

**Description**

The untransduced T cells are produced by mock lentiviral transduction of human primary CD4+CD8+ T cells. These cells are subjected to comparable manipulations as CAR-T cells: activation, spinoculation (without lentivirus), and expansion. These T cells are meant to be negative controls in experiments using lentivirus-transduced primary CAR-T cells.

**Application**

Negative control for lentivirus-transduced CAR-T cells

**Materials Provided**

Components	Format
One vial of frozen cells	Each vial contains $2 \times 10^6$ cells in 1 ml of CryoStor® CS10)

**Mycoplasma Testing**

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

**Storage Conditions**

Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a  $-80^{\circ}\text{C}$  freezer for long term storage. Contact technical support at support@bpsbioscience.com if the cells are not frozen in dry ice upon arrival.

**Recommended T Cell Medium:** StemSpan SFEM (Stemcell Technologies, #09650) supplemented with 10% heat-inactivated FBS (Life Technologies, #10082147), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01), plus 10 ng/ml IL-2 (BPS Bioscience, #90184)

**Cell Thawing and Culture Protocol:**

1. Swirl the vial of frozen cells for approximately 60 seconds in a  $37^{\circ}\text{C}$  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 T cell growth medium.

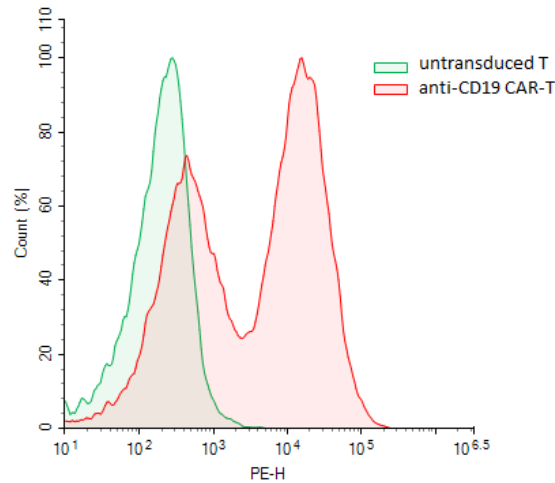
**Leaving the cells in the water bath at  $37^{\circ}\text{C}$  for too long will result in rapid loss of viability.**

2. Immediately spin down the cells at  $300 \times g$  for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed T cell growth medium.
3. Transfer the resuspended cells to a T25 flask and add T cell activation reagents. Activate the cells at  $37^{\circ}\text{C}$  with 5%  $\text{CO}_2$  for 24 – 48 hours.
4. Centrifuged the cells gently at  $300 \times g$  for 5 minutes and resuspend in fresh T cell medium. Continue to culture the cells at  $37^{\circ}\text{C}$  with 5%  $\text{CO}_2$ . Do not allow the cell density to exceed  $2.0 \times 10^6$  cells/ml. Transfer the cells in larger culture vessels and add fresh medium when the density reaches  $2.0 \times 10^6$  cells.



It is recommended to activate the untransduced T cells for expansion after thawing. Since these are primary cells, the extent of expansion is not predictable. Perform the cytotoxicity assay as soon as possible to avoid exhaustion. The untransduced T cells should not be in culture for more than 8-10 days. It is not recommended to freeze the cells again once they have been activated and expanded.

## Validation



*Figure 2: The expression of anti-CD19 CAR in anti-CD19 CAR-T cells, with untransduced T cells as negative control. Anti-CD19 CAR-T cells (BPS Bioscience #78171) and untransduced T cells were thawed and activated for 48 hours. Anti-CD19 CAR-T and untransduced T cells were then expanded for another 4 days, and ~50,000 cells were analyzed by flow cytometry using PE-anti-FMC63 ScFv (Acrobiosystems, #FM3-HPY53-25tests). Untransduced T cells were used as negative control.*

## Warnings

Donors have been screened and determined negative for:

- Hepatitis B (anti-HBc EIA, HBsAg EIA)
- Hepatitis C (anti-HCV EIA)
- Human Immunodeficiency Virus (HIV-1/HIV-2 plus O)
- Human T-Lymphotropic Virus (HTLV-I/II)
- HIV-1/HCV/HBV
- West Nile Virus
- Trypanosoma cruzi

**Note:** Testing cannot guarantee that any sample is completely virus-free. These cells should be treated as potentially infectious and appropriate biological safety level 2 precautions should be used.

## Troubleshooting Guide

For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

**Related Products**

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
Anti-CD19 CAR-T Cells	78171	1 vial
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ)	78600	50 μl
Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-CD28-CD3ζ)	78603	50 μl
Anti-CD20 CAR Lentivirus (Clone Leu-16 ScFv-CD8-4-1BB-CD3ζ)	78606	50 μl
Anti-CD22 CAR Lentivirus (Clone m971 ScFv-CD8-4-1BB-CD3ζ)	78608	50 μl
Anti-CD19/CD22 Bispecific CAR Lentivirus (Clones FMC63/m971 ScFv-CD8-4-1BB-CD3ζ)	78609	50 μl