

Description

The TCR Knockout Jurkat cell line was generated by CRISPR/Cas9 genome editing to remove the TRAC (T-Cell Receptor Alpha Constant) and TRBC1 (T-Cell Receptor Beta Constant 1) domains of the TCR α and β chains.

Background

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. Engagement of the TCR initiates downstream NFAT signaling resulting in T cell activation. The TCR consists of a heterodimer of two different protein chains, of which the alpha (α) and beta (β) chains are the predominant chains. CRISPR/Cas9 genome editing was utilized to remove the TRAC (T-Cell Receptor Alpha Constant) and TRBC1 (T-Cell Receptor Beta Constant 1) regions of the α and β chains, resulting in a loss of TCR expression.

Application

Use as a control when generating or characterizing CAR-T Cells

Materials Provided

Components	Format
2 vials of frozen cells	Each vial contains 2×10^6 cells in 1 ml of cell freezing medium (BPS Bioscience #79796)

Parental Cell Line

Jurkat (clone E6-1), human T lymphoblast, suspension

Mycoplasma Testing

The cell line has been screened to confirm the absence of Mycoplasma species.

Materials Required but Not Supplied

These materials are not supplied with the cell line but are necessary for cell culture and cellular assays. BPS Bioscience's reagents are validated and optimized for use with this cell line and are highly recommended for best results. Media components are provided in the Media Formulations section below.

Media Required for Cell Culture

Name	Ordering Information
Thaw Medium 2	BPS Bioscience #60184

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°C freezer for long term storage. Contact technical support at support@bpsbioscience.com if the cells are not frozen in dry ice upon arrival.

Media Formulations

For best results, it is *highly recommended* to use these validated and optimized media from BPS Bioscience. Other preparations or formulations of media may result in suboptimal performance.



Note: Thaw Media do not contain selective antibiotics. Cells should be grown at 37°C with 5% CO_2 . BPS Bioscience's cell lines are stable for at least 15 passages when grown under proper conditions.

Media Required for Cell Culture

Thaw Medium 2 (BPS Bioscience #60184):

RPMI 1640 medium supplemented with 10% FBS, 1% Penicillin/Streptomycin.

Cell Culture Protocol

Cell Thawing

1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°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 Thaw Medium 2.
Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.
2. Immediately spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed Thaw Medium 2.
3. Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO₂ incubator.
4. After 24 hours of culture, check for cell viability. For a T25 flask, add 3-4 ml of Thaw Medium 2, and continue growing in a 5% CO₂ incubator at 37°C until the cells are ready to passage.
5. Cells should be passaged before they reach a density of 2 x 10⁶ cells/ml. At first passage and subsequent passages, use Thaw Medium 2.

Cell Passage

Dilute the cell suspension into new culture vessels before they reach a density of 2 x 10⁶ cells/ml, at no less than 0.2 x 10⁶ cells/ml of Thaw Medium 2. The sub-cultivation ratio should maintain the cells between 0.2 x 10⁶ cells/ml and 2 x 10⁶ cells/ml.

Cell Freezing

1. Spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cell pellet in 4°C Freezing Medium (BPS Bioscience #79796, or 10% DMSO + 90% FBS) at a density of ~2 x 10⁶ cells/ml.
2. Dispense 1 ml of cell aliquots into cryogenic vials. Place the vials in an insulated container for slow cooling and store at -80°C overnight.
3. Transfer the vials to liquid nitrogen the next day for storage.



Note: It is recommended to expand the cells and freeze at least 10 vials at an early passage for future use.

Validation Data

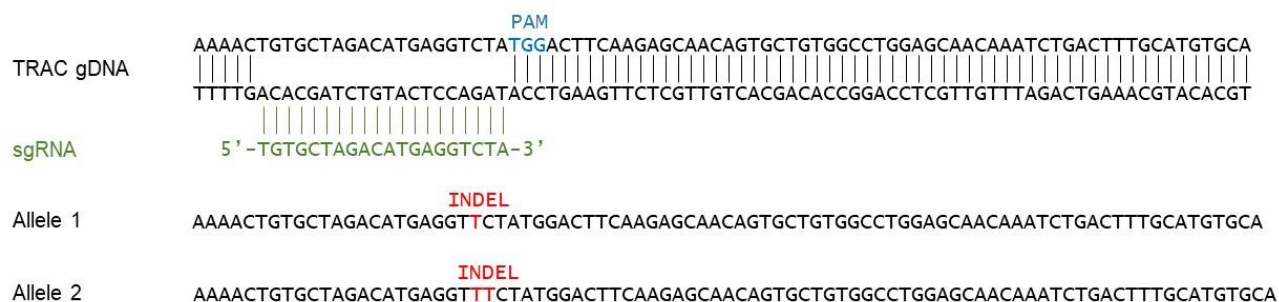


Figure 1. Genomic Sequencing of TRAC in the TCR Knockout Jurkat cells.

Genomic DNA from the TCR Knockout Jurkat cells was isolated and sequenced. The PAM (Protospacer Adjacent Motif) is shown in blue, the sgRNA (synthetic guide RNA) in green, and the Indels (Insertions / Deletions) in the two TRAC alleles are indicated in red.



