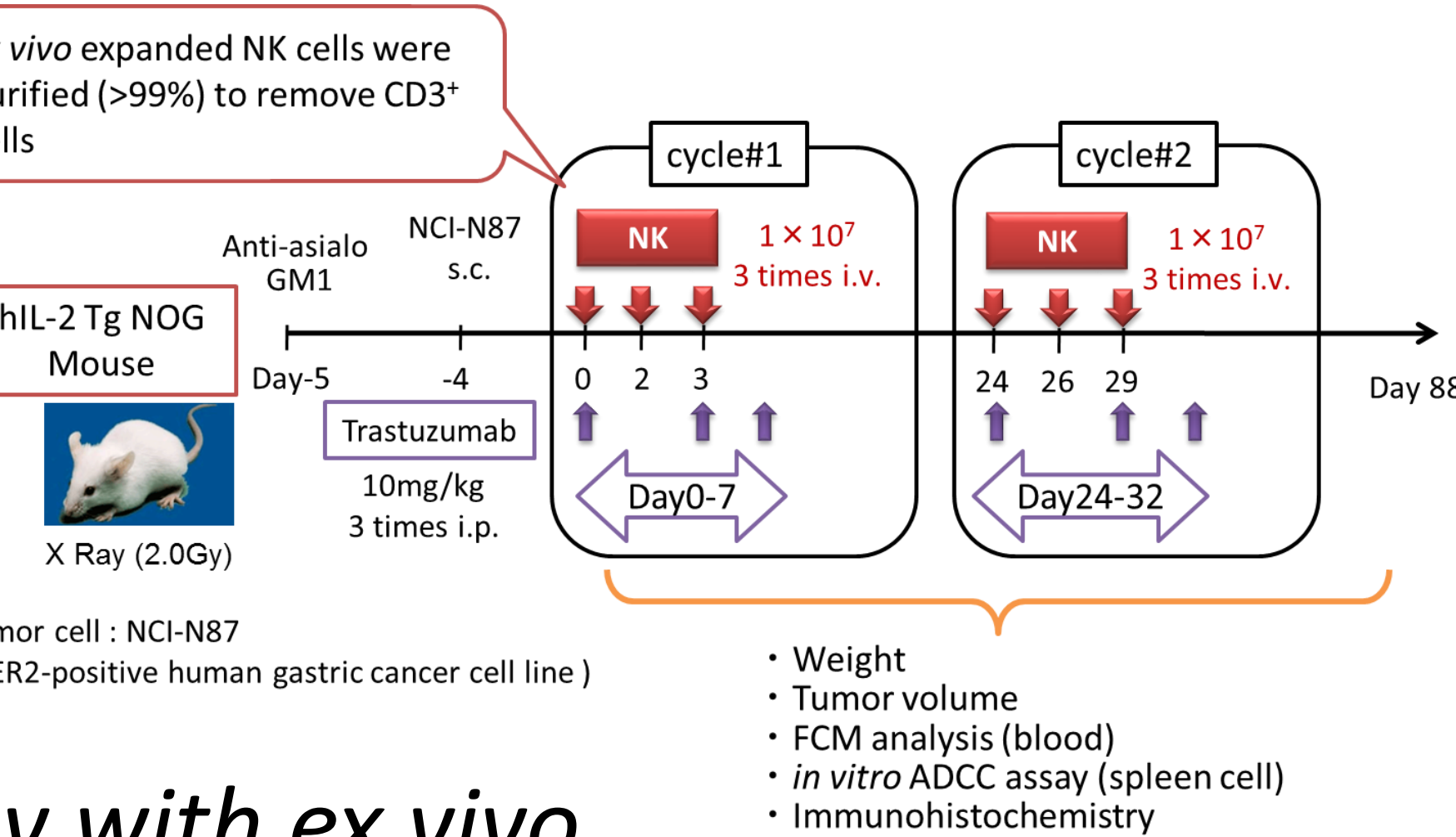
hIL-2 Tg NOG mice: NOG sub-strain expressing human IL-2
(Central Institute for Experimental Animals)



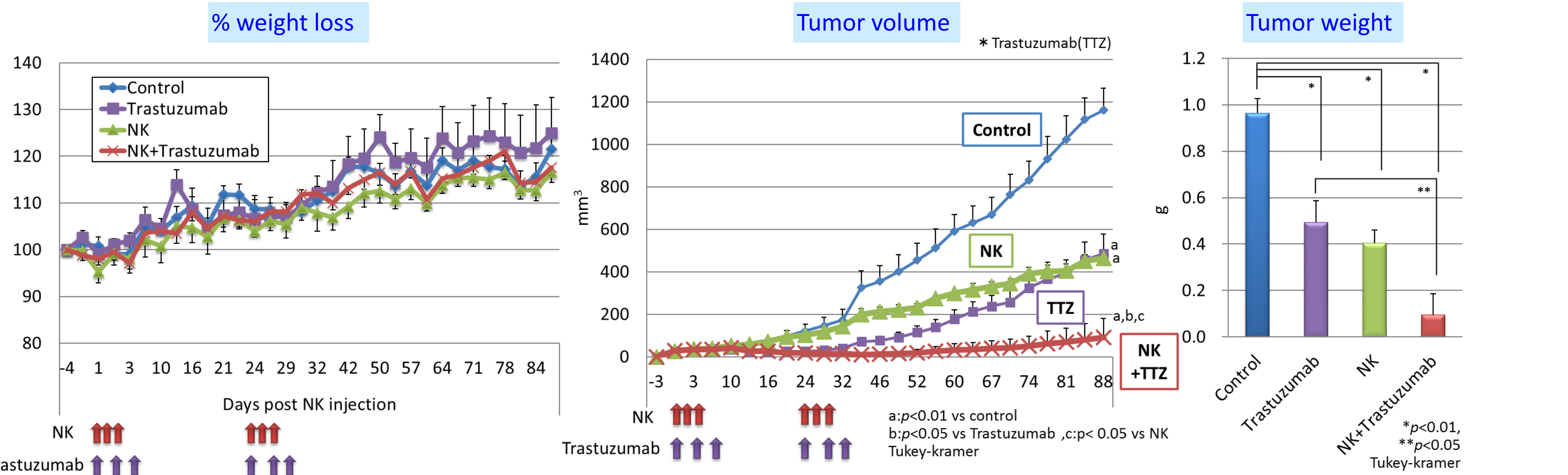
The persistence of *ex vivo* expanded NK cells was remarkably improved by using hIL-2 Tg NOG mice. CD3⁺ cell proportion also increased at day21 and it caused the development of GVHD.

Study design

Group	NK				mAb	n
A	-	-	-	-	-	4
B	-	-	-	-	Trastuzumab (3x2)	4
C	NK(>99%)	Fresh	1x10 ⁷	3x2	-	4
D	NK(>99%)	Fresh	1x10 ⁷	3x2	Trastuzumab (3x2)	4

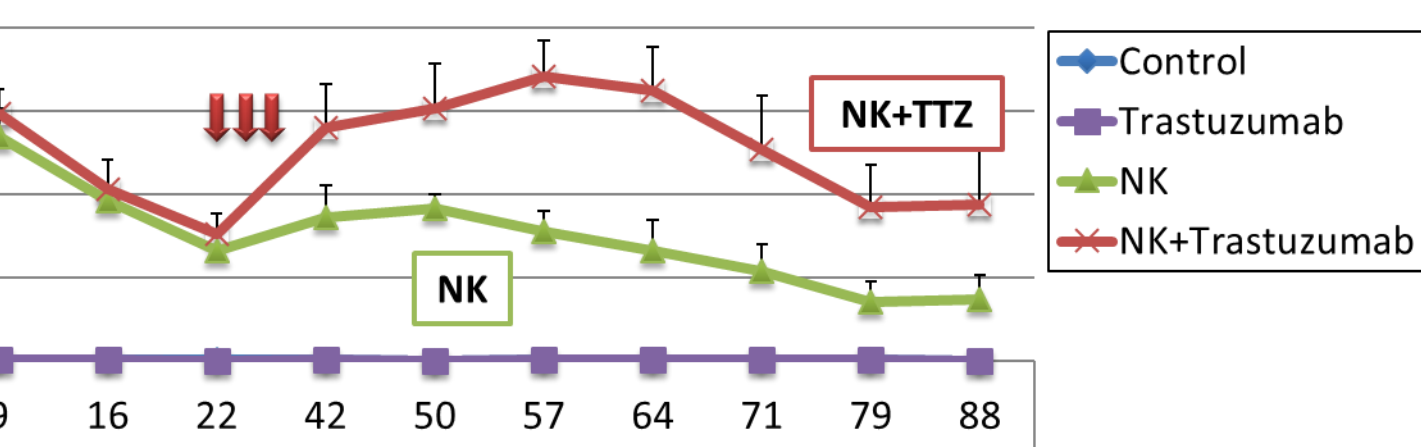


Anti-tumor effect of combination therapy with *ex vivo* expanded human NK cells and Trastuzumab in hIL-2 Tg NOG mice

