

# ProFection® Mammalian Transfection System

INSTRUCTIONS FOR USE OF PRODUCT E1200.

Quick  
PROTOCOL

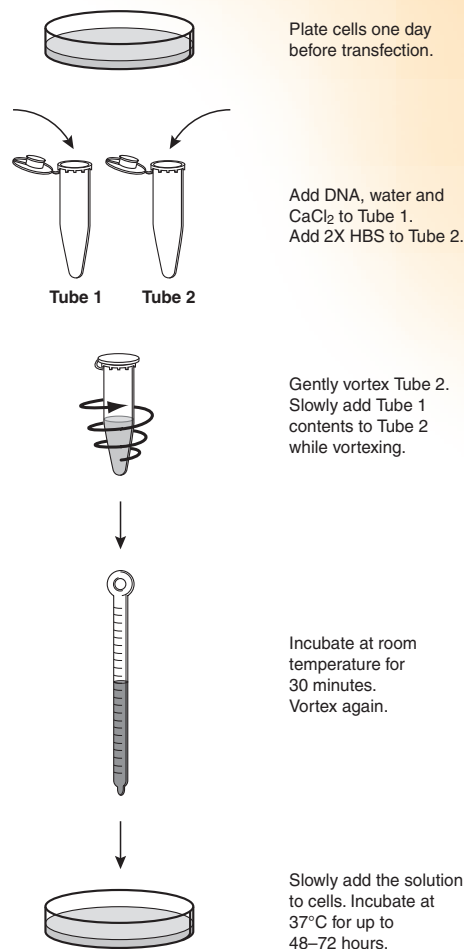
## Calcium-Phosphate Transfection Protocol

1. Plate cells one day before the transfection. A general guideline for plating density is  $8 \times 10^5$  cells per 100mm plate.
2. Three hours before the transfection, replace cell culture medium with fresh growth medium.
3. Thaw all transfection reagents. Warm to room temperature and mix thoroughly.
4. For each transfection, prepare two tubes. Add DNA and water, followed by  $\text{CaCl}_2$ , to Tube 1. Add 2X HBS to Tube 2 (see table below).

	Per 60mm Plate	Per 100mm Plate
<b>Tube 1</b>		
DNA	6–12 $\mu\text{g}$	10–20 $\mu\text{g}$
2M $\text{CaCl}_2$	37 $\mu\text{l}$	62 $\mu\text{l}$
sterile, deionized water to a final volume of	300 $\mu\text{l}$	500 $\mu\text{l}$
<b>Tube 2</b>		
2X HBS	300 $\mu\text{l}$	500 $\mu\text{l}$

5. In a tissue culture hood, gently vortex the 2X HBS. Slowly add the DNA solution in Tube 1 dropwise to the HBS in Tube 2 while vortexing.
6. Incubate the combined solution at room temperature for 30 minutes.
7. Vortex again, then immediately add the solution, dropwise, to cells. Swirl plate to distribute. Incubate at 37°C with  $\text{CO}_2$  for up to 48–72 hours.
8. After incubation cells may be harvested or treated with selective medium.

For additional protocol information, see Technical Manual #TM012, available online at: [www.promega.com/tbs](http://www.promega.com/tbs)



3089MA10\_0A

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA. Revised 3/09.  
Part #9FB049

©2000–2009 Promega Corporation. All Rights Reserved.