

**Introduction**

This protocol shows optimum transfection condition using HilyMax in U937 cells. To tranfect U937 cells in 24-well plate, follow "Optimum Condition for Transfection" and "Transfection Procedure". When using the other vessel, refer to Table 2 and adjust the amounts of cells, medium, DNA and HilyMax in proportion to the relative surface area.

※Important Note※

Optimum Transfection condition is possibly changed by passage number and culture condition. If transfection efficiency is low by followed this protocol, refer to "Transfected Result by HilyMax" and "Troubleshooting".

**Optimum Condition for Transfection (for 24-well plate)**

Table 1 Optimum condition for transfection to U937 cells

Number of cells	3.0x10 <sup>5</sup> cells/well	
DNA-HilyMax complex formation	Serum-free medium	30 μl
	DNA	1 μg
	HilyMax	7.0-9.0 μl
	Incubation time	15 min
Medium change after transfection	Not Determined	

**Transfection Procedure (for 24-well plate)**

Cell preparation

- ↓ Adjust the concentration of cells to be 3.0x10<sup>5</sup> cells/well confluent in 0.5 ml of growth medium.
- ↓ Inoculate the cell suspension onto the 24-well plate.
- ↓ Incubate cells in CO<sub>2</sub> incubator for 24 hr.

Transfection

- ↓ Form the DNA-HilyMax complex
  - Add the serum-free medium(without antibiotics) 30 μl/well in a sterile plastic tube
  - Add plasmid DNA 1.0 μg/well and mix by gentle pipetting
  - Add HilyMax 7.0-9.0 μl/well and mix by gentle pipetting
  - Incubate the mixture of DNA and HilyMax solution at room temperature for 15 minutes
- ↓ Add DNA-HilyMax complex to cells in each well and mix by gentle shaking the plate
- ↓ Incubate cells in CO<sub>2</sub> incubator for 18-48 hr

Assay

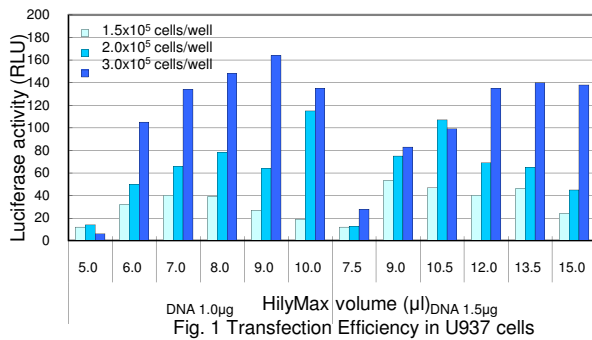
- ↓ Measure protein expression

**Transfection in Various Vessels**

Table 2 Transfection condition in various vessels

Culture Vessel	Culture of Cells		Formation of DNA-HilyMax complex		
	Surface Area	Plating Medium	Serum-free Medium	DNA	HilyMax
96 -well	0.3 cm <sup>2</sup>	0.1 ml	10 μl	0.2 μg	0.8-1.2 μl
24 -well	1.9 cm <sup>2</sup>	0.5 ml	30 μl	1.0 μg	7.0-9.0 μl
12 -well	3.8 cm <sup>2</sup>	1.0 ml	60 μl	2.0 μg	14.0-18.0 μl
6 -well	9.2 cm <sup>2</sup>	2.0 ml	120 μl	4.0 μg	28.0-36.0 μl
35 -mm	8.0 cm <sup>2</sup>	2.0 ml	120 μl	4.0 μg	28.0-36.0 μl
60 -mm	21.0 cm <sup>2</sup>	5.0 ml	300 μl	10.0 μg	70.0-90.0 μl
100 -mm	58.0 cm <sup>2</sup>	15.0 ml	900 μl	30.0 μg	210.0-270.0 μl

**Transfected result by HilyMax**



U937 cells were incubated for 24 hr and transfected pGL3 control vector (Promega) using HilyMax in each conditions. Transfection efficiency (Luciferase activity) was measured in 24 hr after transfection.

U937 cells were cultured in RPMI1640 medium(Gibco) containing 10%FBS(Gibco) for about 2 weeks after thawing.

**Troubleshooting**

**-Low Transfection Efficiency-**

- Change the DNA(μg):HilyMax(μl) ratio to 1:10-1:12.
- Increase the mass of DNA up to 1.5-2.0 times and change the DNA(μg):HilyMax(μl) ratio to 1:7-1:9.

**-High cellular Toxicity-**

- Change the DNA(μg):HilyMax(μl) ratio to 1:3-1:6.
- Decrease the mass of DNA down to half and change the DNA (μg):HilyMax(μl) ratio to 1:2-1:7.

**-Check the Material and Condition-**

- Was HilyMax Reagent dissolved completely when HilyMax was Prepared?
- Was incubation time of cells after transfection optimum for cells and plasmid?
- Was DNA-HilyMax complex formed in medium without serum and antibiotics?