

General Information

HilyMax is a newly developed gene transfection reagent. The transfection reagent forms a liposome to be used on highly efficient gene transfection to a wide variety of cells. HilyMax can be utilized to siRNA transfection as well. Since serum in the growth medium does not interfere with the transfection using HilyMax, no exchange of the media during the transfection is required. HilyMax does not contain any biological components that might interfere with the transfection.

Kit Contents

- HilyMax Reagent x 1
- Lipoform Buffer 1.0 ml x 1

Storage Condition

Kit: Store at 0-5°C.

Reconstituted HilyMax solution: Store at -20°C. The reconstituted solution is stable at -20°C for 6 months.

**For frequent usage of the reconstituted HilyMax solution, store at 0-5°C and use the solution within 1 month.*

Preparation of HilyMax

Add Lipoform Buffer 1.0 ml to HilyMax Reagent and mix the solution well by vortex for 30 seconds or pipetting until the reagent is dissolved completely.

⚠ Check if HilyMax reagent is completely dissolved. If the insoluble form remains in the solution, continue the vortex or pipetting until the reagent is completely dissolved.

Precautions

Transfection efficiency varies significantly with types and densities of cells. We strongly recommend to search or seek a transfection condition for your best performance with HilyMax (see "Optimization Protocol" on reverse page).

The optimized protocol for following 18 cell lines are available at www.dojindo.com.

A549 cell, Caco2 cell, CHO cell, COS-7 cell, HEK293 cell, HeLa cell, HepG2 cell, K562 cell, L6 cell, LNCap cell, MCF7 cell, MDCK cell, Neuro2a cell, NIH3T3 cell, PC3 cell, PC12 cell, U937 cell, Vero cell are available. (Jan. 10, 2011)

General Protocol

Transfection procedure for a 24-well plate

1. Cell preparation

Adherent cells

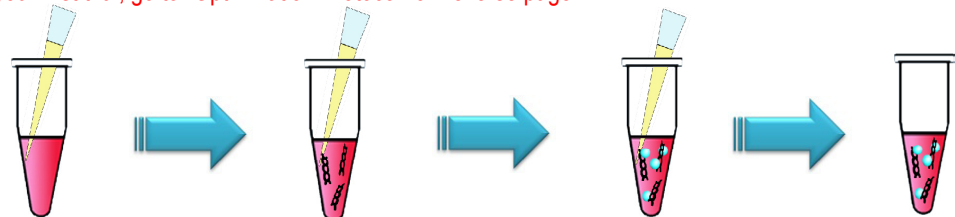
Prepare cell culture plate with 40-90% cell density.^{a)}

Non-adherent cells

Prepare $0.2-3.2 \times 10^6$ cells/ml cell suspension with cell culture medium.^{b)} Add 0.5 ml of cell suspension to each well.

2. DNA-HilyMax complex formation^{c)}

⚠ For your first trial, go to "Optimization Protocol" on reverse page.



Step 1.

Add 30 μ l serum-free medium^{d)} to a plastic tube.

Step 2.

Add plasmid DNA (0.5-1.5 μ g) to the tube and mix by pipetting.

Step 3.

Add HilyMax to prepare DNA (μ g):HilyMax (μ l)=1:2-1:7 and mix by pipetting.

Step 4.

Incubate the tube at room temperature for 15 minutes.^{e)}

3. Addition of complex to cells

Add DNA-HilyMax complex to cell culture well.

4. Transfection

Incubate the plate at 37°C in a CO₂ incubator for 4 hours.^{f)} Change the medium and continue incubation for 24-72 hours.

5. Downstream experiment

Reporter gene assay, protein isolation, protein analysis, etc.

a) 40-90% cell density means that 40-90% of the total surface area is covered with cells.

b) Serum does not interfere with the transfection.

c) The appropriate amount of DNA and HilyMax for various cell densities are indicated in Table 1.

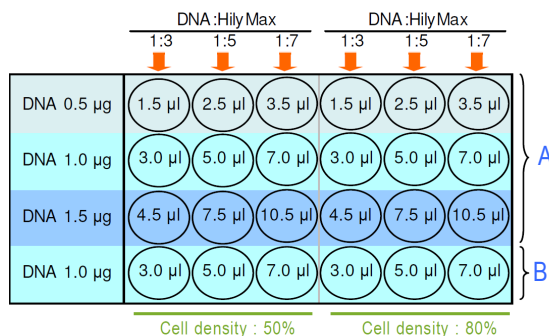
d) Serum and antibiotics in the medium interfere with the DNA-HilyMax complex formation. However Opti-MEM, DMEM, and MEM can be used for transfection.

e) Excess incubation time may cause a lower transfection efficiency.

f) A medium change increases the transfection efficiency and decreases the cytotoxicity for most of cell lines.

