

**Description**

This cell line is a clonal derivative from the Cas9-Expressing Jurkat Cell Pool (BPS Bioscience, #78070). It was generated by limiting dilution of the original pool and isolation of individual clones, which were screened based on Cas9 expression to obtain a high-expressing cell line. The expressed Cas9 protein includes a C-terminal FLAG tag.

**Background**

Cas9 (*Streptococcus pyogenes* CRISPR associated protein 9) is an endonuclease enzyme that, when recruited to a specific DNA sequence by the sgRNA (single guide RNA), introduces a double stranded break into the DNA. This double stranded break is repaired in mammalian cells either through Non-Homologous End Joining or Homologous Recombination. Non-Homologous End Joining often results in the deletion or insertion of several base pairs at the cut site, which, when resulting in a frameshift, causes the functional inactivation of the targeted gene. Cas9-expressing Jurkat cell lines can be transduced or electroporated with sgRNA targeting a gene of interest to quickly generate knock-out cell pools or cell lines.

**Application**

1. Quickly generating knock-out cell pools or cell lines in Jurkat cells.
2. Implementing sgRNA screens in Cas9 expressing Jurkat cells.

**Materials Provided**

Components	Format
2 vials of frozen cells	Each vial contains $2 \times 10^6$ cells in 1 ml of 10% DMSO

**Host Cell**

Jurkat is a human leukemia cell line. Non-adherent T lymphocytes.

**Mycoplasma Testing**

This cell line has been screened using the MycoAlert™ Mycoplasma Detection Kit (Lonza, #LT07-118) to confirm the absence of Mycoplasma contamination. MycoAlert Assay Control Set (Lonza, #LT07-518) was used as a positive control.

**Materials Required but Not Supplied**

These materials are not supplied with this cell line but are necessary for cell culture and cellular assays. BPS Bioscience reagents systems 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.

**Materials Required for Cell Line Culture**

Name	Ordering Information
Thaw Medium 2	BPS Bioscience #60184
Growth Medium 2K	BPS Bioscience #78078

**Storage Conditions**

Cells will arrive upon 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](mailto:support@bpsbioscience.com) if the cells are not frozen in dry ice upon arrival.

**License Disclosure**

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

**Troubleshooting Guide**

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

**Media Formulations**

For best results, it is *highly recommended* to use these validated and optimized media from BPS Bioscience. To formulate a comparable but not BPS validated media, formulation components can be found below.



Note: Thaw Media does *not* contain selective antibiotics. However, Growth Media *does* contain selective antibiotics, which are used for maintaining cell lines over many passages. Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 2K.

**Media Required for Cell Line Culture****Thaw Medium 2 (BPS Bioscience #60184):**

RPMI 1640 medium (Thermo Fisher, #A1049101) supplemented with 10% FBS (Thermo Fisher, #26140079), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01).

**Growth Medium 2K (BPS Bioscience #78078):**

Thaw Medium 2 (BPS Bioscience, #60184) plus 0.25 µg/ml of Puromycin (Invivogen, #ant-pr-1) to ensure recombinant expression.

Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 2K to ensure recombinant expression is maintained.

