

Transfection Reagent References

siLentFect™ Lipid Reagent for RNAi

Cell Line	Reference
184htrt (human breast epithelial) cells	Troussard AA et al., Preferential dependence of breast cancer cells versus normal cells on integrin-linked kinase for protein kinase B/Akt activation and cell survival, <i>Cancer Res</i> 66, 393–403 (2006)
293T cells	Goto Y et al., A new melanoma antigen fatty acid-binding protein 7, involved in proliferation and invasion, is a potential target for immunotherapy and molecular target therapy, <i>Cancer Res</i> 66, 4443–4449 (2006)
293T cells	Kostenko EV et al., Ccpg1, a novel scaffold protein that regulates the activity of the Rho guanine nucleotide exchange factor Dbs, <i>Mol Cell Biol</i> 26, 8964–8975 (2006)
2FTGH cells	Komuro A and Horvath CM, RNA and virus-independent inhibition of antiviral signaling by RNA helicase LGP2, <i>J Virol</i> 80, 12332–12342 (2006)
2FTGH cells	Ulane CM et al., Composition and assembly of STAT-targeting ubiquitin ligase complexes: paramyxovirus V protein carboxyl terminus is an oligomerization domain, <i>J Virol</i> 79, 10180–10189 (2005)
4T1 cells	Gonzalez RR et al., Leptin signaling promotes the growth of mammary tumors and increases the expression of vascular endothelial growth factor (VEGF) and its receptor type two (VEGF-R2), <i>J Biol Chem</i> 281, 26320–26328 (2006)
4T1 cells	Nam JS et al., Bone sialoprotein mediates the tumor cell-targeted prometastatic activity of transforming growth factor β in a mouse model of breast cancer, <i>Cancer Res</i> 66, 6327–6335 (2006)
A549 cells	Bartz SR et al., Small interfering RNA screens reveal enhanced cisplatin cytotoxicity in tumor cells having both BRCA network and TP53 disruptions, <i>Mol Cell Biol</i> 26, 9377–9386 (2006)
A549 cells	Dohadwala M et al., Cyclooxygenase-2-dependent regulation of E-cadherin: prostaglandin E ₂ induces transcriptional repressors ZEB1 and snail in non-small cell lung cancer, <i>Cancer Res</i> 66, 5338–5345 (2006)
Caco-2 BBe cells	Clayburgh DR et al., A differentiation-dependent splice variant of myosin light chain kinase, MLCK1, regulates epithelial tight junction permeability, <i>J Biol Chem</i> 279, 55506–55513 (2004)
Caco-2 cells	Hu Z et al., MAPKAPK-2 is a critical signaling intermediate in NHE3 activation following Na ⁺ -glucose cotransport, <i>J Biol Chem</i> 281, 24247–24253 (2006)
COS-7 cells	Evans K et al., Interaction of two hereditary spastic paraplegia gene products, spastin and atlastin, suggests a common pathway for axonal maintenance, <i>PNAS</i> 103, 10666–10671 (2006)
DU145 cells	Li X et al., Endogenous inhibition of histone deacetylase 1 by tumor-suppressive maspin, <i>Cancer Res</i> 66, 9323–9329 (2006)
F9 cells	Gao Y and Wang H, Casein kinase 2 is activated and essential for Wnt/ β -catenin signaling, <i>J Biol Chem</i> 281, 18394–18400 (2006)
HCT-15 cells	Ludwig JA et al., Selective toxicity of NSC73306 in MDR1-positive cells as a new strategy to circumvent multidrug resistance in cancer, <i>Cancer Res</i> 66, 4808–4815 (2006)
HCT 116 cells	Wang Y and Li G, ING3 promotes UV-induced apoptosis via Fas/caspase-8 pathway in melanoma cells, <i>J Biol Chem</i> 281, 11887–11893 (2006)
HEK 293 cells	Green RA et al., APC and EB1 function together in mitosis to regulate spindle dynamics and chromosome alignment, <i>Mol Biol Cell</i> 16, 4609–4622 (2005)
HEK 293T cells	Komuro A and Horvath CM, RNA and virus-independent inhibition of antiviral signaling by RNA helicase LGP2, <i>J Virol</i> 80, 12332–12342 (2006)
HeLa cells	Evans K et al., Interaction of two hereditary spastic paraplegia gene products, spastin and atlastin, suggests a common pathway for axonal maintenance, <i>PNAS</i> 103, 10666–10671 (2006)
HeLa cells	Komuro A and Horvath CM, RNA and virus-independent inhibition of antiviral signaling by RNA helicase LGP2, <i>J Virol</i> 80, 12332–12342 (2006)
HeLa cells	Olabisi OO et al., Bcr interacts with components of the endosomal sorting complex required for transport-I and is required for epidermal growth factor receptor turnover, <i>Cancer Res</i> 66, 6250–6257 (2006)
HepG2 cells	Simonsson M et al., The balance between acetylation and deacetylation controls Smad7 stability, <i>J Biol Chem</i> 280, 21797–21803 (2005)
Huh-7 cells	Ishii N et al., Diverse effects of cyclosporine on hepatitis C virus strain replication, <i>J Virol</i> 80, 4510–4520 (2006)
HUVECs (human umbilical vein endothelial cells)	Hattori Y et al., Metformin inhibits cytokine-induced nuclear factor κ B activation via AMP-activated protein kinase activation in vascular endothelial cells, <i>Hypertension</i> 47, 1183–1188 (2006)

