

Reverse Transfection (RTF) Optimization Plates

This protocol is for the Dharmacon RTF Optimization Plates.

Introduction:

Optimizing RNA interference (RNAi) experiments is a critical component to successful gene silencing experimentation. The Reverse Transfection Format (RTF) Optimization plates are designed for the optimization of siRNA experiments prior to using Dharmacon™ siGENOME™ or Dharmacon™ ON-TARGETplus™ RTF siRNA Libraries.

Product Description:

Once the optimum Dharmacon™ DharmaFECT™ siRNA transfection reagent (or other transfection reagent of choice) has been determined for your cell line, the RTF Optimization Plates allow simple and flexible determination of (a) the best negative control siRNA, (b) optimal cell density and (c) DharmaFECT transfection reagent amount for the cells and assay being used. The pre-plated controls provided at 6.25 pmol per well in triplicate plate sets are described in the accompanying Product Insert.

Each RTF Optimization Plate is provided with pre-plated controls as described below (Figure 1).

- Rows A-D contain four different Non-targeting siRNAs.
- Rows E contains a Non-targeting siRNA pool.
- Row F contains a validated positive control siRNA pool targeting the housekeeping gene Cyclophilin B (PPIB).
 - a. Row G is recommended for mock transfected (or transfection reagent alone) to determine if there is any effect of the transfection reagent on cell viability and/or mRNA expression.
 - b. Row H is recommended for untreated cells and serves as 100% viability control and/or 100% mRNA level control.

Handling Precautions:

Oligonucleotides are susceptible to enzymatic degradation by nucleases and to chemical degradation by extreme pH and temperatures. Wearing gloves and maintaining nuclease-free conditions when handling the oligonucleotides is highly recommended.

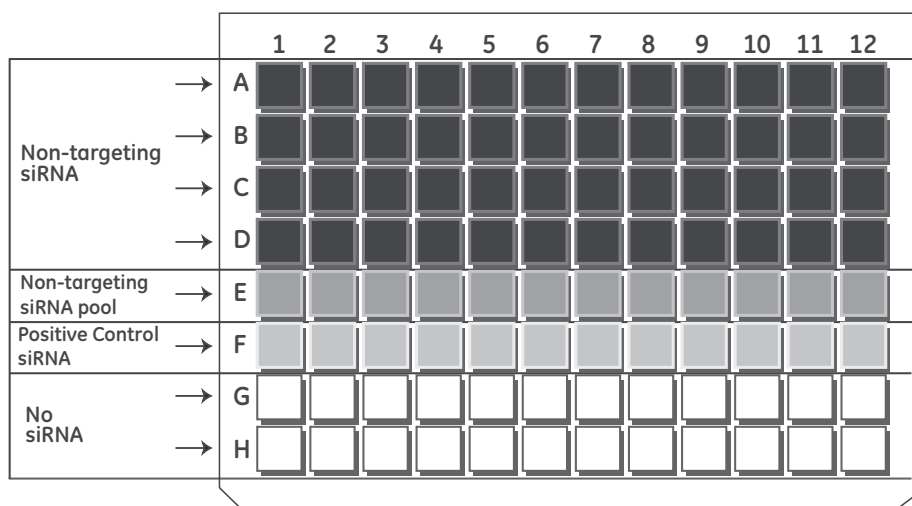


Figure 1. RTF Optimization Plate Layout.



Additional Materials Required for Use:

- DharmaFECT Transfection Reagent 1, 2, 3 or 4 (Cat #T-2001, T-2002, T-2003, T-2004)
- DharmaFECT Cell Culture Reagent (DCCR; 100 mL bottle, Cat #B-004500-100), DCCR is provided with Optimization Plates with catalog numbers ending in -D
- Equipment, reagents and supplies needed for cell culture
- Equipment, reagents and supplies needed for transfection efficiency assessment
- Deep-well plate for preparation of the Master Dilution Plate [for example Nunc™ 96 DeepWell Plate, 1.0 mL, Polystyrene, Sterile, Transparent, from Fisher Scientific (Cat #12-565-553)]

Protocol:

The recommended experimental format for RTF Optimization Plates is shown in Figure 2. The Dharmacon DharmaFECT transfection reagent that has been predetermined to be optimum for the cells (from use in conventional transfection experiments) is tested at two volumes with two cell numbers per well. An alternate method could assess four DharmaFECT reagent volumes with a single cell number per well.

The instructions below describe how to construct the plate format shown in Figure 2. All manipulations should be performed under sterile, RNase-free conditions. The final siRNA concentration during transfection is 50 nM when using this protocol.

1. Allow RTF Optimization Plates and DharmaFECT Cell Culture Reagent (DCCR) to equilibrate to room temperature.
2. Prepare the Rehydration Solution in two 15 mL tubes as described below and in Figure 2. Tube X will prepare reagent for a final volume of 0.15 μL DharmaFECT reagent per well and Tube Y will prepare reagent for a final volume of 0.075 μL DharmaFECT per well.
 - a. Dispense 3519 μL DCCR into Tube X and 1180 μL DCCR into Tube Y.
 - b. Add 21 μL DharmaFECT Transfection Reagent to Tube X and mix well by pipetting up and down.
 - c. Transfer 1180 μL from Tube X to Tube Y. Mix well.

Tubes	Volume of DCCR (μL)	Volume of DharmaFECT (μL)
X	3519	21
Y	1180	None

- d. Prepare Master Dilution Plate by dispensing 330 μL from each Rehydration Solution tube to the appropriate wells (shown below) in a deep-well plate, and 330 μL of DCCR into wells H1 and H2. Following the recommended protocol will facilitate convenient and accurate dispensing of the Rehydration Solution to the RTF Optimization Plate. The Rehydration Solution may be stored under sterile conditions at room temperature for up to 2 hours prior to use.
3. Rehydrate wells of RTF Optimization Plate by transferring 25 μL of DCCR or Rehydration Solution from the appropriate wells of the Master Dilution Plate as described below and in Figure 2.

Tubes	Dispense into Wells of Master Dilution Plate
X	A1 – G1
Y	A2 – G2
DCCR	H1 and H2

