

Description

The T Cell Receptor (TCR) is found on the surface of T-cells and is responsible for recognizing antigens bound to MHC (Major Histocompatibility Complex) molecules. Activation of the TCR results in activation of downstream NFAT signaling. The TCR consists of a heterodimer of two different protein chains, of which the alpha (α) and beta (β) chains are the predominant chains.

The TCR CRISPR Lentiviruses are replication incompetent, HIV-based VSV-G pseudo-typed lentiviral particles that are ready to be transduced into almost all types of mammalian cells, including primary and non-dividing cells. The particles contain a CRISPR/Cas9 gene driven by an EF1A promoter, along with 4 sgRNA (single guide RNA) targeting human TRAC (T-Cell Receptor Alpha Constant) and human TRBC1 (T-Cell Receptor Beta Constant 1) regions of the α and β chains.

The DNA transduced by the integrating lentivirus integrates randomly into the cellular genome to express both Cas9 and sgRNA. Puromycin selection increases the knockout efficiency by forcing high expression levels of both Cas9 and the sgRNA, and can be used with the integrating lentivirus to quickly and easily achieve high knockdown efficiencies in a cell pool. Efficiencies also depend on the cell type and the gene of interest.

Applications

1. Transient knock-down of TCR in target cells.
2. Generation of a stable TCR knock-out cell line following puromycin selection.

Formulation

The lentivirus particles were produced from HEK293T cells. They are supplied in cell culture medium containing 90% DMEM + 10% FBS.

Titer

Two vials of lentivirus (500 μ l x 2) at a titer $\geq 1 \times 10^6$ TU/ml. The titer will vary with each lot; the exact value is provided with each shipment.

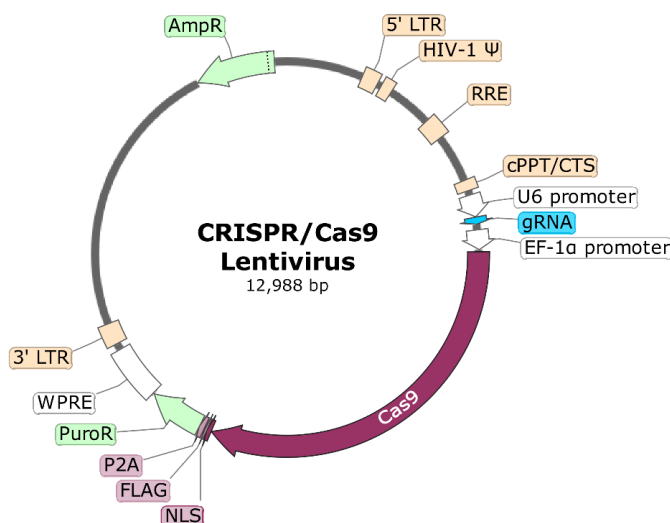


Figure 1: Schematic of the Lenti-vector used to generate the TCR CRISPR/Cas9 Lentivirus.

Gene Target:	Primer ID:	sgRNA Sequence:
TRAC	TCR-1	AGAGTCTCTCAGCTGGTACA
TRAC	TCR-2	TGTGCTAGACATGAGGTCTA
TRBC1	TCR-3	GGAGAATGACGAGTGGACCC
TRBC1	TCR-4	GCAGTATCTGGAGTCATTGA

Figure 2: List of sgRNA Sequences in the TCR CRISPR/Cas9 Lentivirus.

Storage Conditions



Lentiviruses are shipped with dry ice. For long term storage, it is recommended to store the lentiviruses at -80°C. Avoid repeated freeze-thaw cycles. Titers can drop significantly with each freeze-thaw cycle.

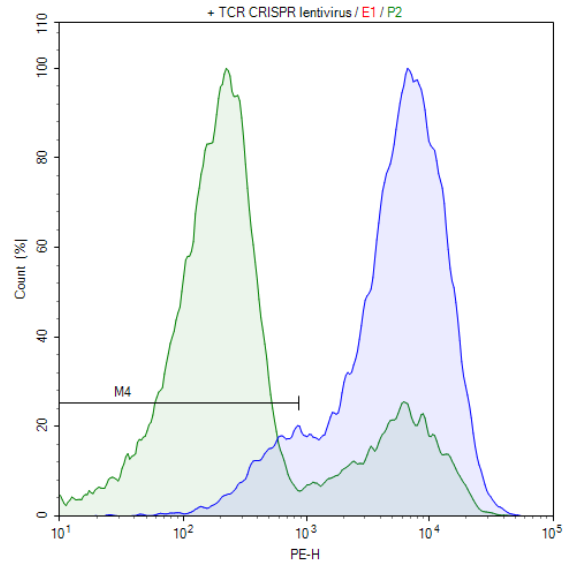
Biosafety



The lentiviruses are produced with the SIN (self-inactivation) lentivector which ensures self-inactivation of the lentiviral construct after transduction and after integration into the genomic DNA of the target cells. None of the HIV genes (*gag*, *pol*, *rev*) will be expressed in the transduced cells, as they are expressed from packaging plasmids lacking the packing signal and are not present in the lentivirus particle. Although the pseudotyped lentiviruses are replication-incompetent, they require the use of a Biosafety Level 2 facility. BPS Bioscience recommends following all local federal, state, and institutional regulations and using all appropriate safety precautions.

Validation Results

A.

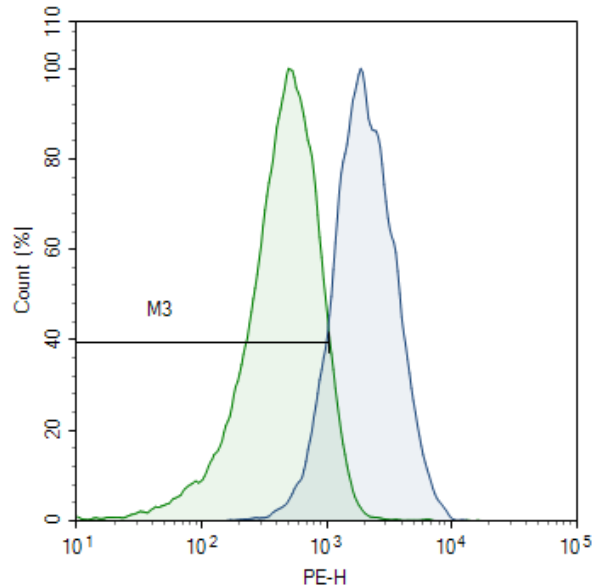


B.

Sample	Gate #	% of Cell Population
Jurkat Parental cells	M4	9.79%
Jurkat cells transduced with TCR CRISPR/Cas9 Lentivirus	M4	75.30%

Figure 3. Knock-down of TCR in Jurkat cells. **A.** Jurkat cells were transduced via spinoculation with 5,000,000 TU/well of TCR CRISPR/Cas9 lentivirus, corresponding to an MOI of approximately 5-10. 72 hours after transduction, cells were stained with PE anti-human TCR antibody (BioLegend, #306708) and analyzed by flow cytometry. Parental Jurkat cells are shown in blue, and the transduced cells are shown in green. **B.** Percentages of cell populations captured in Gate M4.

A.



B.

Sample	Gate #	% of Cell Population
Primary T cells	M3	10.94%
Primary T transduced with TCR CRISPR/Cas9 Lentivirus	M3	90.47%

Figure 4: Knock-down of TCR in primary T cells using TCR CRISPR/Cas9 Lentivirus. A. Primary T cells (1 million cells per well) were transduced via spinoculation with 5,700,000 TU/well of TCR CRISPR/Cas9 lentivirus, corresponding to an MOI ≥ 50 . 48 hours after transduction, cells were stained with PE-labeled anti-human TCR antibody (BioLegend, #306708) and analyzed by flow cytometry. Non-transduced primary T cells are shown in blue, and the transduced cells are shown in green. **B.** Percentages of cell populations captured in Gate M3.

Notes

The CRISPR/CAS9 technology is covered under numerous patents, including U.S. Patent Nos. 8,697,359 and 8,771,945, as well as corresponding foreign patents applications, and patent rights.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
TCR CRISPR/Cas9 Lentivirus (Non-Integrating)	78062	500 µl x 2
PD-1 CRISPR/Cas9 Lentivirus (Integrating)	78052	500 µl x 2
PD-1 CRISPR/Cas9 Lentivirus (Non-Integrating)	78059	500 µl x 2
Cas9, His-tag (<i>S. pyogenes</i>)	100206	50 µg
TCR Knockout NFAT-Luciferase Reporter Jurkat Recombinant Cell Line	79887	2 vials
TCR Activator - Raji Cell Line	60556	2 vials
TCR Activator – CHO Recombinant Cell Line	60539	2 vials