(continues)

siLentFect Lipid Reagent for RNAi (cont.)

Cell Line	Reference
LNCaP cells	Lee DY et al., Histone H3 lysine 9 methyltransferase G9a is a transcriptional coactivator for nuclear receptors, <i>J Biol Chem</i> 281, 8476–8485 (2006)
LNCap-HP cells	Teyssier C et al., Transcriptional intermediary factor 1 α mediates physical interaction and functional synergy between the coactivator-associated arginine methyltransferase 1 and glucocorticoid receptor-interacting protein 1 nuclear receptor coactivators, <i>Mol Endocrinol</i> 20, 1276–1286 (2006)
MCF-7 cells	Schultz-Norton JR et al., Protein disulfide isomerase serves as a molecular chaperone to maintain estrogen receptor α structure and function, <i>Mol Endocrinol</i> 20, 1982–1995 (2006)
MDA-MB-231 (human breast cancer) cells	Troussard AA et al., Preferential dependence of breast cancer cells versus normal cells on integrin-linked kinase for protein kinase B/Akt activation and cell survival, <i>Cancer Res</i> 66, 393–403 (2006)
Melanoma cells	Goto Y et al., A new melanoma antigen fatty acid-binding protein 7, involved in proliferation and invasion, is a potential target for immunotherapy and molecular target therapy, <i>Cancer Res</i> 66, 4443–4449 (2006)
MEWO cells	Wang Y and Li G, ING3 promotes UV-induced apoptosis via Fas/caspase-8 pathway in melanoma cells, <i>J Biol Chem</i> 281, 11887–11893 (2006)
MMRU cells	Wang J et al., The novel tumor suppressor p33ING2 enhances nucleotide excision repair via induction of histone H4 acetylation and chromatin relaxation, <i>Cancer Res</i> 66, 1906–1911 (2006)
MMRU cells	Wang Y and Li G, ING3 promotes UV-induced apoptosis via Fas/caspase-8 pathway in melanoma cells, <i>J Biol Chem</i> 281, 11887–11893 (2006)
Mouse endothelial cells (murine EC)	Isakson BE et al., Heterocellular contact at the myoendothelial junction influences gap junction organization, <i>Circ Res</i> 97, 44–51 (2005)
NIH 3T3 cells	Liu Z et al., Transformation by the Rho-specific guanine nucleotide exchange factor Dbs requires ROCK I-mediated phosphorylation of myosin light chain, <i>J Biol Chem</i> 281, 16043–16051 (2006)
PC-3 cells	Li X et al., Endogenous inhibition of histone deacetylase 1 by tumor-suppressive maspin, <i>Cancer Res</i> 66, 9323–9329 (2006)
PC-3 cells	Yin S et al., Tumor suppressive maspin regulates cell response to oxidative stress by direct interaction with glutathione S-transferase, <i>J Biol Chem</i> 280, 34985–34996 (2005)
Primary HTM (human trabecular meshwork) cells	Husain S et al., Mechanisms linking adenosine A ₁ receptors and ERK1/2 activation in human trabecular meshwork cells, <i>J Pharmacol Exp Ther</i> , 10.1124/jpet.106.110981 (2006)
Primary ovarian epithelial cells	Cheung LWT et al., Pigment epithelium-derived factor is estrogen sensitive and inhibits the growth of human ovarian cancer and ovarian surface epithelial cells, <i>Endocrinology</i> 147, 4179–4191 (2006)
ST2 stromal cells	Deregowski V et al., Notch 1 overexpression inhibits osteoblastogenesis by suppressing Wnt/ β -catenin but not bone morphogenetic protein signaling, <i>J Biol Chem</i> 281, 6203–6210 (2006)
SVEC4 cells	Hattori Y et al., Metformin inhibits cytokine-induced nuclear factor κ B activation via AMP-activated protein kinase activation in vascular endothelial cells, <i>Hypertension</i> 47, 1183–1188 (2006)
TOV21G cells	Bartz SR et al., Small interfering RNA screens reveal enhanced cisplatin cytotoxicity in tumor cells having both BRCA network and TP53 disruptions, <i>Mol Cell Biol</i> 26, 9377–9386 (2006)
Vascular smooth muscle cells (VSMC)	Isakson BE et al., Heterocellular contact at the myoendothelial junction influences gap junction organization, <i>Circ Res</i> 97, 44–51 (2005)
YAMC cells	Tao Y et al., Soluble factors from <i>Lactobacillus</i> GG activate MAPKs and induce cytoprotective heat shock proteins in intestinal epithelial cells, <i>Am J Physiol Cell Physiol</i> 290, C1018–C1030 (2006)

TransFectin™ Reagent

Cell Line	Reference
293 cells	Nakerakanti SS et al., Fli1 and Ets1 have distinct roles in connective tissue growth factor/CCN2 gene regulation and induction of the profibrotic gene program, <i>J Biol Chem</i> 281, 25259–25269 (2006)
293A cells	Lerner-Marmarosh N et al., Human biliverdin reductase: a member of the insulin receptor substrate family with serine/threonine/tyrosine kinase activity, <i>Proc Natl Acad Sci USA</i> 102, 7109–7114 (2005)
293A cells	Miralem T et al., Small interference RNA-mediated gene silencing of human biliverdin reductase, but not that of heme oxygenase-1, attenuates arsenite-mediated induction of the oxygenase and increases apoptosis in 293A kidney cells, <i>J Biol Chem</i> 280, 17084–17092 (2005)
293FT cells	Gööz M et al., 5-HT _{2A} receptor induces ERK phosphorylation and proliferation through ADAM-17 tumor necrosis factor- α -converting enzyme (TACE) activation and heparin-bound epidermal growth factor-like growth factor (HB-EGF) shedding in mesangial cells, <i>J Biol Chem</i> 281, 21004–21012 (2006)
293T cells	Billelo JP et al., A genetic system for rhesus monkey rhadinovirus: use of recombinant virus to quantitate antibody-mediated neutralization, <i>J Virol</i> 80, 1549–1562 (2006)
293T cells	Sakaguchi T et al., AIP1/Alix is a binding partner of Sendai virus C protein and facilitates virus budding, <i>J Virol</i> 79, 8933–8941 (2005)
293T cells	Yuste E et al., Simian immunodeficiency virus engrafted with human immunodeficiency virus type 1 (HIV-1)-specific epitopes: replication, neutralization, and survey of HIV-1-positive plasma, <i>J Virol</i> 80, 3030–3041 (2006)
A375 cells	Reu FJ et al., Overcoming resistance to interferon-induced apoptosis of renal carcinoma and melanoma cells by DNA demethylation, <i>J Clin Oncol</i> 24, 3771–3779 (2006)
ACHN cells	Reu FJ et al., Overcoming resistance to interferon-induced apoptosis of renal carcinoma and melanoma cells by DNA demethylation, <i>J Clin Oncol</i> 24, 3771–3779 (2006)

(continues)

TransFectin Reagent (cont.)