Optimization of transfection for a 24-well plate

The transfection efficiency may increase 2-5 times by executing the following optimization step. The amount of DNA and HilyMax, cell density, and medium change are the most significant factors to optimize your transfection efficiency. We strongly recommend you carry out optimization protocol for your first trial.



Transfection Procedure (for a 24-well plate)

1. Prepare cell culture plate with 50% and 80% cell density.
2. Add serum-free medium*¹ (30 µl) to a plastic tube.
3. Add plasmid DNA*² (0.5-1.5 µg) to the tube and mix.
4. Add HilyMax*² (1.5-10.5 µl) to the tube and mix.
5. Incubate the tube at room temperature for 15 minutes.
6. Add DNA-HilyMax complex to cell culture well (step 1).
7. Incubate at 37°C in CO₂ incubator in 24-48 hours incubation.
*Change the medium for the row B after 4 hours incubation.
8. Measure the transfection efficiency.

Figure. Example of plate arrangement for optimization

Value in each well indicates the volume of HilyMax.
 A: Without Medium Change.
 B: With Medium Change 4 hours after Transfection.

If you are using other size of vessels, change the amount of medium, DNA, and HilyMax.

*¹ Refer to Table 2 (Serum-Free Medium).

*² Multiply the coefficient in Table 1 by the amount of DNA and HilyMax.

Table 1 Coefficient for other vessels

96-well plate : x0.2	12-well plate : x2	6-well plate (35-mm dish) : x4
60-mm dish : x8	100-mm dish : x24	

Transfection Condition for Various Vessels

Table 2 shows suitable amount of medium, DNA, and mixing ratio of DNA and HilyMax solution against each vessel.

Table 2 Transfection condition for various vessels

Culture Vessel	Surface Area	Plating Medium	Serum-Free Medium	DNA	DNA(µg) : HilyMax(µl)
96 -well	0.3 cm ²	0.1 ml	10 µl	0.1-0.3 µg	1:2-1:7
24 -well	1.9 cm ²	0.5 ml	30 µl	0.5-1.5 µg	1:2-1:7
12 -well	3.8 cm ²	1.0 ml	60 µl	1.0-3.0 µg	1:2-1:7
6 -well	9.2 cm ²	2.0 ml	120 µl	2.0-6.0 µg	1:2-1:7
35 -mm	8.0 cm ²	2.0 ml	120 µl	2.0-6.0 µg	1:2-1:7
60 -mm	21.0 cm ²	5.0 ml	300 µl	5.0-15.0 µg	1:2-1:7
100 -mm	58.0 cm ²	15.0 ml	900 µl	15.0-45.0 µg	1:2-1:7

Troubleshooting

Transfection efficiency is low.

1. The volume of HilyMax is not enough. Increase the HilyMax volume.
2. Cell density is too high. Reduce the cell density. The appropriate cell density for transfection is about 40-90%.
3. HilyMax reagent is not be dissolved completely. Please check that the HilyMax solution is homogeneous.
4. Incubation time for the preparation of HilyMax and DNA complex is too long.
5. Your culture medium for DNA-HilyMax complex formation contain serum and/or antibiotics. Please use serum and antibiotics free medium for the complex formation.

Toxicity is high. Most cells died.

1. Reduce the amount of DNA and/or HilyMax prepare the complex.
2. Cell density is too low. Increase the cell density. The appropriate cell density for transfection is about 40-90%.

FAQ

Are there transfection conditions for specific cell lines available?

Please contact our customer service at 1-877-987-2667 or info@dojindo.com.

Does frequent freeze and thaw of HilyMax solution affect transfection efficiency?

No. Freeze/thaw process is repeatable without reducing the efficiency.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories
 2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto
 861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525
 E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.
 Tel: +1-301-987-2667 Web: http://www.dojindo.com/
 Dojindo EU GmbH
 Tel: +49-89-3540-4805 Web: http://www.dojindo.eu.com/
 Dojindo China Co., Ltd
 Tel: +86-21-6427-2302 Web: http://www.dojindo.cn/