

293-Free[™] Transfection Reagent

Table of Contents

| | |
|---------------------------------|---|
| About the Kits | 2 |
| Description | 2 |
| Storage | 2 |
| General Considerations | 3 |
| Cell Transfection Protocol..... | 4 |
| Transfection Procedure | 4 |

© 2011 EMD Chemicals Inc., an affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. Novagen[®] is a registered trademark Merck KGaA, Darmstadt, Germany. 293-Free[™] is trademark of Merck KGaA, Darmstadt, Germany. Opti-Pro[™] is a trademark of Life Technologies, Corporation.

USA and Canada

Tel (800) 628-8470
bioscienceshelp@
emdchemicals.com

France
Freephone
0800 126 461

Germany
Freecall
0800 100 3496

Ireland
Toll Free
1800 409 445

United Kingdom
Freephone
0800 622 935

**All other
European Countries**
+44 115 943 0840

All Other Countries

Contact Your Local Distributor
www.merck4biosciences.com
bioscienceshelp@
emdchemicals.com

techservice@merckbio.eu

www.merck4biosciences.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.

About the Kits

| | | |
|--------------------------------|-----------|---------|
| 293-Free™ Transfection Reagent | 1 x 1 ml | 72181-3 |
| | 5 x 1 ml | 72181-4 |
| | 10 x 1 ml | 72181-5 |

Description

293-Free™ Transfection Reagent has a unique polycationic liposomal formulation and is designed expressly for transfection of HEK293 cells grown in suspension culture. It is ideal for mammalian protein production. Derived from non-animal sources, 293-Free Transfection Reagent gives minimal cellular toxicity. It is provided as a sterile, ready-to-use solution. Benefits of this proprietary polycationic liposomal formulation include:

- Optimized for transient transfections of HEK293 suspension cultures
- Minimal cellular toxicity
- Derived from non-animal sources
- Protocol easily scales up for production
- Compatible with both serum-containing and serum-free media
- Eliminates the need for media changes

Each 1 ml of 293-Free™ Transfection Reagent is sufficient to transfect a 1-liter culture.

Storage

Store tightly capped at 4°C.

General Considerations

- Use only high quality, endotoxin-free DNA. DNA should be at a concentration of 0.5–1 µg/µl. If the DNA concentration is lower, decrease the volume of serum-free medium or PBS in the transfection protocol to compensate for the larger volume of DNA.
- Propagate and maintain cell cultures on an orbital shaker at 125 rpm at 37°C (8% CO₂).
- Passage cells regularly (e.g., every 2–3 days). Avoid high density. Use only rapidly proliferating cells for transfection. To ensure reproducibility, keep cell growth conditions and density consistent.
- 293-Free™ Transfection Reagent is compatible with both serum-containing and serum-free media.

Note: Do not include serum and antibiotics during the formation of the transfection reagent/DNA complex.

Cell Transfection Protocol

The following protocol is optimized for transfection of 25 ml HEK-293 cells in suspension in a 125-ml polycarbonate Erlenmeyer flask, however, the amounts of reagent and DNA can be varied if necessary. For other culture sizes, the reagent and DNA amounts can be scaled up or down proportionately. See Table 1 below.

Transfection Procedure

1. The day before transfection, passage HEK293 suspension cells at approximately 0.5×10^6 cells/ml. Incubate at ~125 rpm at 37°C (8% CO₂) overnight. The cell density should be approximately $1.0\text{--}1.5 \times 10^6$ cells/ml at the time of transfection.
2. On the day of transfection, dilute the cells to 1×10^6 cells/ml and place 25 ml in a 125 ml shake flask.
3. Place 1 ml serum-free MEM, Opti-Pro™ SFM, or PBS into a sterile tube. Add 12.5 µg DNA. Mix.
Note: Do not use the serum-free culture medium in which HEK293 cells were grown.
4. Add 25 µl 293-Free Transfection Reagent. Mix.
5. Incubate transfection mixture at room temperature for 15 minutes to allow 293-Free Transfection Reagent/DNA complex formation.
6. Add entire volume of transfection mixture drop wise to prepared cells.
7. Incubate cultures for 48–168 hours, at ~125 rpm at 37°C (8% CO₂).
8. Harvest cells for protein purification, characterization, or reporter assays.

Table 1. Preparation of Transfection Mixture for Various Culture Sizes

| Transfection of Plasmid DNA | Culture Size | | | |
|---|--------------|-----|-----|------|
| | 25 | 250 | 500 | 1000 |
| Culture Volume (ml) | 25 | 250 | 500 | 1000 |
| Number of cells ($\times 10^6$) | 25 | 250 | 500 | 1000 |
| Volume of MEM or PBS in the transfection mixture (ml) | 1 | 10 | 20 | 40 |
| Volume of 293-Free Transfection Reagent (µl) | 25 | 250 | 500 | 1000 |
| Amount of plasmid DNA (µg) | 12.5 | 125 | 250 | 500 |