

In vitro DNA Transfection using Thermo Scientific TurboFect™ Transfection Reagent

This protocol is for the *In vitro* DNA Transfection using Thermo Scientific TurboFect™ Transfection Reagent.

Reagents to be Supplied by the User:

Serum free DMEM, RPMI or other growth medium. The presence of antibiotics in the medium has no effect on transfection efficiency. The protocol below is provided for 24-well plate. Quantities and volumes should be scaled up or down according to the number of cells/or wells to be transfected (see Table below for guidelines).

1. In each well, seed $\sim 5 \times 10^4$ adherent cells or $\sim 5 \times 10^5$ suspension cells 24 hours prior to transfection.

Note

- The recommended confluency for adherent cells on the day of transfection is 50-70%.
 - Suspension cells should be in logarithmic growth phase at the time of transfection.
2. Dilute 1 μ g of DNA in 100 μ L of serum-free DMEM or other growth medium.
 3. Add 2 μ L of TurboFect™ to the diluted DNA and mix the solution by pipetting.
 4. Incubate 15-20 minutes at room temperature.

Note

- Prepare immediately prior to transfection.
 - We recommend starting with 1 μ g of DNA and 2 μ L of TurboFect per well in a 24-well plate (see Table below).
 - Subsequent optimization may further increase transfection efficiency depending on the cell line and transgene used.
5. Dilute 1 μ g of DNA in 100 μ L of serum-free DMEM or other growth medium.
 6. Gently rock the plate to achieve even distribution of the complexes.
 7. Incubate at 37 °C in a CO₂ incubator.
 8. Analyze transgene expression 24-48 hours.
For stable transfection cells should be grown in selective medium for 10-15 days. Plates can be centrifuged for 2-5 minutes at 200 \times g to facilitate sedimentation of cells to the bottom of the plate.

Table. Scale-up /down ratios for transfection with TurboFect™ Transfection Reagent.

Tissue Culture Vessel	Growth Area, cm ² /well	Media, mL	Adherent cells to seed the day before transfection*	DNA, µg** (µL***)	Volume of TurboFect (µL)**	
					Recommended	Range
96-well plate	0.3	0.2	0.5-1.20 × 10 ⁴	0.2 (20)	0.4	0.3-0.6
48-well plate	0.7	0.5	1.0-3.0 × 10 ⁴	0.5 (50)	1.0	0.5-1.4
24-well plate	2.0	1.0	2.0-6.0 × 10 ⁴	1.0 (100)	2.0	1.0-2.8
12-well plate	4.0	2.0	0.4-1.2 × 10 ⁵	2.0 (200)	4.0	2.0-6.0
6-well plate	9.5	4.0	0.8-2.4 × 10 ⁵	4.0 (400)	6.0	4.0-8.0
60 mm plate	20.0	6.0	2.0-6.3 × 10 ⁵	6.0 (600)	12.0	8.0-16.0

*These numbers were determined using HeLa cells. Actual value depends on the cell type.

** Amount of DNA and TurboFect™ Transfection Reagent used may require optimization.

*** The volume of the DNA solution should represent 1/10 of the total volume of the culture medium.

www.thermoscientific.com/onebio

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Europe

Customer Service
cs.molbio.eu@thermofisher.com

Technical Support
ts.molbio.eu@thermofisher.com

Tel 00800 222 00 888
Fax 00800 222 00 888

United States

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 235 9880
Fax 800 292 6088

Canada

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 340 9026
Fax 800 472 8322

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific