

Description

This cell line is a clonal derivative from the Cas9-Expressing Neuro2A Cell Pool ([BPS Bioscience, #78087](#)). 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 low-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 Neuro2A 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 Neuro2A cells.
2. Implementing sgRNA screens in Cas9 expressing Neuro2A cells.

Materials Provided

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

Host Cell

Neuro2A is a mouse neuroblastoma cell line. Adherent neuroblast cells.

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 1	BPS Bioscience #60187
Growth Medium 1Q	BPS Bioscience #78096

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 if the cells are not frozen in dry ice upon arrival.

License Disclosure

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

Troubleshooting Guide

For all questions, please email 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% CO₂ using Growth Medium 1Q.

Media Required for Cell Line Culture

Thaw Medium 1 (BPS Bioscience #60187):

MEM medium (Gibco, #11095-080) supplemented with 10% FBS (Gibco, #26140-079), 1X MEM Non-essential Amino Acids (Corning, #25-025-Cl), 1% Sodium Pyruvate (Corning, #25-000-Cl), and 1% Penicillin/Streptomycin (Gibco, #15140-122).

Growth Medium 1Q (BPS Bioscience #78096):

Thaw Medium 1 (BPS Bioscience, #60187) plus 1 µg/ml Puromycin (Invivogen, #ant-pr-1) to ensure recombinant expression.

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 1Q to ensure recombinant expression is maintained.

Recommended Culture Conditions

Frozen Cells:

1. Prepare a 15 mL conical tube with 10 ml of pre-warmed Thaw Medium 1 (**no Puromycin**).
2. Quickly thaw cells in a 37°C water bath with constant and slow agitation.
3. Clean the outside of the vial with 70% ethanol and immediately transfer the entire contents to the conical tube. Spin cells at 200 x g for 5 minutes, and remove all medium from the pellet.
4. Resuspend in 15 ml Thaw Medium 1 (**no Puromycin**) and transfer to a T-75 flask.
5. Gently rock the flask to distribute the cells. Incubate the cells in a humidified 37°C incubator with 5% CO₂.
6. 24 hours after incubation, change culture to fresh Thaw Medium 1 (**no Puromycin**); avoid disturbing the attached cells.
7. Continue to monitor growth for 2-3 days and change the media to remove dead cell debris, if necessary.
8. Begin adding Growth Medium 1Q (contains Puromycin) after multiple cell colonies (in clumps) start to appear (indicative of healthy cell division).

Subculture:

1. When cells have reached 90% confluency, remove Growth Medium 1Q and gently wash cells twice with PBS (without Magnesium or Calcium).
2. Treat cells with 2 ml of 0.25% Trypsin/EDTA and incubate for 2-3 minutes at 37°C.
3. Dispense 10 ml of pre-warmed Growth Medium 1Q to the trypsinized cells and gently pipette up and down to neutralize the trypsin and break apart any cell clumps.
4. Transfer cells to a conical tube and centrifuge at 200 x g for 5 minutes.
5. Remove the supernatant and re-suspend the cell pellet in 10-14 ml of prewarmed Growth Medium 1Q.
6. Dispense 2 ml of cell suspension into a new T-75 flask containing prewarmed 15 ml of Growth Medium 1Q.
7. Incubate cells in a humidified 37°C incubator with 5% CO₂.

Cryopreservation:

1. When cells have reached 90% confluency, use trypsin to remove cells from plate as above, 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

