

Pub. No. MAN0010803 Rev D.0

	<b>Package Contents</b>	<b>Catalog Number</b>	<b>Size</b>
		<ul style="list-style-type: none"> <li>LMRNA001</li> <li>LMRNA003</li> <li>LMRNA008</li> <li>LMRNA015</li> <li>LMRNA150</li> </ul>	<ul style="list-style-type: none"> <li>0.1 mL vial</li> <li>0.3 mL vial</li> <li>0.75 mL vial</li> <li>1.5 mL vial</li> <li>15 mL vial</li> </ul>
	<b>Storage Conditions</b>	Store at 4°C (do not freeze).	
	<b>Required Materials</b>	<ul style="list-style-type: none"> <li>mRNA (0.5–5 µg/µL stock)</li> <li>mRNA positive control</li> <li>Opti-MEM™ Reduced Serum Medium</li> <li>Eppendorf tubes</li> </ul>	
	<b>Timing</b>	Preparation: 10 minutes Incubation: 10 minutes, 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™ MessengerMAX™ mRNA transfection reagent is a proprietary formulation that is optimized to deliver the highest amount of mRNA possible into neurons and a range of difficult-to-transfect primary cells.</li> </ul>	
	<b>Important Guidelines</b>	<ul style="list-style-type: none"> <li>mRNA-Lipofectamine™ MessengerMAX™ complexes must be made in serum-free medium, but can be added directly to cells in culture medium with/without serum/antibiotic.</li> <li>It is not necessary to remove complexes or change/add medium after transfection.</li> <li>The amount of Lipofectamine™ MessengerMAX™ reagent required for successful transfection varies depending on the cell type and passage number. Start any new transfection by testing the two recommended concentrations of Lipofectamine™ MessengerMAX™ Reagent to determine an optimum amount.</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> .	

## Protocol Outline

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

## Transfection Amounts

Component	96-well	24-well	6-well
Final mRNA per well*	100 ng	500 ng	2500 ng
Final Lipofectamine™ MessengerMAX™ Reagent per well	0.15 and 0.3 µL	0.75 and 1.5 µL	3.75 and 7.5 µL

\* To help ensure validity of transfection results, we strongly recommend the use of a positive control mRNA available from various third party vendors including TriLink BioTechnologies (e.g. EGFP mRNA).

## mRNA Synthesis

The Ambion™ mMESSAGE mMACHINE™ T7 Ultra Kit (Cat. no. AM1345) is recommended for synthesis of mRNA transcripts incorporating a 5' ARCA cap and 3' poly(a) tail.

Any plasmid with the gene of interest regulated by a T7 polymerase promoter can be used as the reaction template.

A PCR product of the gene of interest may also be used. See below for details:

### Forward Primer

GC TAATACGACTCACTATAGGG ACAG GCCACC ATG (gene specific sequence)

1 2 3 4 5 6

### Reverse Primer

Use the reverse complement sequence of the gene of interest.

## Primer Characteristics

## Genome Editing Applications

Lipofectamine™ MessengerMAX™ reagent increases the likelihood of successful cleavage and recombination with GeneArt™ CRISPR Nuclease mRNA (Cat. no. A25640) through highly efficient transfection, and ultimately maximizes the efficiency of genetic modifications and simplifies downstream processes.

## Scaling Up or Down Transfections

## Limited Product Warranty and Disclaimer Details

# Lipofectamine™ MessengerMAX™ Reagent mRNA Transfection Protocol

Transfect cells according to the following table. Use the indicated volume of transfection reagent at the two recommended doses as a starting point for optimization.

Volumes in each column are for a single well. Scale the volumes proportionally for additional wells.

Each reaction mix volume is for one well and accounts for pipetting variations.

Timeline		Steps
Day 0	1	 <b>Seed cells to be 70–90% confluent at transfection</b>
	2	 Diluted MessengerMAX™ Reagent Vortex 2–3 sec <b>Dilute MessengerMAX™ Reagent in Opti-MEM™ Medium (2 tubes) – Mix well</b>
Day 1	3	 <b>Incubate</b>
	4	 Diluted mRNA <b>Prepare Diluted mRNA master mix by adding mRNA to Opti-MEM™ Medium – Mix well</b>
	5	 <b>Add Diluted mRNA to each tube of Diluted MessengerMAX™ Reagent (1:1 ratio)</b>
	6	 <b>Incubate</b>
	7	 <b>Add mRNA-lipid complex to cells</b>
	8	 <b>Visualize/analyze transfected cells</b>

Procedure Details (Two Reaction Optimization)			
Component	96-well	24-well	6-well
Adherent cells	1–4 × 10 <sup>4</sup>	0.5–2 × 10 <sup>5</sup>	0.25–1 × 10 <sup>6</sup>
Opti-MEM™ Medium	5 µL × 2	25 µL × 2	125 µL × 2
Lipofectamine™ MessengerMAX™ Reagent	0.15 and 0.3 µL	0.75 and 1.5 µL	3.75 and 7.5 µL
Incubate diluted MessengerMAX™ Reagent in Opti-MEM™ Medium for 10 minutes at room temperature.			
Opti-MEM™ Medium	10 µL	50 µL	250 µL
mRNA (0.5–5 µg/µL)	0.2 µg	1 µg	5 µg
Diluted mRNA	5 µL	25 µL	125 µL
Diluted Lipofectamine™ MessengerMAX™ Reagent	5 µL	25 µL	125 µL
Incubate for 5 minutes at room temperature.			
Component (per well)	96-well	24-well	6-well
mRNA-lipid complex	10 µL	50 µL	250 µL
mRNA	100 ng	500 ng	2500 ng
Lipofectamine™ MessengerMAX™ Reagent	0.15 and 0.3 µL	0.75 and 1.5 µL	3.75 and 7.5 µL
Incubate cells for 1–2 days at 37°C. Then, analyze transfected cells.			