Cell Line	Reference
BC3 cells	Keller SA et al., NF- κ B is essential for progression of KSHV- and EBV-infected lymphomas in vivo, <i>Blood</i> 107, 3295–3302 (2006)
C2C12 cells	Mortusewicz O et al., Differential recruitment of DNA ligase I and III to DNA repair sites, <i>Nucleic Acids Res</i> 34, 3523–3532 (2006)
C2C12 cells	Mortusewicz O et al., Recruitment of DNA methyltransferase I to DNA repair sites, <i>Proc Natl Acad Sci USA</i> 102, 8905–8909 (2005)
C2C12 myoblasts (mouse)	Perucca P et al., Spatiotemporal dynamics of p21 ^{CDKN1A} protein recruitment to DNA-damage sites and interaction with proliferating cell nuclear antigen, <i>J Cell Sci</i> 119, 1517–1527 (2006)
C6 cells	De Vuyst E et al., Connexin hemichannels and gap junction channels are differentially influenced by lipopolysaccharide and bFGF, <i>Mol Biol Cell</i> , 10.1091/mbc.E06-03-0182 (2006)
CHP-212 cells	Zhong N et al., DJ-1 transcriptionally up-regulates the human tyrosine hydroxylase by inhibiting the sumoylation of pyrimidine tract-binding protein-associated splicing factor, <i>J Biol Chem</i> 281, 20940–20948 (2006)
COS-7 cells	Colletti KS et al., Human cytomegalovirus UL84 oligomerization and heterodimerization domains act as transdominant inhibitors of <i>ori</i> Lyt-dependent DNA replication: evidence that IE2-UL84 and UL84-UL84 interactions are required for lytic DNA replication, <i>J Virol</i> 78, 9203–9214 (2004)
CV-1 cells	Chen Y et al., Unraveling protein-protein interactions in living cells with fluorescence fluctuation brightness analysis, <i>Biophys J</i> 88, 4366–4377 (2005)
CV-1 cells	Sanchez-Andres A et al., Molecular brightness determined from a generalized form of Mandel's Q-parameter, <i>Biophys J</i> 89, 3531–3547 (2005)
HEK 293 cells	Derler I et al., Dynamic but not constitutive association of calmodulin with rat TRPV6 channels enables fine tuning of Ca ²⁺ -dependent inactivation, <i>J Physiol</i> 577, 31–44 (2006)
HEK 293 cells	De Vuyst E et al., Connexin hemichannels and gap junction channels are differentially influenced by lipopolysaccharide and bFGF, <i>Mol Biol Cell</i> , 10.1091/mbc.E06-03-0182 (2006)
HEK 293 cells	Gavin MA et al., Single-cell analysis of normal and FOXP3-mutant human T cells: FOXP3 expression without regulatory T cell development, <i>PNAS</i> 103, 6659–6664 (2006)
HEK 293 cells	Zhang C et al., Toll-like receptor 2 mediates alveolar macrophage response to <i>Pneumocystis murina</i> , <i>Infect Immun</i> 74, 1857–1864 (2006)
HEK 293T cells	An HJ et al., Activation of Ras up-regulates pro-apoptotic BNIP3 in nitric oxide-induced cell death, <i>J Biol Chem</i> 281, 33939–33948 (2006)
HEK 293T cells	Brown L et al., In vitro analysis of partial loss-of-function ZIC2 mutations in holoprosencephaly: alanine tract expansion modulates DNA binding and transactivation, <i>Hum Mol Genet</i> 14, 411–420 (2005)
HEK 293T cells	Pohjoismäki JLO et al., Alterations to the expression level of mitochondrial transcription factor A, TFAM, modify the mode of mitochondrial DNA replication in cultured human cells, <i>Nucleic Acids Res</i> 34, 5815–5828 (2006)
HeLa cells	De Vuyst E et al., Connexin hemichannels and gap junction channels are differentially influenced by lipopolysaccharide and bFGF, <i>Mol Biol Cell</i> , 10.1091/mbc.E06-03-0182 (2006)
HeLa cells	Mortusewicz O et al., Differential recruitment of DNA ligase I and III to DNA repair sites, <i>Nucleic Acids Res</i> 34, 3523–3532 (2006)
HeLa cells	Mortusewicz O et al., Recruitment of DNA methyltransferase I to DNA repair sites, <i>Proc Natl Acad Sci USA</i> 102, 8905–8909 (2005)
HepG2 cells	Kern MA et al., Cyclooxygenase-2 inhibition induces apoptosis signaling via death receptors and mitochondria in hepatocellular carcinoma, <i>Cancer Res</i> 66, 7059–7066 (2006)
Huh-7 cells	Kern MA et al., Cyclooxygenase-2 inhibition induces apoptosis signaling via death receptors and mitochondria in hepatocellular carcinoma, <i>Cancer Res</i> 66, 7059–7066 (2006)
InR1G9 cells	Mamin A and Philippe J, Activin A decreases <i>glucagon</i> and <i>arx</i> gene expression in α -cell lines, <i>Mol Endocrinol</i> , 10.1210/me.2005-0530 (2006)
Jurkat cells	Niinuma A et al., Aberrant activation of the interleukin-2 autocrine loop through the nuclear factor of activated T cells by nonleukemogenic human T-cell leukemia virus type 2 but not by leukemogenic type 1 virus, <i>J Virol</i> 79, 11925–11934 (2005)
Jurkat cells	Satou Y et al., HTLV-I basic leucine zipper factor gene mRNA supports proliferation of adult T cell leukemia cells, <i>Proc Natl Acad Sci USA</i> 103, 720–725 (2006)
LLC-MK2 cells	Cseke G et al., Human metapneumovirus fusion protein vaccines that are immunogenic and protective in cotton rats, <i>J Virol</i> , 10.1128/JVI.00844-06 (2006)
LNCaP _{nan} or LNCaP-HP cells	Unni E et al., Changes in androgen receptor nongenotropic signaling correlate with transition of LNCaP cells to androgen independence, <i>Cancer Res</i> 64, 7156–7168 (2004)
MEF (mouse embryonic fibroblast) cells	Alirol E et al., The mitochondrial fission protein hFis1 requires the endoplasmic reticulum gateway to induce apoptosis, <i>Mol Biol Cell</i> 17, 4593–4605 (2006)
MEF (mouse embryonic fibroblast) cells	Goulimari P et al., α 12/13 is essential for directed cell migration and localized Rho-Dia1 function, <i>J Biol Chem</i> 280, 42242–42251 (2005)
MIN6 β cells	Oh E and Thurmond D, The stimulus-induced tyrosine phosphorylation of Munc18c facilitates vesicle exocytosis, <i>J Biol Chem</i> 281, 17624–17634 (2006)
NIH 3T3 cells	Evans KJ et al., Linking axonal degeneration to microtubule remodeling by Spastin-mediated microtubule severing, <i>J Cell Biol</i> 168, 599–606 (2005)
NIH 3T3 cells	Haritunians T et al., Functional analysis of a recurrent missense mutation in Notch3 in CADASIL, <i>J Neurol Neurosurg Psychiatry</i> 76, 1242–1248 (2005)

