

# Transfection of neural stem cells with Lipofectamine Stem Transfection Reagent in StemPro medium

## NSC media, passaging reagents, and complexation medium

Component	Cat. No.
StemPro NSC SFM	A1050901
CTS GlutaMAX-I Supplement	A1286001
Geltrex LDEV-Free, hESC-Qualified, Reduced Growth Factor Basement Membrane Matrix	A1413302
Dulbecco's Phosphate Buffered Saline (DPBS) without calcium and magnesium	14190144
StemPro Accutase Cell Dissociation Reagent	A1110501
Opti-MEM I Reduced Serum Medium	31985062

Starting with undifferentiated human primary neural stem cells (NSCs) or pluripotent stem cell (PSC)-derived NSCs, expanded in a defined culture system such as Gibco™ StemPro™ NSC Serum-Free Medium (SFM) on Gibco™ Geltrex™ matrix, is ideal for efficient transfection.

### Passaging

- Maintain NSCs in the format of your choice, such as 6-well plates, 60 cm dishes, T-25 flasks, or T-75 flasks coated with Geltrex matrix, in StemPro SFM. Propagating NSCs in T-25 flasks and transfecting in 24-well plates are convenient formats used in this protocol.
- Passage NSCs every 3 to 5 days at 90–100% confluence.
- Use Gibco™ StemPro™ Accutase™ Cell Dissociation Reagent to generate a single-cell suspension of NSCs for both expansion and seeding for transfection.

### Precoating 24-well plates with Geltrex matrix for transfection

1. Prepare a 1:100 dilution of Geltrex matrix in cold Gibco™ DMEM/F-12 with Gibco™ GlutaMAX™ Supplement (Cat. No. 10565).
  2. Add 300 µL of diluted Geltrex matrix to each well of a 24-well plate and incubate at 37°C for ≥1 hour, before use.
- **Tip:** Geltrex matrix-coated plates can be prepared ahead of time and stored for up to 2 weeks at 4°C. Equilibrate at room temperature for 1 hour before plating cells.

### Seeding cells for transfections

1. When NSC cultures are ~90–100% confluent, remove the StemPro NSC SFM.
2. Wash NSCs once with 10 mL of DPBS without calcium and magnesium; aspirate the medium and discard.
3. Add 1 mL of room-temperature StemPro Accutase reagent to each T-25 flask, swirl to evenly coat the NSCs, and incubate for 2–5 minutes at room temperature.
  - **Important:** To maximize transfection efficiency, seeding a single-cell suspension of NSCs prepared with StemPro Accutase reagent is recommended.
4. Observe cells on an inverted microscope to confirm that NSCs are detached; firmly tap the flask to aid in the detachment of NSCs, as necessary.
5. Add 9 mL of StemPro NSC SFM to inactivate the StemPro Accutase reagent.
6. Gently triturate and rinse the flask to generate a single-cell suspension, and transfer the cell suspension into a 15 mL conical tube.
7. Centrifuge the NSC cell suspension at 200 x g for 4 minutes.
8. Aspirate the supernatant and resuspend the pellet to a single-cell suspension in 3 mL of StemPro NSC SFM.
9. Perform a total viable cell count with the Invitrogen™ Countess™ II Automated Cell Counter or another method.
10. Dilute with additional StemPro NSC SFM to a final concentration of 150,000 cells/mL.
11. Aspirate the Geltrex matrix from the wells of a precoated 24-well plate.
  - **Important:** Proliferating NSC cultures need room to expand during transfection, so plate the recommended starting number of cells (step 12) to achieve 30–60% confluence on the day of transfection.
12. Add 0.5 mL of the NSC suspension in StemPro NSC SFM to plate 75,000 cells/well in the precoated 24-well plate.
13. Return the plate to the incubator and culture at 37°C with 5% CO<sub>2</sub>, overnight.
  - **Important:** You do not need to change the medium on the day of transfection.