**Recommended Culture Conditions****Frozen cells:**

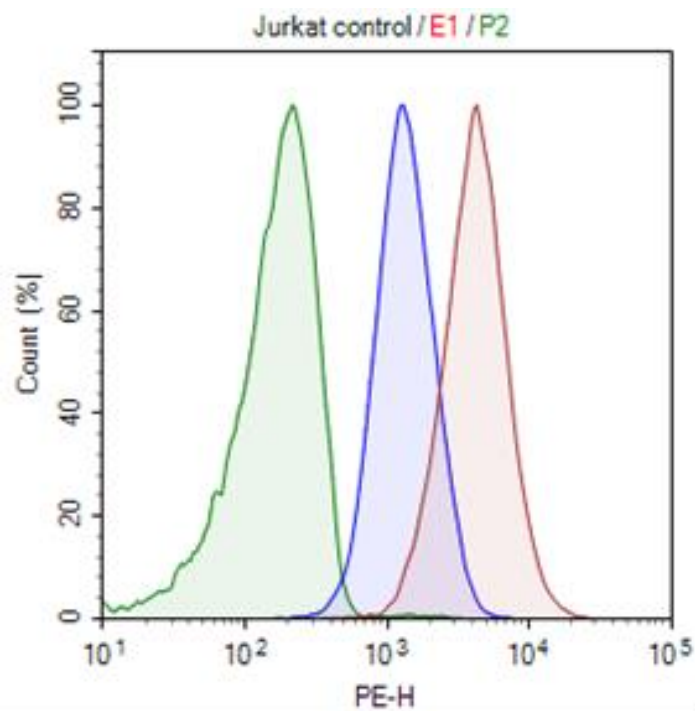
1. It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, and then transfer the entire contents of the vial to a tube containing 10 ml of pre-warmed Thaw Medium 2 (**no Puromycin**).
2. Then spin the cells down, remove the supernatant, and resuspend the cells in 5 ml of pre-warmed Thaw Medium 2 (**no Puromycin**).
3. Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 24 hours of culture, add an additional 3-4 ml of Thaw Medium 2 (**no Puromycin**).
5. At first passage, switch to Growth Medium 2K (contains Puromycin).
6. Cells should be split before they reach 2 x 10<sup>6</sup> cells/ml.

**Cryopreservation:**

1. When cells have reached 90% confluency, spin cells and remove medium from the pellet.
2. Resuspend the cells in freezing medium (10% DMSO in FBS).
3. Freeze cells using a reduced rate freezing box (-0.5°C to -1°C per minute) down to -80°C, then move cells to liquid nitrogen for long term storage.



Cells have been demonstrated to be stable for at least 15 passages; BPS recommends preparing frozen stocks so cells are not used beyond passage 20.

**Validation Data**

*Figure 1. Flow cytometry analysis of intracellular expression of Cas9 in Jurkat cells.*

*Cells were stained with PE-labeled anti-FLAG antibody (BioLegend, #637309) and analyzed by flow cytometry. The parental Jurkat cells are shown in green, the Cas9-expressing Jurkat High expression cell line (BPS Bioscience, #78136-H) is shown in red, and the Cas9-expressing Jurkat Low expression cell line (BPS Bioscience, #78136-L) is shown in blue.*

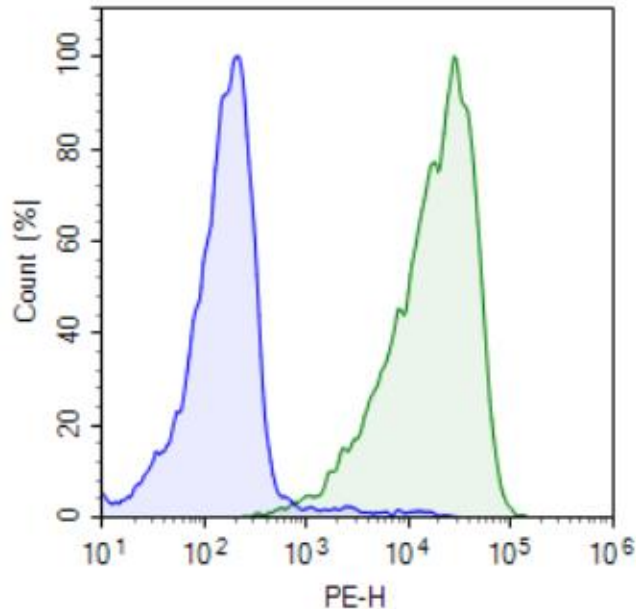


Figure 2. Flow cytometry analysis of TCR expression following knock-out in Cas9-expressing Jurkat cells. Cas9-expressing cells (BPS Bioscience, #78136-H) were electroporated with 0.1 nmol TCR sgRNA, and TCR expression was analyzed by flow cytometry 72 hours later. The parental Cas9-expressing Jurkat High expression cell line (BPS Bioscience, #78136-H) is shown in green, and the TCR knock-out cell pool is shown in blue.

### Vector and Sequence

*Streptococcus pyogenes* Cas9, including a C-terminal FLAG tag, was transduced via lentivirus (BPS Bioscience, #78066).

MDKKYSIGLDIGTNSVGVAVITDEYKVPSSKFKVLGNTDRHSIKKNLIGALLFDSGETAEATRLKRTARRRYTRRKNRICYLQEIFSN  
 EMAKVDDSFHRLSEESFLVEEDKKHERHPIFGNIVDEVAYHEKYPTIYHLRKKLVDSTDKADLRILIYALAHMIKFRGHFLIEGDLNP  
 DNSDVKLFIQLVQTYNQLFEENPINASGVDAKAILSARLSKSRLENLIAQLPGEKKNLFGNLIASLGLTPNFKSNFDLAEDAKL  
 QLSKDYDDDLNLLAQIGDQYADLFLAAKNLSDAILLSDLRVNTEITKAPLSASMIKRYDEHHQDLTLLKALVRQQLPKEYKEIFF  
 DQSKNGYAGYIDGGASQEEFYKFIKPILEKMDGTEELLVKNLREDLLRKQRTFDNGSIPHQIHLGELHAILRRQEDFYFPLKDNREKI  
 EKILTFRIPYYVGPLARGNSRFAWMTRKSEETITPWNFEVVDKGASAQSFIERMTNFDKNLPNEKVLPKHSLLYEYFTVYNELTKV  
 KYVTEGMRKPAFLSGEQKKAIVDLLFKTNRKVTVKQLKEDYFKKIECFDSVEISGVEDRFNASLGTYHDLLKIIKDKDFLDNEENEDIL  
 EDIVLTLTLFEDREMIEERLKYAHLFDDKVMKQLKRRRYTGWGRLSRKLINGIRDKQSGKTILDFLKSDFANRNFMLIHDDSLT  
 FKEDIQKAQVSGQGDSLHEHIANLAGSPAIKKILQTVKVVDELVKVMGRHKPENIVIEMARENQTTQKGQKNSRERMKRIEEGI  
 KELGSQILKEHPVENTQLQNEKLYLYLQNGRDMYVDQELDINRLSDYDVDHIVPQSFLKDDSIDNKVLRSDKNRKGSDNVPSEE  
 VVKMKNYWRQLLNAKLITQRKFDNLTKAERGGSELDKAGFIKRLVETRQITKHVAQILDSRMNTKYDENDKLIREVKVITLKS  
 KLVSDFRKDFQFYKREINNYHHAHDAYLNAVVGTAIIKKYKPLESEFVYGDYKVVYDVRKMIKAKSEQEIGKATAKYFFYSNIMNFFK  
 TEITLANGEIRKPLIETNGETGEIVWDKGRDFATVRKVLSPQVNIKKTEVQTGGFSKESILPKRNSDKLIARKKDWDPKYYGGF  
 DSPTVAYSVLVAKVEKGSKLLKSVKELLGITIMERSSEFKNPIDFLEAKGYKEVKKDLIILPKYSLELENGRKRMLASAGELQKG  
 NELALPSKYVNFYLAHYEKLKGPEDNEQKQLFVEQHKHYLDEIIIEQISEFSKRVLADANLDKVL SAYNKHRDKPIREQAENIIHL  
 FTLNLGAPAAFKYFDTTIDRKRYTSTKEVLDTLIHQISITGLYETRIDLSQLGGDKRPAATKKAGQAKKKKDYKDDDDK

**Related Products**

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
Cas9-Expressing Jurkat cell pool	<a href="#">78070</a>	2 vials
Cas9 Lentivirus (puromycin selection)	<a href="#">78066</a>	500 µl x 2
Cas9, His-tag ( <i>S. pyogenes</i> )	<a href="#">100206-1</a>	50 µg
Thaw Medium 2	<a href="#">60184-1</a>	100 ml
Growth Medium 2K	<a href="#">78078</a>	500 ml

**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.*