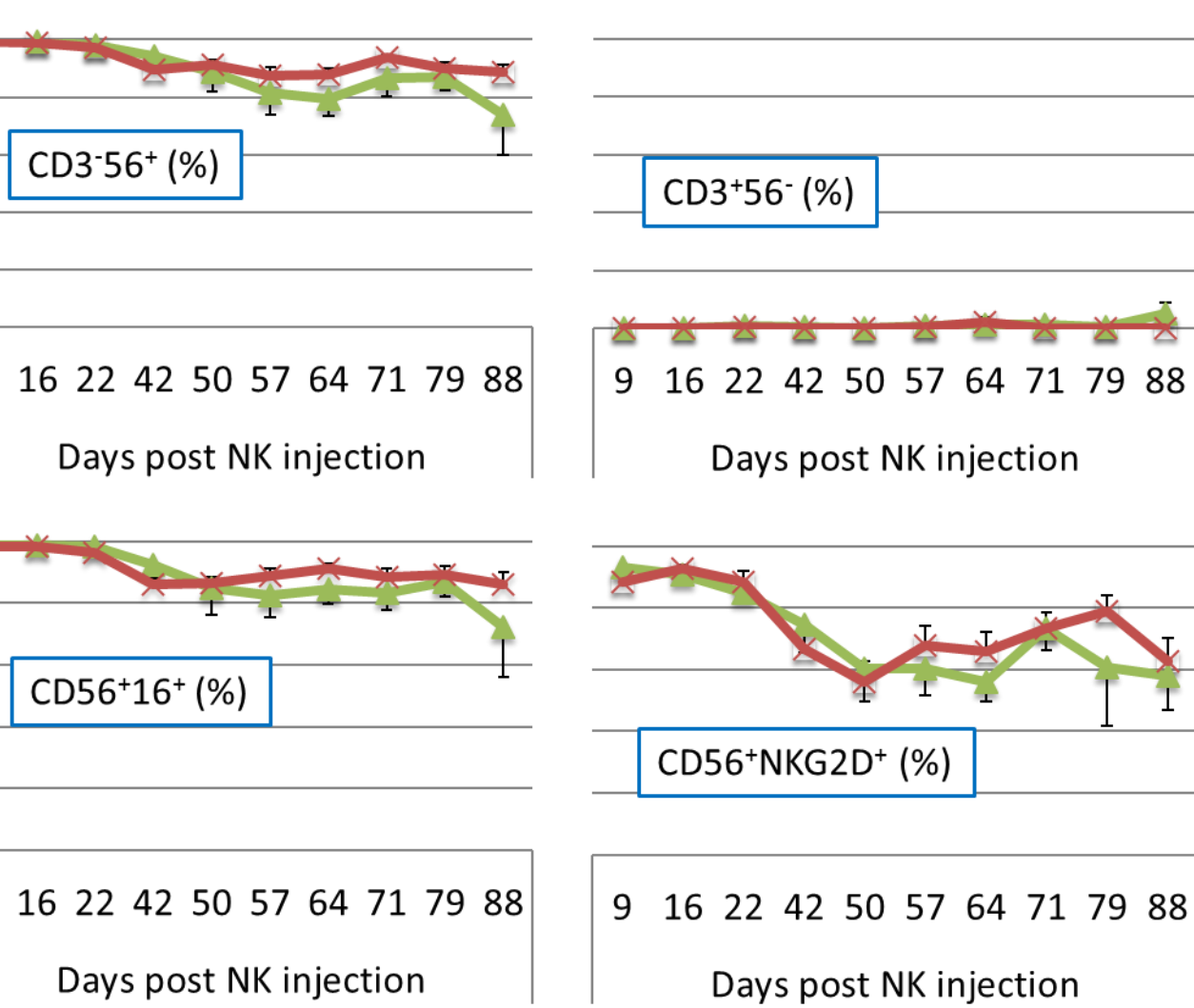


No significant differences in weight loss were observed among the groups.

