

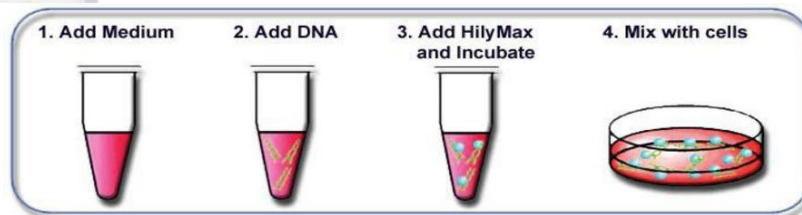
HilyMax

FEATURES

- ◇ Good Transfection Efficiency in Various Cell Lines
- ◇ Low Cytotoxicity
- ◇ Cost Effective



H357-10, 1x1mL
 H357-15, 5x1mL
 H357-20, 10x1mL



HilyMax is a newly developed gene transfection reagent that forms a liposome to be used for highly efficient gene transfection to a wide variety of cells. Serum in growth medium does not interfere with transfection performances by using HilyMax, HilyMax does not contain any biological components that might interfere with transfections.

TRANSFECTION EFFICIENCY IN VARIOUS CELL LINES

Cell line	Efficiency	Cell line	Efficiency
3T3-L1	30%	L6	30%
A549	50%	LNCap	30%
CHO	90%	MCF-7	70%
COS7	40%	MDCK	20%
HC	50%	MG63	20%
HEK293	60%	Neuro2a	70%
HeLa	70%	NIH3T3	70%
HepG2	10%	PC3	70%
Jurkat	3%	U1SMC	10%
K562	30%	Vero	40%

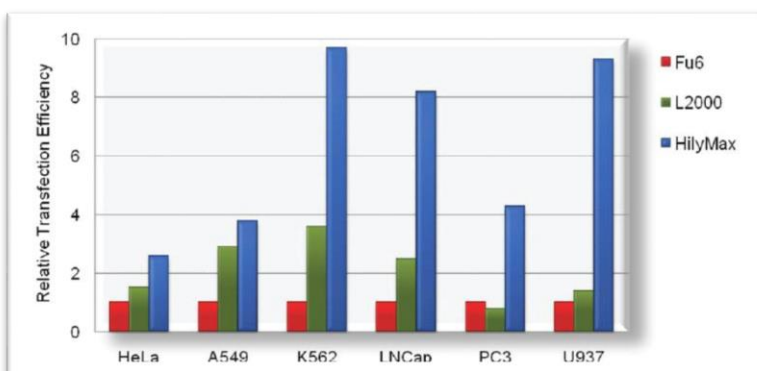
OPTIMIZED PROTOCOLS ARE AVAILABLE FOR BELOW CELL LINES!

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		

To get protocol, please visit us online at www.dojindo.eu.com

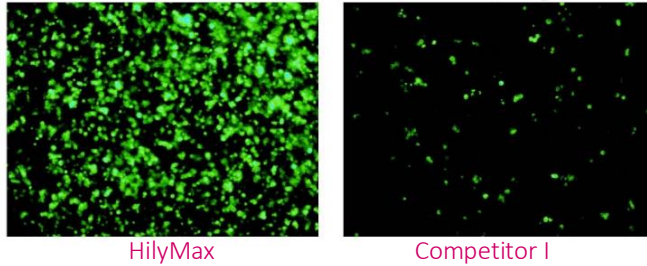


COMPARISON WITH OTHER SUPPLIERS



The transfection efficiency of HilyMax is higher than commercially available reagents. GFP-expressed DNA was transfected using HilyMax and the other transfection reagents. The highest transfection efficiencies were observed when HilyMax was used.

COMPARISON IN TRANSFECTION USING S2 CELLS



Culture Conditions

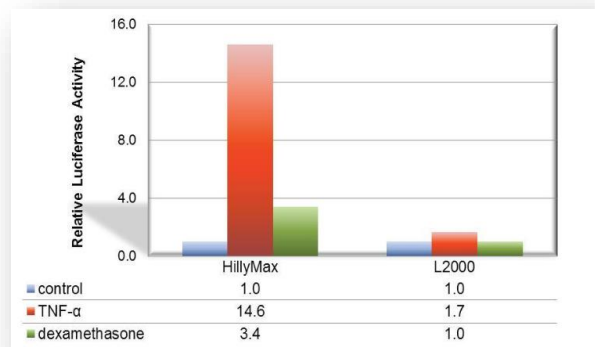
Cell	S2(Schneider 2) cell, 200,000 cells/well
Media	Schneider's Drosophila medium with 10% FCS
Antibiotics	50 units Penicillin/ml, 20 µg Streptomycin/ml
Microplate	24-well plate

Transfection Conditions

Vector	1 µg/well <pAct Gal4(6 kb), pUAS-mCD8::GFP(10 kb)>
Reagent	5 µl/well <HilyMax or Cellfectin>
	Medium was changed in 4 hours after transfection.
	Schneider's Drosophila medium (without serum and antibiotics) was used for the complex preparation.

*Data was kindly provided by Dr. Takashi Suzuki at Max Planck Institute of Neurobiology.

CELL SIGNALING RESEARCH



Signal transduction from A549 cell was examined with the TNF-α stimulation. For detection of cellular response, IL-8 dependent luciferase expression vector was transfected with HilyMax or Lipo2000. Good signal response was observed at HilyMax experiment.

IMPOTANT NOTES

To achieve good transfection efficiency, **optimization of cell density, DNA amount, HilyMax reagent amount are required.** Please try to find the best conditions by following DOJINDO' general protocol or optimized protocols.

Once Lipoform buffer is added to HilyMax reagent, store the mixture at -20°C for longer storage (Max 6 months). Mix the mixture well by vortex or pipetting before use. For frequent use, you can store the mixture at 0-5°C for 1 month.



EUROPEAN HEADQUARTERS
 DOJINDO EU GMBH
 Leopoldstr. 254, 80807 Munich, Germany
 Phone +49 89 3540-4806
 Fax +49 89 3540-4806
 email info@dojindo.eu.com
 www.dojindo.eu.com

DISTRIBUTED BY