Lipofectamine™ Stem Transfection Reagent

USER GUIDE

Pub. No. MAN0017080 Rev. A.0



Contents

Catalog Number Size • STEM00001 $0.1 \, \mathrm{mL}$ STEM00003 $0.3 \, \mathrm{mL}$ STEM00008 $0.75 \, \text{mL}$ STEM00015 1.5 mL



Storage

Store at 4°C (do not freeze).



- Opti-MEM™ I Reduced Serum Medium
- Microcentrifuge tubes



Timing

Preparation: 10 minutes Incubation: 10 minutes Final incubation: 1–2 days



Product description

- Lipofectamine[™] Stem Reagent is a proprietary formulation for transfecting nucleic acids into a wide range of human stem cells.
- Lipofectamine[™] Stem Reagent should be made in Opti-MEM™ I Reduced Serum Medium and can be added directly to cells in culture growing with or without serum.
- It is not necessary to remove transfection complexes. Additional medium may be added the following day to stem cell cultures without interfering with the transfection.



Important quidelines

- The amount of Lipofectamine™ Stem Reagent required for optimal transfection depends upon the amounts of stem cells plated and DNA used (See recommended amounts).
 - □ Proliferating stem cell culture need room to expand, so plate the recommended starting cell numbers to achieve 30-60% confluency on the day of transfection.
 - □ Using lower amounts of DNA, mRNA, or ribonucleoprotein (RNP) complexes can yield improved expression, while minimizing risks of cytotoxicity from excess foreign constructs.



Online resources

Visit our product page for additional information and cell specific protocols. For support, visit thermofisher.com/support.

Protocol outline

- A. Plate cells so they will be 30–60% confluent at the time of transfection.
- B. Prepare transfection complexes.
- C. Add transfection complexes to cells.

Cell seeding recommendations for various stem cell models

Stem Cell Type	96-well	24-well	6-well
Human Pluripotent	$1-1.5 \times 10^4 \text{ cells}$	$5-7.5 \times 10^4 \text{ cells}$	$2.5-3.75 \times 10^5$ cells
Human Neural	1.5×10^4 cells	7.5×10^4 cells	3.75×10^5 cells
Human Mesenchymal	0.5×10^4 cells	2.5×10^4 cells	1.25×10^5 cells

DNA transfection

See page 2 to view a typical DNA transfection procedure.

Component	96-well	24-well	6-well
Final DNA per well	50–100 ng	250–500 ng	1.25–2.5 µg
Final Lipofectamine™ Stem Reagent per well	0.2–0.4 μL	1.0-2.0 μL	5.0–10 μL

mRNA Transfection

mRNA can be transfected in a 24-well plate with 1 or 2 µL Lipofectamine™ Stem Reagent complexed with 250 ng to 500 ng of mRNA per well.

Ribonucleoprotein (RNP) Transfection

RNP complexes can be transfected into stem cells in a 24-well plate by combining 0.5 to 1.5 µg of GeneArt™ Platinum™ Cas9 nuclease with 125 to 375 ng of gRNA and then adding 1 to 2 µL of Lipofectamine Stem[™] Transfection Reagent per reaction for each well.

Limited Product Warranty and Disclaimer Details



Lipofectamine™Stem Transfection Reagent Protocol:

To optimize transfection of your stem cells in your culture conditions, we recommend starting with two amounts of DNA and two amounts of Lipofectamine™ Stem Transfection Reagent following the chart below. Volumes are listed on a per well basis in a 24-well format. Adjust the final amounts of components according to your experimental design and cell culture format. For additional information on setting up mRNA or RNP transfections see thermofisher.com/lipofectaminestem.

Timeline		neline	Step	Procedure details					
Day 0	1	 	Seed cells to be 30-60% confluent at transfection	Adherent stem cells: 2.5–7.5 x 10 ⁴ per well in 4 wells of a 24-well plate					
Day 1	2			Well number					
			Dilute 2 amounts of Lipofectamine™ Stem Reagent in Opti-MEM™ I Medium	Component	1	2	3	4	
	_	UUUU		Opti-MEM™ I Medium	25 µL	25 µL	25 μL	25 µL	
			Lipofectamine™ Stem Reagent	1 μL	1 μL	2 μL	2 μL		
	3	Dilute 2 amounts of DNA in Opti-MEM™ I Medium	Opti-MEM™ I Medium	25 µL	25 µL	25 μL	25 µL		
			DNA	250 ng	500 ng	250 ng	500 ng		
			Add diluted DNA to diluted	Diluted DNA volume	25 µL	25 µL	25 μL	25 µL	
	4	Lipofectamine™ Stem Reagent (1:1 ratio)	Diluted Lipofectamine™ Stem Reagent volume	25 μL	25 μL	25 μL	25 μL		
	5	10	Incubate	Incubate for 10 minutes at room temperature.					
	6			Well number					
				Component	1	2	3	4	
		Add DNA-lipid complex to cells	DNA-lipid complex per well	50 μL	50 μL	50 μL	50 µL		
				Final DNA used per well	250 ng	500 ng	250 ng	500 ng	
			Final Lipofectamine™ Stem Reagent used per well	1 μL	1 µL	2 μL	2 µL		
Day 2-3	7		Visualize/analyze transfected cells	Incubate and monitor transfected stem cells at 37°C for 1–2 days.					