Chimerism of human CD45⁺ cells in PBL

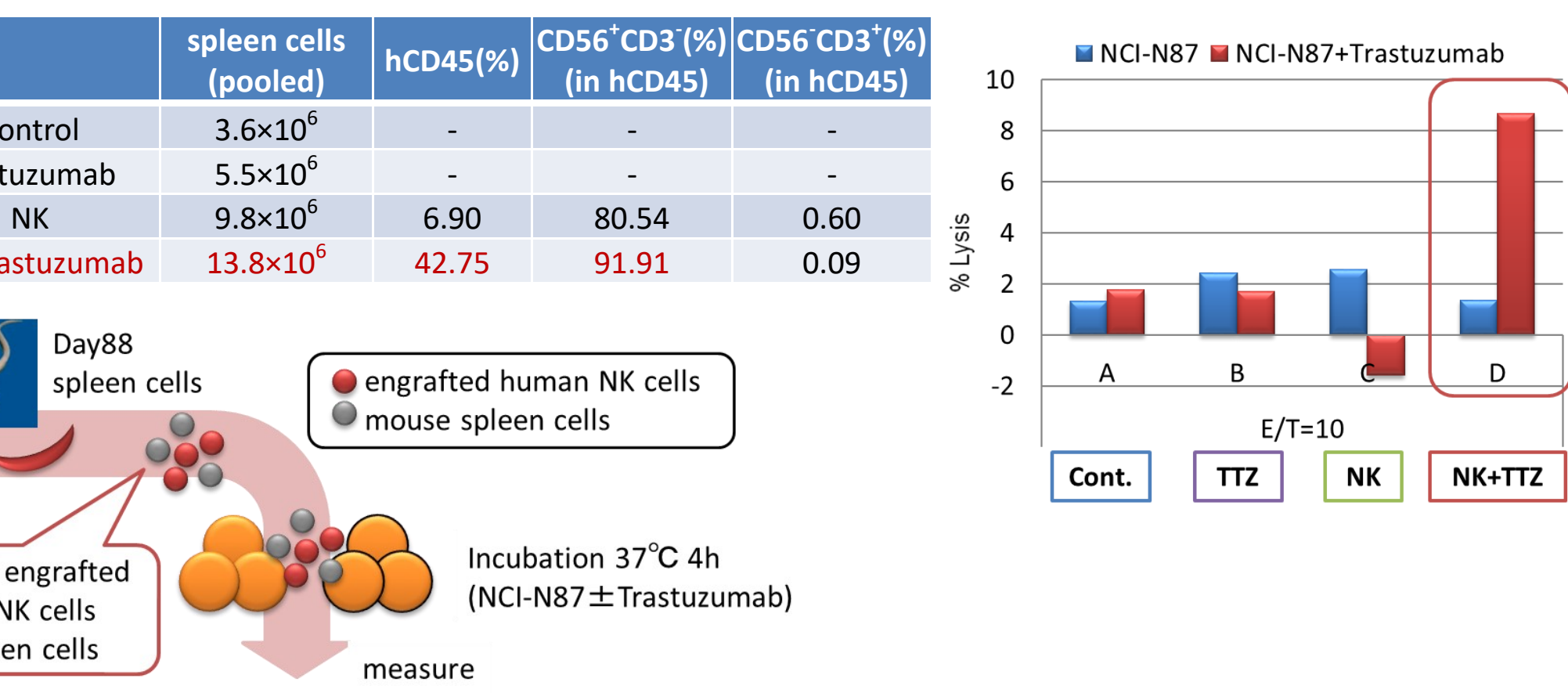


NK-cell surface maker in human CD45⁺ gate (%)



