

I. Introduction

This protocol is provided for transfection with **Xfect Single Shots (Maxi)** (Cat. Nos. 631437, 631447), single-transfection aliquots of lyophilized Xfect Transfection Reagent supplied in individual tubes. Use this procedure to transfect DNA into mammalian cells **in a 10 cm dish format**. The amount of reagent in each tube is optimized for a single transfection in a 10 cm dish. Transfections can be carried out entirely in the presence of serum.

II. General Considerations

A. Storage and Handling

- Store Xfect Single Shots at -20°C in the supplied foil pouch containing the desiccant sachet.
- Make sure to return any unused Xfect Single Shots to the supplied foil pouch containing the desiccant sachet, and store at -20°C .

B. Mock Transfections

Use a plasmid that does not contain your gene of interest. You should include a source of nucleic acids to assemble with the Xfect polymer.

III. Transfection Protocol

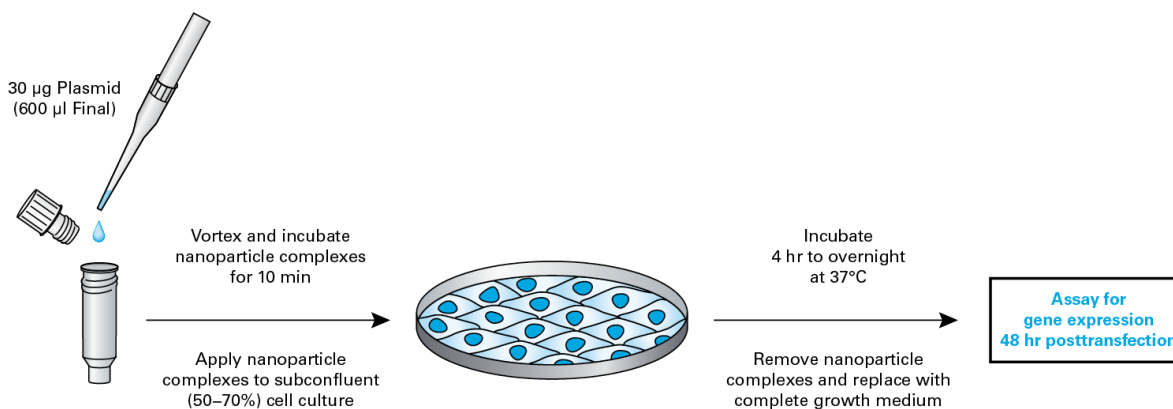


Figure 1. The Xfect Single Shots (Maxi) transfection protocol.

1. Prepare cells for transfection **in a 10 cm dish** as follows:

Adherent cells: One day prior to the transfection, plate cells in 8 ml of complete growth medium so that the cells will be 50–70% confluent at the time of transfection.

NOTE: The 8 ml volume of growth medium used here is important, since it affects the concentrations of the transfection reaction components.

Xfect™ Single Shots (Maxi) Protocol-At-A-Glance

- In a sterile microfuge tube, dilute 25 µg of your plasmid DNA with sterile water to a final volume of 600 µl. Mix thoroughly by vortexing.

NOTES:

- Always dilute your DNA in water prior to mixing with Xfect— do not add water and DNA separately to the Xfect Single Shot.
- Do not use less than 12.5 µg of DNA per 10 cm dish—25 µg of plasmid DNA works best for most cell lines.

- Add the 600 µl of diluted DNA to a tube of Xfect Single Shots, replace the cap, and vortex well at a high speed for 10 sec. The pellet should dissolve completely.
- Incubate the samples for 10 min at room temperature to allow nanoparticle complexes to form. After the 10 min incubation, centrifuge the tube for 2 sec to collect the contents at the bottom of the tube.

NOTE: Sample tubes can be inserted into 1.5 ml microfuge tubes for centrifugation.

- Add the entire 600 µl of nanoparticle complex solution dropwise to the cell culture prepared in Step 1. Rock the plate gently back and forth to mix.

NOTE: It is normal for the medium to change color slightly upon addition of nanoparticle complex solution.

- Incubate the cells at 37°C from 4 hr to overnight.

NOTE: A 4 hr incubation with Xfect-DNA nanoparticles is sufficient for optimal transfection. Incubation may be continued overnight for convenience but does not generally increase transfection efficiency. If you have sensitive cells, we recommend incubating for no more than 4 hr.

- Remove the nanoparticle complexes from the cells by aspiration, replace with 10 ml of fresh complete growth medium, and return the cells to the 37°C incubator until the time of analysis. Peak expression is typically reached 48 hr posttransfection.

Contact Us For Assistance	
Customer Service/Ordering	Technical Support
Telephone: 800.662.2566 (toll-free)	Telephone: 800.662.2566 (toll-free)
Fax: 800.424.1350 (toll-free)	Fax: 650.424.1064
Web: www.clontech.com	Web: www.clontech.com
E-mail: orders@clontech.com	E-mail: tech@clontech.com

Notice to Purchaser

Clontech® products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Clontech Laboratories, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech, the Clontech logo, and Xfect are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2014 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.