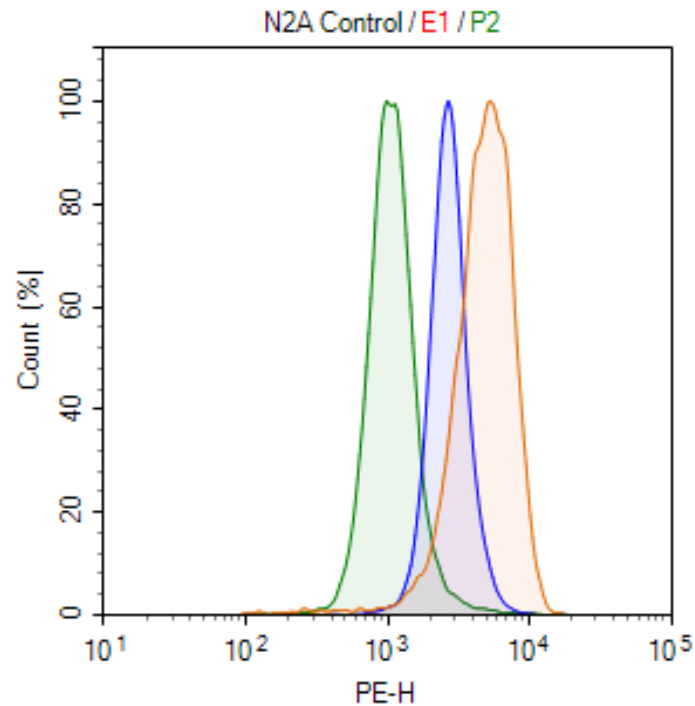


Figure 1. Flow cytometry analysis of intracellular expression of Cas9 in Neuro2A cells. Cells were stained with PE-labeled anti-FLAG antibody (BioLegend, #637309) and analyzed by flow cytometry. The parental Neuro2A cells are shown in green, the Cas9-expressing Neuro2A High expression cell line (BPS Bioscience, #78137-H) is shown in orange, and the Cas9-expressing Neuro2A Low expression cell line (BPS Bioscience, #78137-L) is shown in blue.

Vector and Sequence

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

MDKKYSIGLDIGTNSVGVAVITDEYKVPSPKFKVLGNTDRHSIKKNLIGALLFDSGETAEATRLKRTARRRYTRRKNRICYLQEIFSN
 EMAKVDDSFHRLVESFLVEEDKKHERHPIFGNIVDEVAYHEKYPTIYHLRKKLVDSTDKADLRILIYALAHMIKFRGHFLIEGDLNP
 DNSDVKLFIQLVQTYNQLFEENPINASGVDAKAILSARLSKSRLENLIAQLPGEKKNLFGNLIASLGLTPNFKSNFDLAEDAKL
 QLSKDYDDDLNLLAQIGDQYADLFLLAAKNLSDAILLSDILRVNTEITKAPLSASMIKRYDEHHQDLTLLKALVRQQLPKEYKEIFF
 DQSKNGYAGYIDGGASQEEFYKFIKPILEKMDGTEELLVKNLREDLLRKQRTFDNGSIPHQIHLGELHAILRRQEDFYFPLKDNREKI
 EKILTFRIPYYVGPLARGNSRFAWMTRKSEETITPWNFEVVDKGASAQSFIERMTNFDKNLPNEKVLPKHSLLYEYFTVYNELTKV
 KYVTEGMRKPAFLSGEQKKAIVDLLFKTNRKVTVKQLKEDYFKKIECFDSVEISGVEDRFNASLGTYHDLKIIKDKDFLDNEENEDIL
 EDIVLTLTLFEDREMIEERLKTYAHLFDDKVMKQLKRRRYTGWGRLSRKLINGIRDKQSGKTILDFLKSDGFANRNFQMQLIHDDSLT
 FKEDIQKAQVSGQGDSLHEHIANLAGSPAIKKILQTVKVVDELVKVMGRHKPENIVIEMARENQTTQKGQKNSRERMKRIEEGI
 KELGSQILKEHPVENTQLQNEKLYLYLQNGRDMYVDQELDINRLSDYDVDHIVPQSFLKDDSIDNKVLRSDKNRKGSDNVPSEE
 VVKMKNYWRQLLNAKLITQRKFDNLTKAERGGSELDKAGFIKRLVETRQITKHVAQILDSRMNTKYDENDKLIREVKVITLKS
 KLVSDFRKDFQFYKREINNYHHAHDAYLNAVVGTAIIKKYKPLESEFVYGDYKVVYDVRKMIKAKSEQEIGKATAKYFFYSNIMNFFK
 TEITLANGEIRKRPDIETNGETGEIVWDKGRDFATVRKVLSPQVNIKKTEVQTGGFSKESILPKRNSDKLIARKKDWDPKYYGGF
 DSPTVAYSVLVVAKEKGSKLLKSVKELLGITIMERSSEFNPIDFLEAKGYKEVKDLIIPKYSLELENGRKRMLASAGELQKG
 NELALPSKYVNFYLAHYEKLKSPEDNEQKQLFVEQHKHYLDEIIEQISEFSKRVLADANLDKVL SAYNKHRDKPIREQAENIIHL
 FTLTNLGAPAAFKYFDTTIDRKRYTSTKEVLDTLIHQISITGLYETRIDLSQLGGDKRPAATKKAGQAKKKKDYKDDDDK

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
Cas9-Expressing Neuro2a Cell line – High Expression	78137-H	2 vials
Cas9-Expressing Neuro2a cell pool	78087	2 vials
Cas9 Lentivirus (puromycin selection)	78066	500 µl x 2
Cas9, His-tag (<i>S. pyogenes</i>)	100206-1	50 µg
Thaw Medium 1	60187-1	100 ml
Growth Medium 1Q	78096	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.