

# Rapid, column-free two-step procedure for the enrichment of human Th17 cells from peripheral blood

M. Fairhurst, J. Yu, and T. E. Thomas  
STEMCELL Technologies Inc., Vancouver, BC, Canada

## Introduction

Human interleukin 17 (IL17) producing CD4<sup>+</sup> T helper (Th17) cells have been identified as a distinct effector T cell subset. They act as key drivers of autoimmune diseases such as rheumatoid arthritis, psoriasis, and inflammatory bowel disease. Phenotypically characterized as CD4<sup>+</sup> T cells expressing CCR6, CCR4, CD161, and IL-23R, they also lack expression of CXCR3. Although their cytokine profile can be heterogeneous, they typically produce IL17 cytokines, are IL17<sup>+</sup>IFN- $\gamma$ <sup>-</sup> and express the lineage-specific transcription factor RORC. A major disadvantage of current isolation methods is the requirement for previous *in vitro* stimulation. We have developed a two-step EasySep<sup>™</sup> immunomagnetic column-free method for the enrichment of CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells from fresh peripheral blood nucleated cells (PBNC). Non-CD4<sup>+</sup> T cells and CXCR3<sup>+</sup> cells are first targeted for depletion using a cocktail of antibody complexes and dextran-coated magnetic particles. Labeled cells are separated using an EasySep<sup>™</sup> magnet, and pre-enriched CD4<sup>+</sup>CXCR3<sup>-</sup> T cells are poured off into a new tube. Next, CCR6<sup>+</sup> cells are positively selected from the pre-enriched fraction. The procedure can be automated using RoboSep<sup>™</sup>. Starting with frequencies of  $5 \pm 2\%$  CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells in fresh PBNC, purities of  $91 \pm 3\%$  (n=16) are obtained. Enriched CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells show increased levels of IL17 cytokine production (minimal IFN- $\gamma$ ) as assessed by ELISA and intracellular staining. Increased RORC mRNA expression is found in the enriched CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells compared to total CD4<sup>+</sup> or CD4<sup>+</sup>CXCR3<sup>+</sup> T cells. Dissecting the role of human Th17 cells in the modulation of immune responses is a requirement for the development of future therapies. These studies will be facilitated by the enrichment of non-activated *in vivo* derived Th17 (CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup>) cells using this easy and rapid enrichment strategy.

## Methods

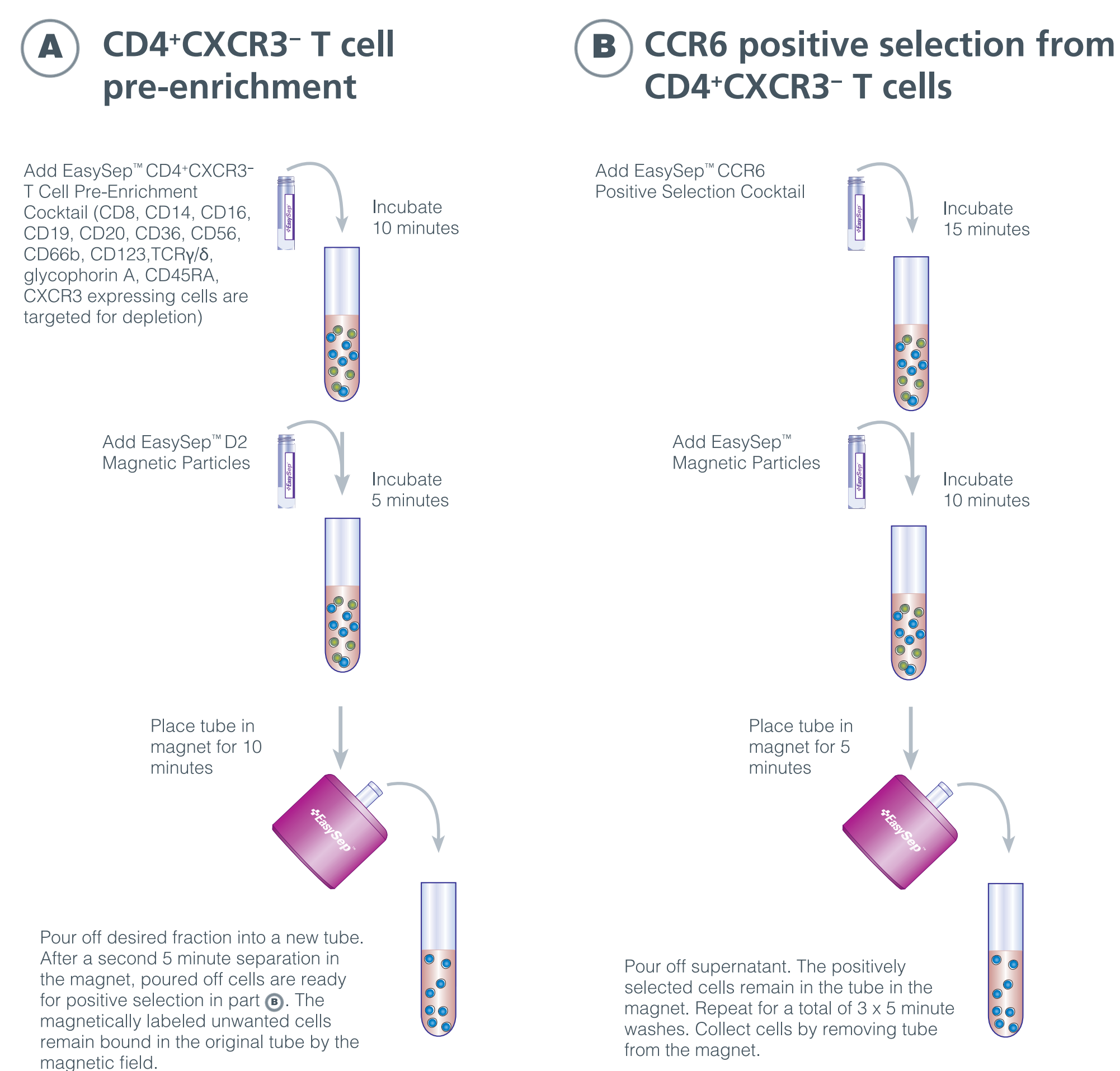
### Preparation of Starting Cell Suspension

A single cell suspension of mononuclear cells (PBMC) was prepared from either fresh whole blood or buffy coat suspensions of peripheral blood using Ficoll-paque PLUS. Alternatively, peripheral blood apheresis (Leucopak PBNC) cells were used following red blood cell lysis and one or more washes to remove platelets. Use of fresh ( $\leq 24$  hours) cells is recommended. Typically start cells were resuspended at  $5 \times 10^7$  cells/mL in PBS + 2% FBS and 1mM EDTA.

### Assessing Enriched Cells

The purity of Th17 (CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup>) T cells can be measured by flow cytometry after staining with fluorochrome-conjugated anti-CD4, anti-CXCR3, and anti-CCR6 antibodies. In addition, intracellular staining of IL-17 cytokine was assessed after stimulation of cells with PMA-Ionomycin. For ELISA, enriched cells were stimulated with anti-CD3/anti-CD28 beads for 24, 48, and 72 hours. Supernatants were collected and analysed for secreted cytokines.

## FIGURE 1: Two-step EasySep<sup>™</sup> procedure for column-free enrichment of Th17 (CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup>) cells from human peripheral blood



## Results

TABLE 1: Purity and recovery of human Th17 (CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup>) cells enriched from peripheral blood by manual EasySep<sup>™</sup> or RoboSep<sup>™</sup>

n	% start	% purity enriched	Typical recovery of CD4 <sup>+</sup> CXCR3 <sup>-</sup> CCR6 <sup>+</sup> T cells from start
16	4.7 $\pm$ 1.8	91.3 $\pm$ 3.7	1.5x10 <sup>6</sup> CD4 <sup>+</sup> CXCR3 <sup>-</sup> CCR6 <sup>+</sup> T cells from 1x10 <sup>8</sup> peripheral blood nucleated cells

Purities determined by flow cytometry. Values are expressed as means  $\pm$  SD.