## DNA transfection protocol

Perform the following steps, which have been optimized for using Invitrogen™ Lipofectamine™ Stem Transfection Reagent with NSCs:

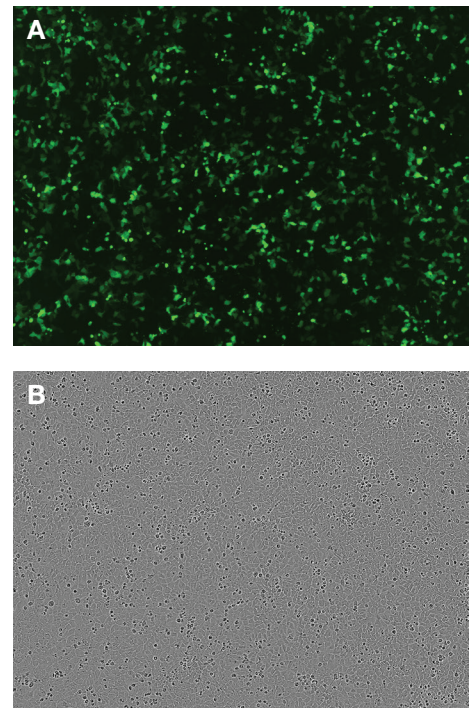
Step	Tube	Complexation component	Amount per well (24-well plate)
1	Tube 1	Opti-MEM I medium	25 $\mu$ L
		Lipofectamine Stem reagent	1 $\mu$ L
2	Tube 2	Opti-MEM I medium	25 $\mu$ L
		DNA (0.5–5 $\mu$ g/ $\mu$ L)	500 ng
3	Add tube 2 solution to tube 1, and mix well.		
4	Incubate mixture from step 3 for 10 minutes at room temperature.		
5	Add 50 $\mu$ L of complex from step 4 to each well; gently swirl plate to ensure even distribution of the complex across the entire well.		
6	Return culture dish to incubator and culture at 37°C with 5% CO <sub>2</sub> , overnight.		
7	The following day, overlay an additional 0.5 mL of StemPro NSC SFM per well, if NSCs are going to be transfected for 48 hours.		

## Analysis of transfection efficiency

Observe NSCs transfected with a GFP reporter construct at 24 and 48 hours posttransfection by fluorescence microscopy or flow cytometry for endpoint analysis (Figure 1).

## Tips and tricks

- The amount of Lipofectamine Stem reagent required for optimal transfection depends on the amount of NSCs plated and the amount of DNA used.
- If cytotoxicity from the DNA preparation is evident, reducing the amount of DNA to 250 ng per well can improve survival while maintaining efficient transfection.
- Using a plasmid with a promoter that is active in human NSCs, such as the EF1 $\alpha$  promoter, is critical for assessing transfection efficiency; some promoters such as the cytomegalovirus (CMV) promoter can be transcriptionally silenced in NSCs.



**Figure 1. Posttransfection analysis of NSCs.** (A) Fluorescence image demonstrating 59% transfection efficiency, and (B) bright-field image. NSCs are shown 24 hours after transfection with 500 ng of a 6 kb EF1 $\alpha$ -GFP plasmid and 1  $\mu$ L of Lipofectamine Stem reagent in StemPro NSC SFM on Geltrex matrix.

## mRNA transfection protocol

Perform the following steps, which have been optimized for using Lipofectamine Stem reagent with NSCs:

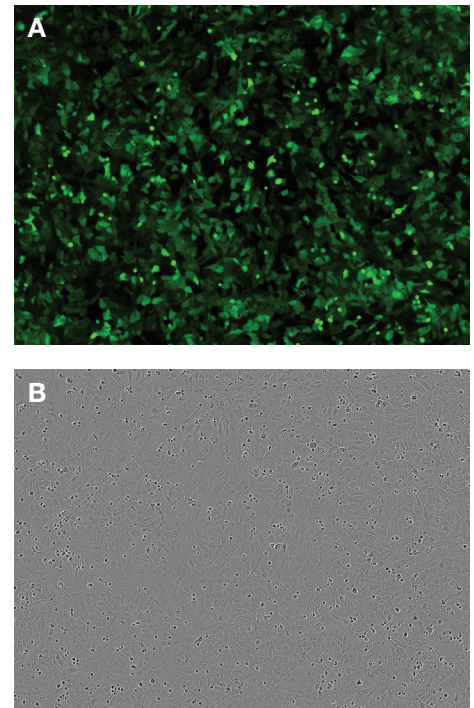
Step	Tube	Complexation component	Amount per well (24-well plate)
1	Tube 1	Opti-MEM I medium	25 $\mu$ L
		Lipofectamine Stem reagent	1 $\mu$ L
2	Tube 2	Opti-MEM I medium	25 $\mu$ L
		mRNA (0.5–5 $\mu$ g/ $\mu$ L)	250 ng
3	Add tube 2 solution to tube 1, and mix well.		
4	Incubate mixture from step 3 for 10 minutes at room temperature.		
5	Add 50 $\mu$ L of complex from step 4 to each well; gently swirl plate to ensure even distribution of the complex across the entire well.		
6	Return culture dish to incubator and culture at 37°C with 5% CO <sub>2</sub> , overnight.		
7	The following day, overlay an additional 0.5 mL of StemPro NSC SFM per well, if NSCs are going to be transfected for 48 hours.		

### Analysis of transfection efficiency

Observe PSCs transfected with a fluorescent mRNA at 24 and 48 hours posttransfection by fluorescence microscopy or flow cytometry for endpoint analysis (Figure 2).

### Tips and tricks

- The amount of mRNA required to generate a specific biological readout will vary by user application; Lipofectamine Stem reagent efficiently delivers mRNA into NSCs across a range of dosages.
- Including an independent GFP mRNA (50 ng) in addition to your transcript of interest allows an independent assessment of transfection efficiency.
- If cytotoxicity from the mRNA preparation is evident, reducing the amount of mRNA to 125 ng per well can improve survival while maintaining efficient transfection.
- The method of generation and purification of *in vitro* transcribed (IVT) mRNA can contribute to toxicity as well as translational repression.
  - An anti-reverse cap analog (ARCA) system, included in the Invitrogen™ mMACHINE™ Kit for *in vitro* transcription, and Invitrogen™ MEGAclear™ columns can be used to eliminate uncapped transcripts and small unincorporated nucleotides that can contribute to cytotoxicity.



**Figure 2. Posttransfection analysis of NSCs.** (A) Fluorescence image demonstrating 70% transfection efficiency, and (B) bright-field image. iPSC-derived NSCs (NCRM1) are shown 36 hours after transfection with 250 ng of GFP mRNA and 1  $\mu$ L of Lipofectamine Stem reagent in StemPro NSC SFM on Geltrex matrix.

## Ribonucleoprotein (RNP) transfection protocol

RNP complex components:

- Invitrogen™ GeneArt™ Platinum™ Cas9 Nuclease (Cat. No. B25641)
- gRNA (see “Designing and generating gRNA by *in vitro* transcription”)

Perform the following steps, which have been optimized for using Lipofectamine Stem reagent with NSCs:

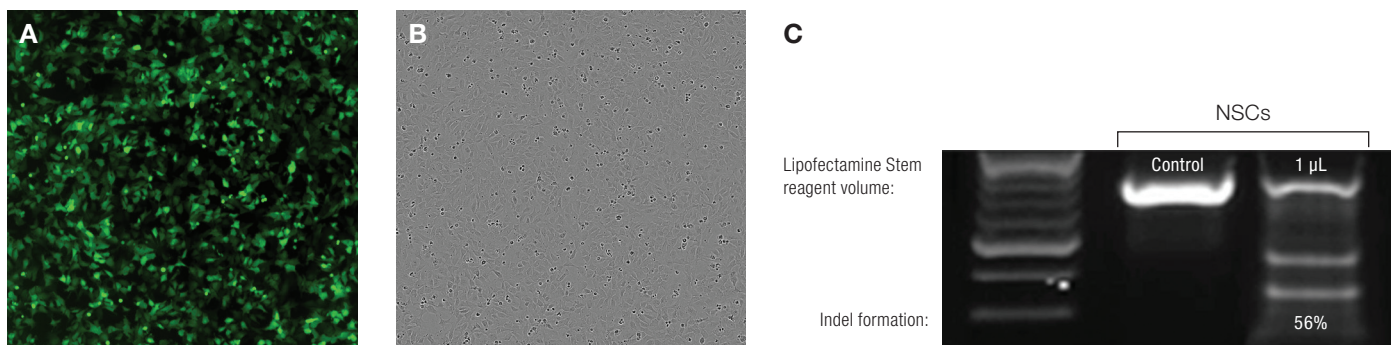
Step	Tube	Complexation component	Amount per well (24-well plate)
1	Tube 1	Opti-MEM I medium	25 $\mu$ L
		Lipofectamine Stem Reagent	1 $\mu$ L
2	Tube 2	Opti-MEM I medium	25 $\mu$ L
		Cas9 nuclease	500 ng
		gRNA (0.1–0.5 $\mu$ g/ $\mu$ L)	125 ng
3	Add tube 2 solution to tube 1, and mix well.		
4	Incubate mixture from step 3 for 10 minutes at room temperature.		
5	Add 50 $\mu$ L of complex from step 4 to each well; gently swirl plate to ensure even distribution of the complex across the entire well.		
6	Return culture dish to incubator and culture at 37°C with 5% CO <sub>2</sub> , overnight.		
7	The following day, overlay an additional 0.5 mL of StemPro NSC SFM per well, if NSCs are going to be transfected for 48 hours.		

## Analysis of transfection efficiency

Observe PSCs transfected with a GFP reporter construct at 24 and 48 hours posttransfection by fluorescence microscopy or flow cytometry, and analyze double-stranded break (DSB) formation using the Invitrogen™ GeneArt™ Genomic Cleavage Detection Kit or a similar assay (Figure 3).

## Tips and tricks

- Adding 50 ng of GFP mRNA to the transfection complex along with the RNP complex can provide an independent measure of transfection efficiency.



**Figure 3. Posttransfection analysis of NSCs.** (A) Fluorescence image demonstrating 60% transfection efficiency, and (B) bright-field image. iPSC-derived NSCs (NCRM1) are shown 24 hours posttransfection with 500 ng of GeneArt Platinum Cas9 Nuclease, 125 ng of gRNA, 50 ng of GFP mRNA, and 1  $\mu$ L of Lipofectamine Stem reagent in StemPro NSC SFM on Geltrex matrix. (C) Genomic cleavage detection analysis of iPSC-derived NSCs 48 hours posttransfection, demonstrating 56% indel formation within the *EMX1* locus.



### **Designing and generating gRNA by *in vitro* transcription**

In addition to RNP transfection efficiency, the efficiency of DSB/indel formation at a given locus can depend on gRNA design. Use the Invitrogen™ GeneArt™ CRISPR Search and Design Tool, available at [thermofisher.com/crisprdesign](https://thermofisher.com/crisprdesign), to search our database of >600,000 predesigned gRNA sequences specific to every gene in the human genome. These predesigned gRNAs are optimized for gene knockout and typically target the first 3 transcribed exons per gene.

Clone and generate your own gRNA using the Invitrogen™ GeneArt™ Precision gRNA Synthesis Kit (Cat. No. A29377). gRNA concentration can be quantified on the Invitrogen™ Qubit™ 3 Fluorometer (Cat. No. Q33216) coupled with the Invitrogen™ Qubit™ RNA BR Assay Kit (Cat. No. Q10210).

Find out more at [thermofisher.com/transfection](https://thermofisher.com/transfection)

**ThermoFisher**  
S C I E N T I F I C