

Introduction

This protocol shows optimum transfection condition using HilyMax in COS-7 cells. To tranfect COS-7 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

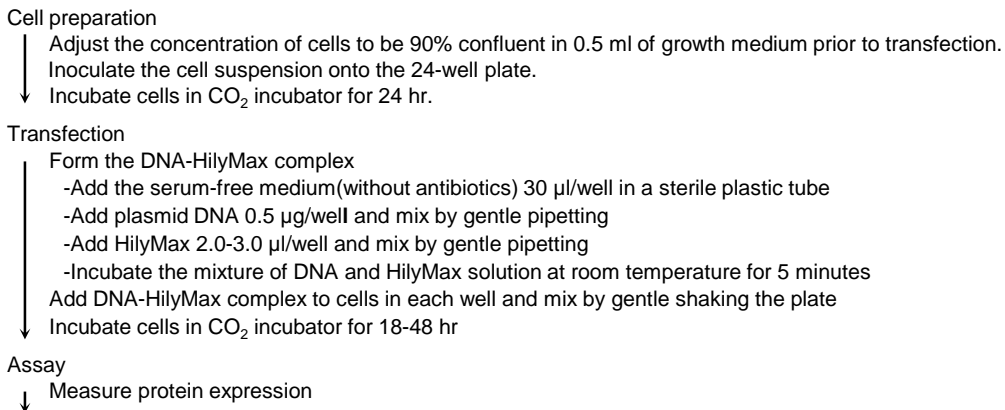
Optimum Transfection condition is possibly changed by passage number and culture condition. If transfection efficiency is low by followed this protocol, refer to "Transfected Result by HilyMax" and "Troubleshooting".

Optimum Condition for Transfection (for 24-well plate)

Table 1 Optimum condition for tranfection to COS-7 cells

Cell Density	90%	
DNA-HilyMax complex formation	Serum-free medium	30 μ l
	DNA	0.5 μ g
	HilyMax	2.0-3.0 μ l
	Incubation time	5 min
Medium change after transfection	Not Necessary	

Transfection Procedure (for 24-well plate)

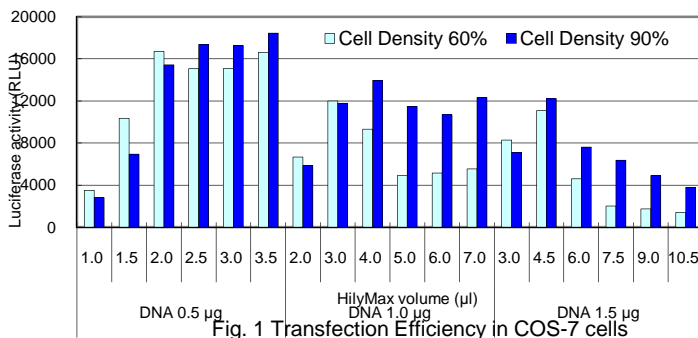


Transfection in Various Vessels

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 ²	0.1 ml	10 μ l	0.1 μ g	0.4-0.6 μ l
24 -well	1.9 cm ²	0.5 ml	30 μ l	0.5 μ g	2.0-3.0 μ l
12 -well	3.8 cm ²	1.0 ml	60 μ l	1.0 μ g	4.0-6.0 μ l
6 -well	9.2 cm ²	2.0 ml	120 μ l	2.0 μ g	8.0-12.0 μ l
35 -mm	8.0 cm ²	2.0 ml	120 μ l	2.0 μ g	8.0-12.0 μ l
60 -mm	21.0 cm ²	5.0 ml	300 μ l	5.0 μ g	20.0-30.0 μ l
100 -mm	58.0 cm ²	15.0 ml	900 μ l	15.0 μ g	60.0-90.0 μ l

Transfected result by HilyMax



COS-7 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. COS-7 cells were cultured in DMEM medium(Gibco) containing 10%FBS(Gibco) for about 2 weeks after thawing.

60% confluent: 0.7×10^5 cells/well 90% confluent: 1.1×10^5 cells/well

Troubleshooting

-Low Transfection Efficiency-
Change the DNA(μ g):HilyMax(μ l) ratio to 1:7-1:9.
Increase the mass of DNA up to 2.0-3.0 times and change the DNA(μ g):HilyMax(μ l) ratio to 1:4-1:6.

-High Cellular Toxicity-
Change the DNA(μ g):HilyMax(μ l) ratio to 1:2-1:3.

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