FIGURE 2: Phenotypic assessment of human Th17 (CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup>) cells enriched using EasySep<sup>™</sup>

Start: 5.4% CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T Cells

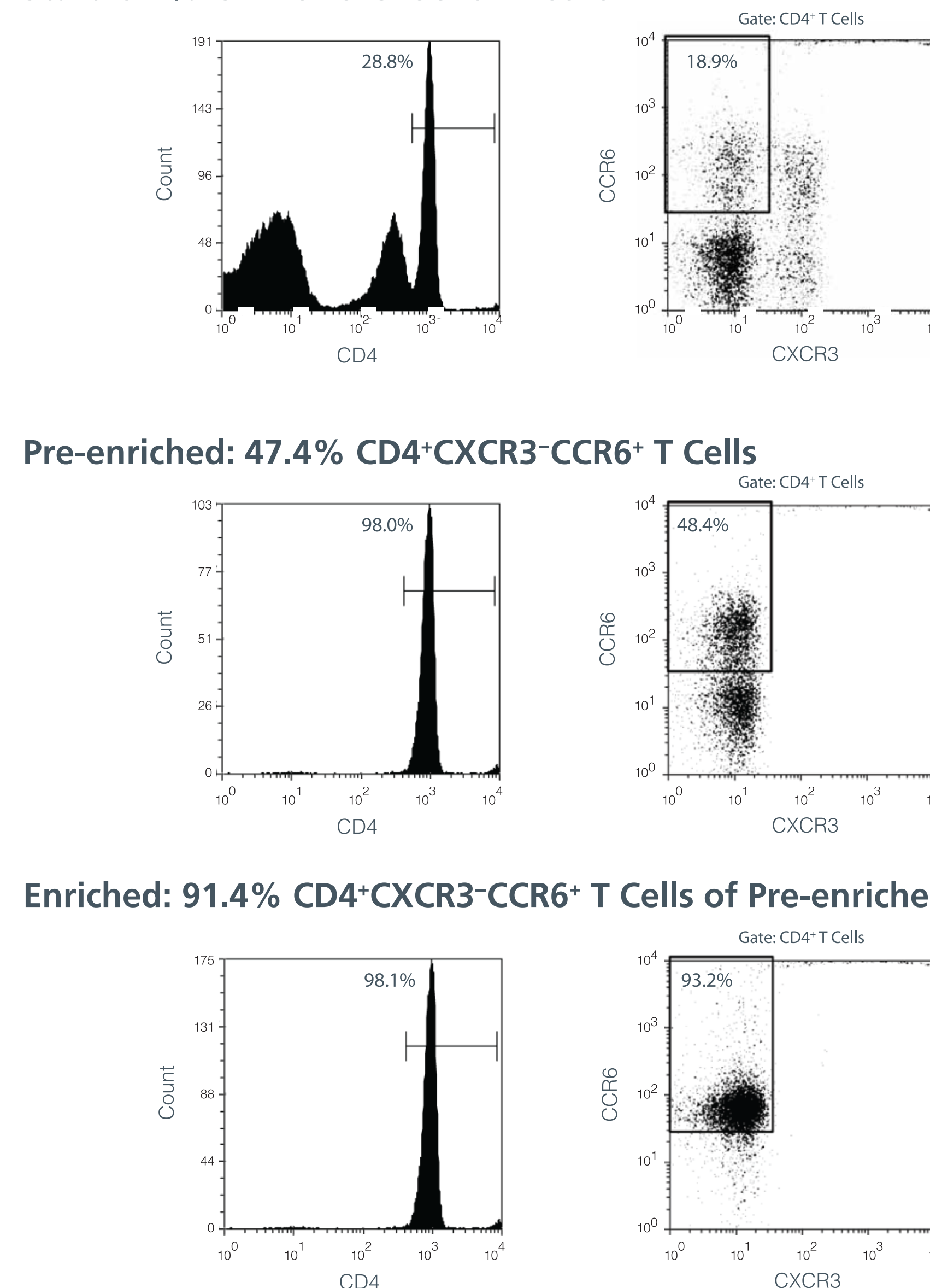
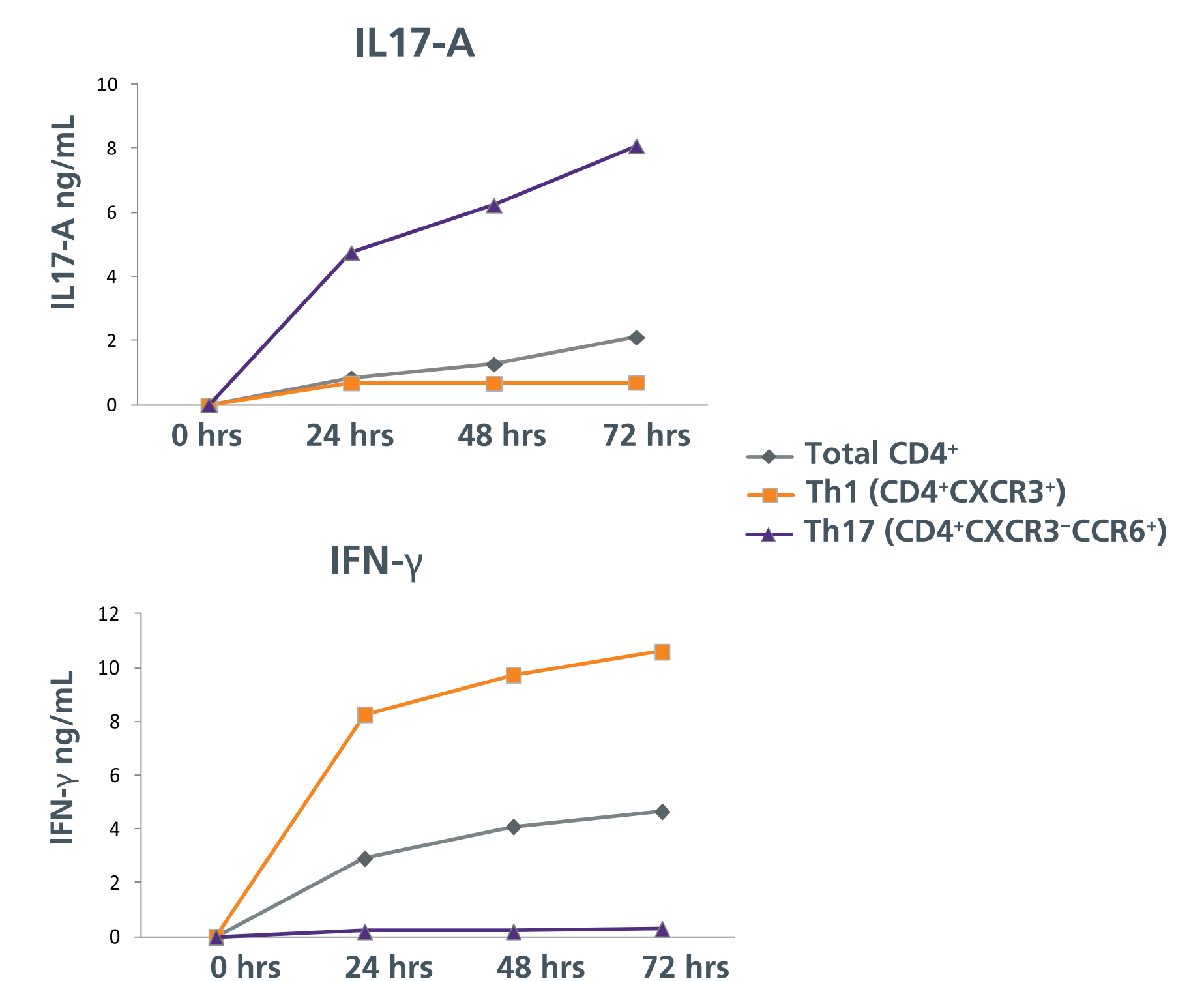
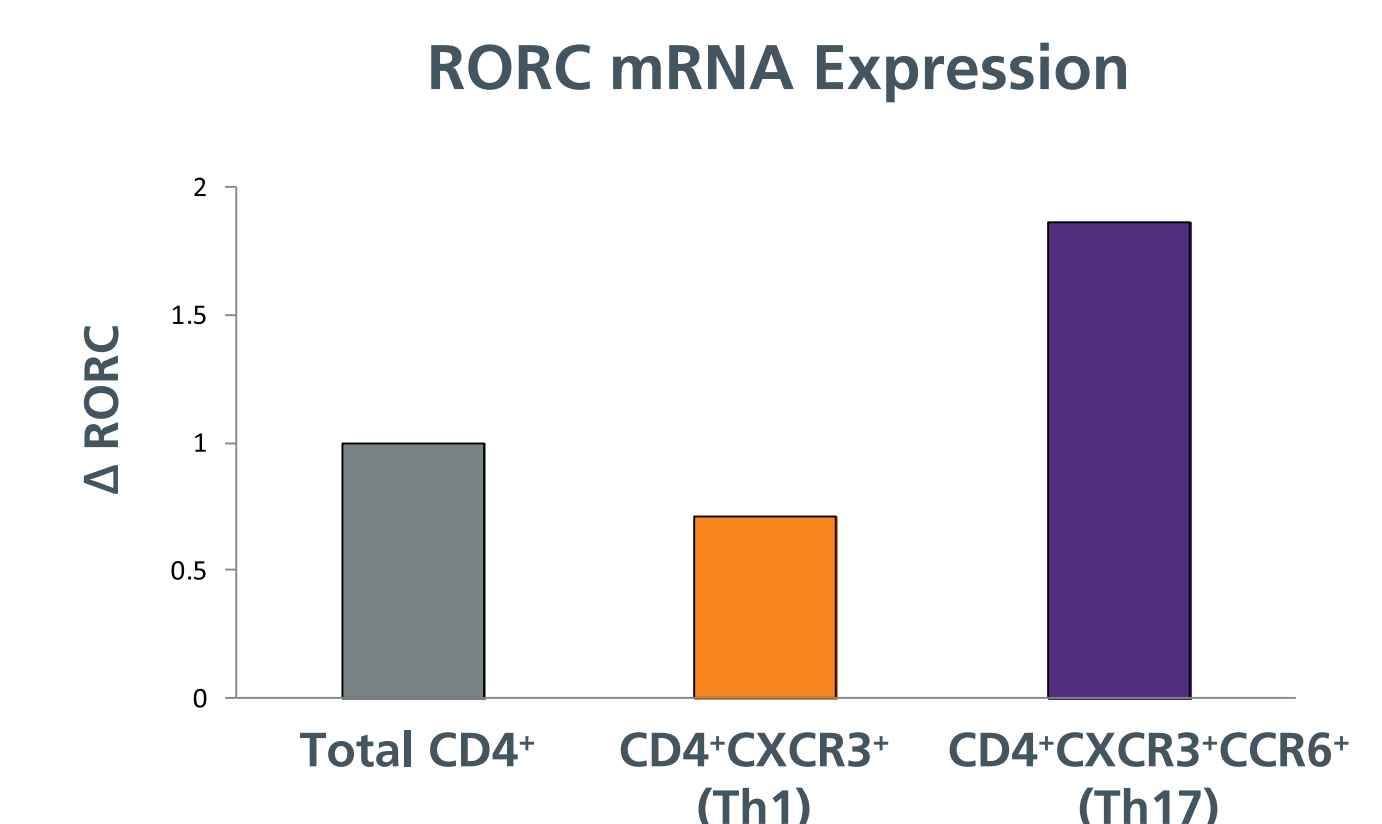


FIGURE 3: Enriched CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells show increased levels of IL17-A secretion by ELISA with minimal IFN- $\gamma$  secretion compared to CD4<sup>+</sup>CXCR3<sup>+</sup> T cells



Freshly isolated CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells were stimulated with anti-CD3/anti-CD28 coated beads at a ratio of 8:1 (cells: beads). Supernatants were collected at 24, 48, and 72 hours and cytokines were assessed by ELISA. Results for one of three comparable experiments are shown. A similar trend was observed with intracellular staining.

FIGURE 4: Enriched CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells show an increased level of RORC mRNA



Freshly isolated CD4<sup>+</sup>CXCR3<sup>-</sup>CCR6<sup>+</sup> T cells were stimulated with anti-CD3/anti-CD28 beads at a ratio of 8:1 (cells: beads) for 48 hours. Expression of RORC relative to that of 18S was analysed by quantitative RT-PCR.