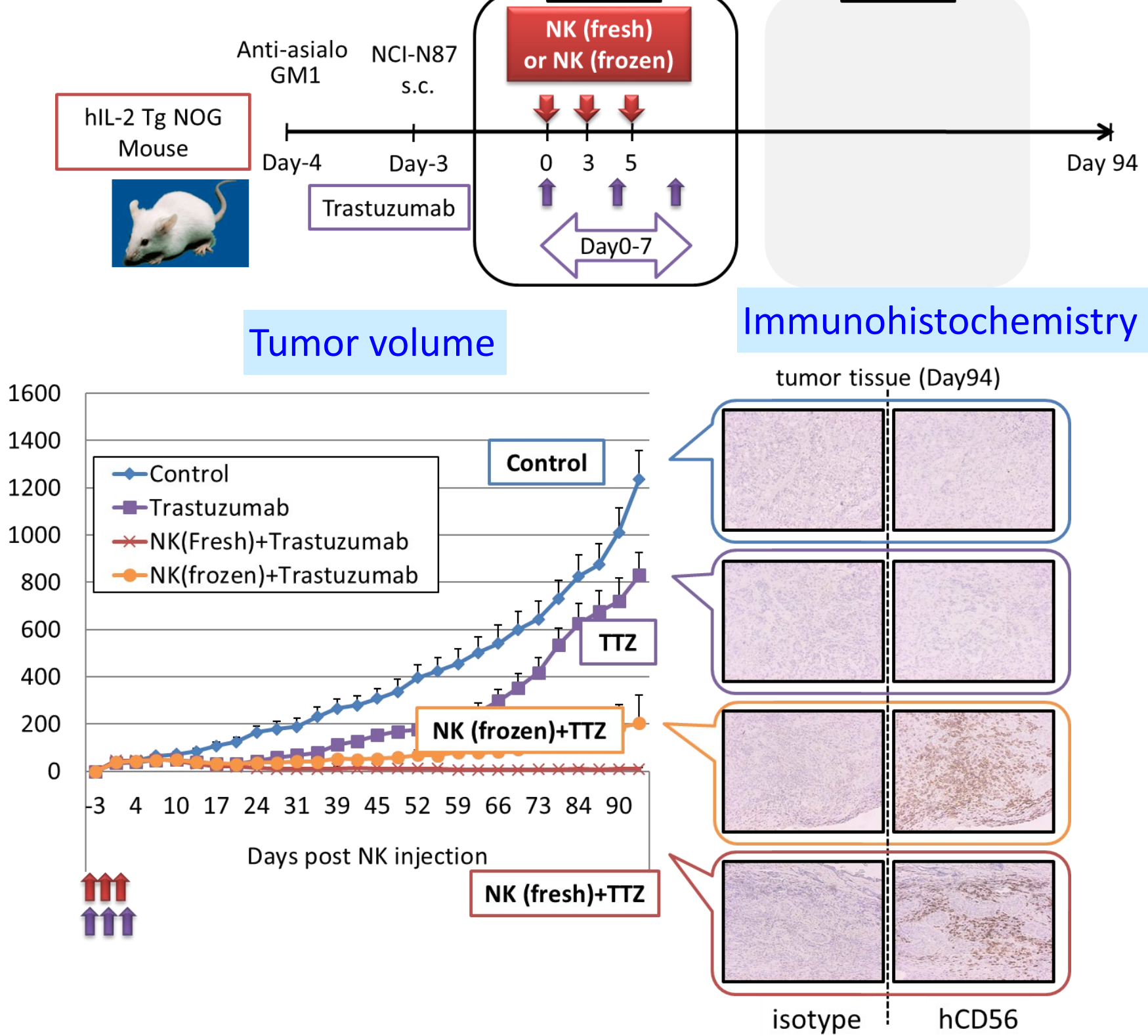
Chimerism of human CD45⁺ in PBL was observed for long period. Functional NK-cell surface markers (CD16, NKG2D) were maintained in human CD45⁺ cells.

in vitro ADCC activity of human NK cells from spleen cells



In combination with Trastuzumab, the infused NK cells were kept potent ADCC activity.

additional experiment



Even in only single cycle of treatment, NK-cell with Trastuzumab significantly reduced tumor volume. Cryopreserved NK cells also showed anti-tumor effect in combination with Trastuzumab. *Ex vivo* expanded human NK cells infiltrated in tumor tissue and they were observed even at 3 months after injection.

Summary

- ✓ High purity of NK cells can be obtained from small amount of peripheral blood by using our expansion system.
- ✓ The expanded NK cells showed *in vivo* as well as *in vitro* anti-tumor activity.
- ✓ Our NK-cell expansion system is highly potent tool to produce a large amount of functional NK cells.