









# Lipofectamine® RNAiMAX Reagent

 <b>Package Contents</b>	<table border="1"> <thead> <tr> <th>Catalog Number</th> <th>Size</th> </tr> </thead> <tbody> <tr> <td>13778-100</td> <td>0.1 mL</td> </tr> <tr> <td>13778-030</td> <td>0.3 mL</td> </tr> <tr> <td>13778-075</td> <td>0.75 mL</td> </tr> <tr> <td>13778-150</td> <td>1.5 mL</td> </tr> <tr> <td>13778-500</td> <td>15 mL</td> </tr> </tbody> </table>	Catalog Number	Size	13778-100	0.1 mL	13778-030	0.3 mL	13778-075	0.75 mL	13778-150	1.5 mL	13778-500	15 mL
Catalog Number	Size												
13778-100	0.1 mL												
13778-030	0.3 mL												
13778-075	0.75 mL												
13778-150	1.5 mL												
13778-500	15 mL												
 <b>Storage Conditions</b>	Store at 4°C (do not freeze).												
 <b>Required Materials</b>	<ul style="list-style-type: none"> <li>siRNA or miRNA (10 µM stock)</li> <li>Opti-MEM® Reduced Serum Medium</li> <li>Eppendorf tubes</li> </ul>												
 <b>Timing</b>	Preparation: 10 minutes Incubation: 5 minutes Final Incubation: 1-3 days												
 <b>Selection Guide</b>	<a href="#">Lipofectamine® Reagents</a> Go online to view related products.												
 <b>Product Description</b>	<ul style="list-style-type: none"> <li>Lipofectamine® RNAiMAX Transfection Reagent is a proprietary formulation for transfecting small RNAs (e.g., siRNA, Silencer® Select siRNA, Stealth® RNAi, mirVana™ miRNA mimics and inhibitors) into a wide range of eukaryotic cells.</li> </ul>												
 <b>Important Guidelines</b>	<ul style="list-style-type: none"> <li>RNA-Lipofectamine® RNAiMAX 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.</li> <li>It is not necessary to remove complexes or change/add medium after transfection.</li> <li>Use 10 nM RNAi duplex as a starting point. BLOCK-iT™ Alexa Fluor® Red Fluorescent Oligo (Cat. no. 14750100) can be used to determine transfection efficiency.</li> </ul>												
 <b>Online Resources</b>	Visit our <a href="#">product page</a> for additional information and protocols. For support, visit <a href="http://www.lifetechnologies.com/support">www.lifetechnologies.com/support</a> .												



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

## Protocol Outline

- Plate cells so they will be 60-80% confluent at the time of transfection.
- Prepare RNA-lipid complexes.
- Add RNA-lipid complexes to cells.

## Lipofectamine® RNAiMAX Transfection Protocol

 See page 2 to view a typical RNAiMAX transfection procedure.

## Transfection Amounts

	96-well	24-well	6-well
Final siRNA used per well	1 pmol	5 pmol	25 pmol
Final Lipofectamine® RNAiMAX used per well	0.3 µL	1.5 µL	7.5 µL

## Reverse Transfection of RNAi

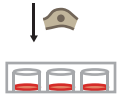


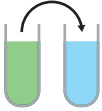

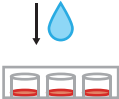

Reverse transfection is faster to perform than forward transfection and is the method of choice for high-throughput transfection. Perform reverse transfection by preparing complexes inside the wells, and then adding cells and medium. Because the cells and siRNA-lipid complexes are prepared on the same day, we recommended using 2.5× more cells than for a regular transfection method.

## Scaling Up or Down Transfections

## Limited Product Warranty and Disclaimer Details

## Typical RNAiMAX Transfection Procedure

Transfect cells according to the following table. The transfection is designed for one RNA amount combined with one amount of Lipofectamine® RNAiMAX. **The prepared mix is enough to have triplicates (96-well), duplicates (24-well), and single well (6-well) transfections, and account for pipetting variations.** For additional information on scaling your transfection reaction, see page 1.

Timeline			Steps	Procedure Details			
Day 0	1		Seed cells to be 60-80% confluent at transfection	Component	96-well	24-well	6-well
	2		Dilute Lipofectamine® RNAiMAX Reagent in Opti-MEM® Medium	Adherent cells	1–4 × 10 <sup>4</sup>	0.5–2 × 10 <sup>5</sup>	0.25–1 × 10 <sup>6</sup>
Day 1	3		Dilute siRNA in Opti-MEM® Medium	Opti-MEM® Medium	25 µL	50 µL	150 µL
	4		Add diluted siRNA to diluted Lipofectamine® RNAiMAX Reagent (1:1 ratio)	Lipofectamine® RNAiMAX Reagent	1.5 µL	3 µL	9 µL
	5		Incubate	Opti-MEM® Medium	25 µL	50 µL	150 µL
	6		Add siRNA-lipid complex to cells	siRNA (10 µM)	0.5 µL (5 pmol)	1 µL (10 pmol)	3 µL (30 pmol)
Day 2–4	7		Visualize/analyze transfected cells	Diluted siRNA	25 µL	50 µL	150 µL
				Diluted Lipofectamine® RNAiMAX Reagent	25 µL	50 µL	150 µL
				Incubate for 5 minutes at room temperature.			
				Component	96-well	24-well	6-well
				siRNA-lipid complex per well	10 µL	50 µL	250 µL
				Final siRNA used per well	1 pmol	5 pmol	25 pmol
				Final Lipofectamine® RNAiMAX used per well	0.3 µL	1.5 µL	7.5 µL
				Incubate cells for 1–3 days at 37°C. Then, analyze transfected cells.			