



Another Successful Block

siLentFect™ Lipid Reagent for RNAi



Effective gene silencing using low amounts of siRNA and lipid.

siLentFect — the first step toward successful gene silencing

Bio-Rad's siLentFect lipid reagent was developed specifically for the delivery of siRNA to cultured mammalian cells. Its high affinity for siRNA and exceptional delivery characteristics will help you achieve robust silencing of your target gene with low lipid volumes and low concentrations of siRNA.

Key Features and Benefits

- **Effective transfection of siRNA to a broad range of cell lines** — use one lipid reagent for effective knockdown in all of your cell lines
- **Low volumes of lipid required per transfection** — minimize cell stress while saving money on each transfection
- **Use 5 nM or less siRNA for effective knockdown** — less siRNA reduces the risk of off-target effects
- **Low cytotoxicity** — achieve robust knockdown while maintaining culture health

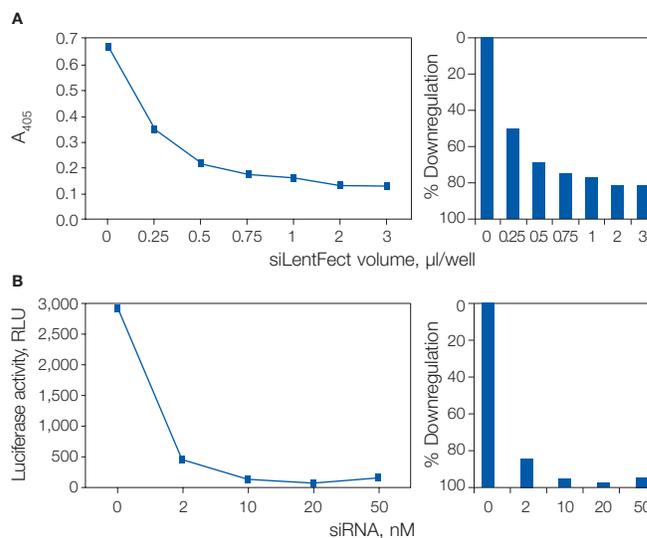
Examples of Successfully Transfected Cell Lines

Cell Line	Cell Type
A549	Human lung carcinoma
C166-GFP	Mouse yolk sac endothelial
Caco-2	Human colon adenocarcinoma
CHO-K1	Hamster ovary epithelial
COS-7	SV40-transformed African green monkey kidney
HEK 293	Human kidney epithelial
HeLa	Human cervical adenocarcinoma
HepG2	Human liver carcinoma
HUVEC	Human umbilical vein endothelial
Jurkat	Human T lymphoblast
LNCap	Human epithelial prostate carcinoma
MCF-7	Human breast adenocarcinoma
NIH-3T3	Mouse embryo fibroblast
Primary fibroblast	Human foreskin fibroblast
Primary keratinocyte	Human keratinocytes
SCC12B2	Human squamous cell carcinoma

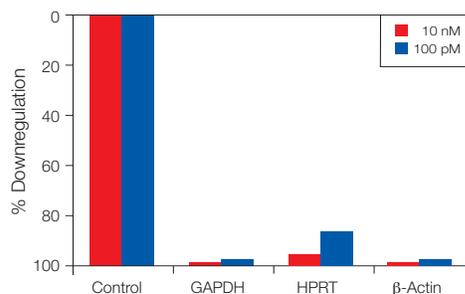
Ordering Information

Catalog #	Description
Reagents for siRNA Delivery	
170-3360	siLentFect Lipid Reagent for RNAi, 0.5 ml
170-3361	siLentFect Lipid Reagent for RNAi, 1.0 ml
170-3362	siLentFect Lipid Reagent for RNAi, 5 x 1.0 ml
Reagents for Plasmid DNA Delivery	
170-3350	TransFectin™ Lipid Reagent, 0.5 ml
170-3351	TransFectin Lipid Reagent, 1.0 ml
170-3352	TransFectin Lipid Reagent, 5 x 1.0 ml

SYBR is a trademark of Invitrogen Corp.



siLentFect mediates efficient transfection with low volumes of lipid and low concentrations of siRNA. A, CHO cells stably expressing the *lacZ* gene were grown in 24-well plates and transfected with an siRNA targeting *lacZ*. Cells were lysed and assayed for β -galactosidase activity 24 hr posttransfection. A significant reduction in expression of the *lacZ* gene product was observed even with low volumes of siLentFect. B, CHO cells stably expressing the luciferase gene were transfected in 96-well plates using 0.3 μ l siLentFect and a 21-mer anti-luciferase siRNA. The addition of siRNA at concentrations >10 nM did not produce any further significant reduction in luciferase activity.



siLentFect allows effective silencing with low concentrations of siRNA. HeLa cells were transfected with siLentFect and either 10 nM or 100 pM of a 27-mer siRNA targeted against three different endogenous genes. RT-qPCR was used 24 hr posttransfection to measure expression relative to a nonsilencing control. For real-time detection, total RNA was isolated using the Aurum™ total RNA mini kit and analyzed by the Experion™ automated electrophoresis system. For RT-qPCR, cDNA was synthesized using the iScript™ cDNA synthesis kit and amplified with iQ™ SYBR® Green supermix and the iCycler iQ® real-time detection system.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 2527 3454 Canada (905) 712-2771 China (86 21) 6426 0808 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-0 Greece 30 210 777 4396 Hong Kong (852) 2789 3300 Hungary 36 1 455 8800 India (91-124)-2398112/3/4, 5018111, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 331 99 99 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-64153188 South Africa 00 27 11 4428508 Spain 34 91 590 52 00 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan (886 2) 2578 7189/2578 7241 United Kingdom 020 8328 2000