

Lipofectamine® Reagent

	Package Contents	Catalog Number	Size
		<ul style="list-style-type: none"> 18324-010 18324-012 18324-020 	<ul style="list-style-type: none"> 0.3 mL 1.0 mL 4 × 1 mL
	Storage Conditions	Store at 4°C (do not freeze).	
	Required Materials	<ul style="list-style-type: none"> Plasmid DNA (0.5–5 µg/µL stock) Opti-MEM® Reduced Serum Medium Eppendorf tubes 	
	Timing	Preparation: 10 minutes Incubation: 5 minutes Final Incubation: 1–3 days	
	Selection Guide	Lipofectamine® Reagents Go online to view related products.	
	Product Description	<ul style="list-style-type: none"> Lipofectamine® Reagent is a proprietary formulation for transfecting nucleic acids into a wide range of eukaryotic cells. 	
	Important Guidelines	<ul style="list-style-type: none"> DNA-Lipofectamine® complexes must be made in serum-free medium such as Opti-MEM® Reduced Serum Medium and can be added directly to cells in culture medium, in the presence or absence of serum/antibiotic. It is not necessary to remove complexes or change/add medium after transfection. The amount of Lipofectamine® Reagent required for successful transfection varies depending on the cell type and passage number. Start any new transfection by testing the recommended four concentrations of Lipofectamine® Reagent to determine an optimum amount. 	
	Online Resources	Visit our product page for additional information and protocols. For support, visit www.lifetechnologies.com/support .	

For Research Use Only. Not for use in diagnostic procedures.



Protocol Outline

- Plate cells so they will be 70–90% confluent at the time of transfection.
- Prepare plasmid DNA-lipid complexes.
- Add DNA-lipid complexes to cells.

Lipofectamine® DNA Transfection Reagent Protocol

i See page 2 to view a typical plasmid transfection procedure.

Transfection Amounts

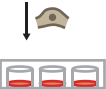


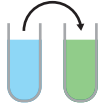

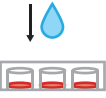

Component	96-well	24-well	6-well
DNA per well	100 ng	500 ng	2500 ng
PLUS™ Reagent (Optional)	0.1 µL	0.5 µL	2.5 µL
Lipofectamine® Reagent per well	0.2–0.5 µL	1–2.5 µL	5–12.5 µL

i **Scaling Up or Down Transfections**

i **Limited Product Warranty and Disclaimer Details**

Lipofectamine® DNA Transfection Reagent Protocol

Transfect cells according to the following table. Use the indicated volume of DNA and PLUS™ Reagent with each of the four volumes of Lipofectamine® Reagent. **Each reaction mix is sufficient for triplicate (96-well), duplicate (24-well), and single well (6-well) transfections, and accounts for pipetting variations.** For even less toxicity, reduce the amount of DNA-lipid complex to the cells, or reduce the amount of DNA used to make complexes.

Timeline		Steps	Procedure Details				
Day 0	1		Seed cells to be 70-90% confluent at transfection	Component	96-well	24-well	6-well
	2		Dilute Lipofectamine® Reagent in Opti-MEM® Medium	Adherent cells	1–4 × 10 ⁴	0.5–2 × 10 ⁵	0.25–1 × 10 ⁶
Day 1	3		Dilute DNA in Opti-MEM® Medium, then add PLUS™ Reagent	Opti-MEM® Medium	25 µL × 4	50 µL × 4	150 µL × 4
	4		Add diluted DNA to diluted Lipofectamine® Reagent (1:1 ratio)	Lipofectamine® Reagent	1, 1.5, 2, 2.5 µL	2, 3, 4, 5 µL	6, 9, 12, 15 µL
	5		Incubate	Opti-MEM® Medium	125 µL	250 µL	700 µL
	6		Add DNA-lipid complex to cells	DNA (0.5–5 µg/µL)	2.5 µg	5 µg	14 µg
Day 2–4	7		Visualize/analyze transfected cells	PLUS™ Reagent (Optional)	2.5 µL	5 µL	14 µL
				Diluted DNA (with PLUS™ Reagent) Total	25 µL	50 µL	150 µL
				Diluted Lipofectamine® Reagent	25 µL	50 µL	150 µL
				Incubate for 5 minutes at room temperature.			
				Component	96-well	24-well	6-well
				DNA-lipid complex per well	10 µL	50 µL	250 µL
				Final DNA used per well	100 ng	500 ng	2500 ng
				Final Lipofectamine® Reagent used per well	0.2–0.5 µL	1.0–2.5 µL	5.0–12.5 µL
				Incubate cells for 1–3 days at 37°C. Then, analyze transfected cells.			