(continues)

TransFectin Reagent (cont.)

Cell Line	Reference
Primary human corneal fibroblasts	Lu Y et al., Inhibitory effect of triptolide on chemokine expression induced by proinflammatory cytokines in human corneal fibroblasts, <i>Invest Ophthalmol Vis Sci</i> 46, 2346–2352 (2005)
RAW264.7 macrophage cells	An HJ et al., Activation of Ras up-regulates pro-apoptotic BNP3 in nitric oxide-induced cell death, <i>J Biol Chem</i> 281, 33939–33948 (2006)
S2 cells (<i>Drosophila</i>)	Paschinger K et al., Molecular basis of anti-horseradish peroxidase staining in <i>Caenorhabditis elegans</i> , <i>J Biol Chem</i> 279, 49588–49598 (2004)
S2 cells (<i>Drosophila</i>)	Rendic D et al., Modulation of neural carbohydrate epitope expression in <i>Drosophila melanogaster</i> cells, <i>J Biol Chem</i> 281, 3343–3353 (2006)
SH-SY5Y cells	Xu J et al., The Parkinson's disease-associated DJ-1 protein is a transcriptional co-activator that protects against neuronal apoptosis, <i>Hum Mol Genet</i> 14, 1231–1241 (2005)
SH-SY5Y cells	Zhong N et al., DJ-1 transcriptionally up-regulates the human tyrosine hydroxylase by inhibiting the sumoylation of pyrimidine tract-binding protein-associated splicing factor, <i>J Biol Chem</i> 281, 20940–20948 (2006)
SW480 (colon cancer) cells	Beck SE et al., Bone morphogenetic protein signaling and growth suppression in colon cancer, <i>Am J Physiol Gastrointest Liver Physiol</i> 291, G135–G145 (2006)
TC7	Evans KJ et al., Linking axonal degeneration to microtubule remodeling by Spastin-mediated microtubule severing, <i>J Cell Biol</i> 168, 599–606 (2005)
THP-1 monocyte cells	Munteanu A et al., Antagonistic effects of oxidized low density lipoprotein and α -tocopherol on CD36 scavenger receptor expression in monocytes: involvement of protein kinase B and peroxisome proliferator-activated receptor- γ , <i>J Biol Chem</i> 10, 6489–6497 (2006)
U-2 OS cells	Giallourakis C et al., A molecular-properties-based approach to understanding PDZ domain proteins and PDZ ligands, <i>Genome Res</i> 16, 1056–1072 (2006)

For direct link to abstracts, go to www.bio-rad.com/transfectionreferences/

Last update: November 2006



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **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 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000