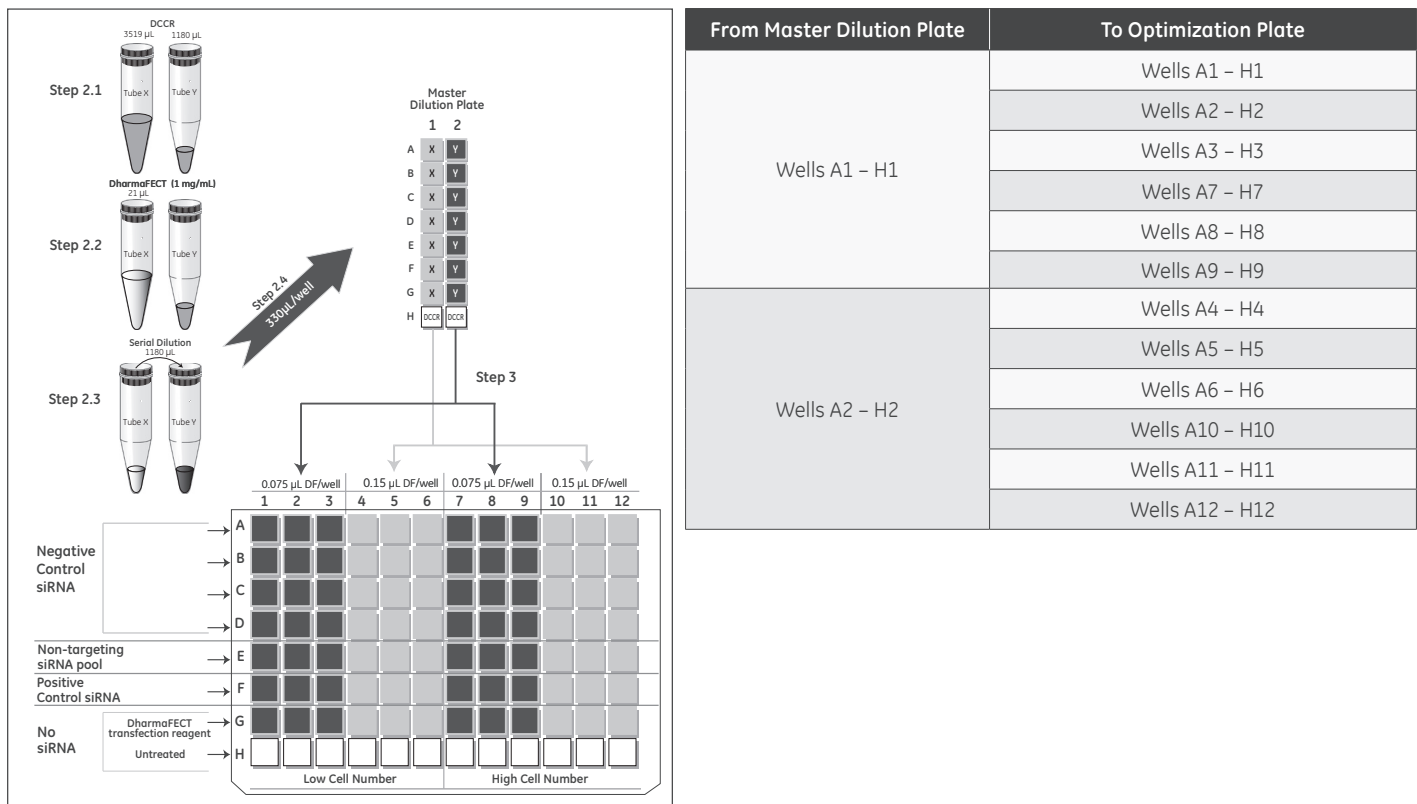


Figure 2. RTF Optimization Plate Rehydration Guide.

Note: The Rehydration Solution may not coat the bottom of the wells but rather “ring” the bottom of the well. This is expected and does not interfere with rehydration of the siRNA.

4. After the transfer of Rehydration Solution to the RTF plate is completed, allow the DharmaFECT transfection reagent to complex with siRNA by incubating the rehydrated RTF Optimization plates at room temperature for 15 to 90 minutes. Optimum incubation time is 30 minutes.
5. Add cells to rehydrated RTF Optimization Plates:
 - a. Prepare at least 5.5 mL of cell suspension in antibiotic-free complete medium. The cell concentration should be such that when 100 μ L is dispensed into each well, the required cell number per well is achieved. For example, to add 1×10^3 cells per well, the cell suspension should have 1×10^4 cells per mL, and to add 2×10^3 cells per well, the cell suspension should have 2×10^4 cells per mL.
 - b. Add 100 μ L of the low cell number cell suspension to Columns 1-6 and add 100 μ L of the high cell number cell suspension to columns 7-12. The final volume in each well will be 125 μ L. The final concentration of siRNA is 50 nM.
6. Incubate the RTF Optimization Plates under appropriate culture conditions (for example 37 °C, 5 % CO₂).
 - For the most consistent results, we recommend that RTF plates be incubated for at least 2 days before performing the transfection efficiency and screening assays of choice.
7. Perform the transfection efficiency assay and the cell viability assay of choice.

Nunc is a trademark of Thermo Fisher Scientific, Inc. GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon is a trademark GE Healthcare companies. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2015 General Electric Company—All rights reserved. Version published January 2015. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK



GE Healthcare

Orders can be placed at:
gelifesciences.com/dharmacon

Customer Support: cs.dharmacon@ge.com
Technical Support: ts.dharmacon@ge.com or
1.800.235.9880; 303.604.9499 if you have any questions.

VI-0115