Figure 2. Genomic Sequencing of TRBC1 in the TCR Knockout Jurkat cells.

Genomic DNA from the TCR Knockout Jurkat cells was isolated and sequenced. The PAM (Protospacer Adjacent Motif) is shown in blue, the sgRNA (synthetic guide RNA) in green, and the Indels (Insertions / Deletions) in the two TRBC1 alleles are indicated in red.

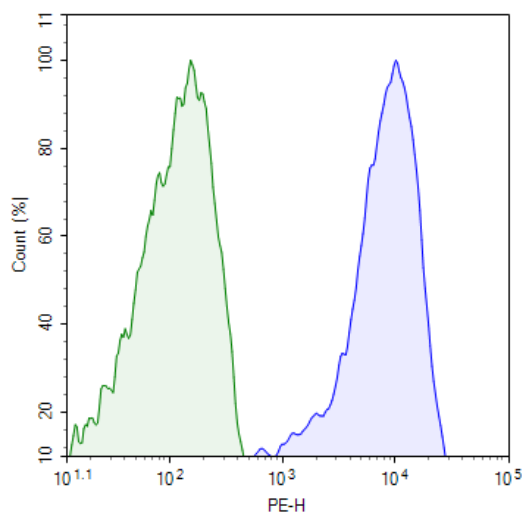


Figure 3. Expression of TCR in the TCR Knockout Jurkat cells.

TCR Knockout Jurkat cells (green) or parental Jurkat cells (blue) were stained with PE-labeled anti-human TCR α/β antibody (BioLegend #306707) and analyzed by flow cytometry. Y-axis is the % cell number. X-axis is the intensity of PE.

Sequence

Human mRNA for T-Cell Receptor Alpha Chain (GenBank Accession #X02592.1), with the sgRNA targeting sequence underlined:

atgctcctgctgctcctcagtgctcgaggtgattttaccctgggaggaaccagagcccagtcggtgaccagcttggcagccagctctgctctgaaggagc
 cctggttctgctgaggtgcaactactcatcgtctgttccacatatctctctggtatgtgcaataccccaaccaaggactccagcttctcctgaagtacacatcagcg
 gccaccctggtaaaggcatcaacggttttagggctgaatttaagaagagtgaacctctccacctgacgaaacctcagcccatatgagcgcgaggctgagt
 acttctgtgctgtagtgatctcgaaccgaacagcagtgctccaagataatcttggatcaggaccagactcagcatccggccaatatccagaacctgacc
 tgcctgtaccagctgagagactctaaatccagtgacaagtctgtctgctattcaccgattttgatttcaaaacaatgtgtcaaaagtaaggattctgatgtga
 tatcacagacaaaactgtgctagacatgaggtctatggacttcaagcaacagtgctgtggcctggagcaaaaatctgactttgcatgtgcaaacgcttcaac
 aacagcattattccagaagacaccttctccccagcccagaaagttcctgtgatgtcaagctggtcgagaaaagcttgaacagatacgaacctaaacttcaaa
 acctgtcagtgattgggtccg aatcctcctcctgaaagtgccgggttaatctgctcatgacgctgaggctgtggtccagctga

Human mRNA for T-Cell Receptor Beta Chain (GenBank Accession #NG_001333), with the sgRNA targeting sequence underlined:

atgggctgcaggctgctctgctgtgctggttctctgctcctgggagcagttccatagacactgaagttaccagacacaaaaacacctggctatgggaatgaaa
 ataagaagtcttgaatgtgaacaacatatggggcagggctatgtattggtacaagcagaaagtaagaagccaccggagctcatgtttgtctacagctatg
 agaaactctataaatgaaagtgtgccaagtgccttctcacctgaatgcccacagctctctctaaaccttcacctacacgcccctgcagccagaagactcagcc
 ctgtatctctgcgcagcagccccctcgggagggaggactgaacactgaagcttcttggacaaggcaccagactcacagttgtagaggacctgaacaaggtg
 ttccacccgaggtcgtgtgtttgagccatcagaagcagagatctccacacccaaaaggccacactggtgtgcttggccacaggcttctccccgaccgctgg
 agctgagctggtgggtgaatgggaaggaggtgcacagtggggtcagcacggaccgagcccctcaaggagcagcccctcaatgactccagatactgcctg
 agcagccgctgaggtctcggccaccttctggcagaacccccgcaaccacttccgctgtcaagtccagttctacgggctctcggagaatgacgagtgaccag
 gataggccaaacctgcaccagatcgtcagcgcggagcctggggtagagcagactgtggtttacctcgggtcctaccagcaaggggtcctgtctgccacc
 atcctctatgagatcctgtaggaaggccacctgtatgctgtgctggtcagcgccttgtgtgtagggccatggtcaagagaaaggatttctga

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.

License Disclosure

Visit bpsbioscience.com/license for the label license and other key information about this product.

Troubleshooting Guide

Visit bpsbioscience.com/cell-line-faq for detailed troubleshooting instructions. For all further questions, please email support@bpsbioscience.com.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
TCR Knockout NFAT-Luciferase Reporter Jurkat Recombinant Cell Line	79887	2 vials
TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78364	2 vials
TCR Activator - Raji Cell Line	60556	2 vials