

## Gene Pulser® siRNA Electroporation References

Cell Line	Reference
33B (rat brain tumor-derived cell line)	Hemmings-Mieszczak M et al., Independent combinatorial effect of antisense oligonucleotides and RNAi-mediated specific inhibition of the recombinant rat P2X <sub>3</sub> receptor, <i>Nucleic Acids Res</i> 31, 2117–2126 (2003)
3T3-F442A (adipocytes)	Mitra P et al., RNAi-based analysis of CAP, Cbl, and Crkl function in the regulation of GLUT4 by insulin, <i>J Biol Chem</i> 279, 37431–37435 (2004)
3T3-L1 (adipocytes)	Jiang ZY et al., Insulin signaling through Akt/protein kinase B analyzed by small interfering RNA-mediated gene silencing, <i>Proc Natl Acad Sci USA</i> 100, 7569–7574 (2003)
3T3-L1 (adipocytes)	Guilherme A et al., Role of EHD1 and EHBP1 in perinuclear sorting and insulin-regulated GLUT4 recycling in 3T3-L1 adipocytes, <i>J Biol Chem</i> 279, 40062–40075 (2004)
3T3-L1 (adipocytes)	Hresko RC and Mueckler M, mTOR/RICTOR is the Ser473 kinase for Akt/PKB in 3T3-L1 adipocytes, <i>J Biol Chem</i> 10.1074/jbc.M508361200 (2005)
3T3-L1 (adipocytes)	Mitra P et al., RNAi-based analysis of CAP, Cbl, and Crkl function in the regulation of GLUT4 by insulin, <i>J Biol Chem</i> 279, 37431–37435 (2004)
3T3-L1 (fibroblasts)	Guilherme A et al., EHD2 and the novel EH domain binding protein EHBP1 couple endocytosis to the actin cytoskeleton, <i>J Biol Chem</i> 279, 10593–10605 (2004)
A7r5 (rat aortic smooth muscle)	Soboloff J et al., Role of endogenous TRPC6 channels in Ca <sup>2+</sup> signal generation in A7r5 smooth muscle cells, <i>J Biol Chem</i> 280, 39786–39794 (2005)
Ba/F3-β	Rasclé A and Lees E, Chromatin acetylation and remodeling at the <i>Cis</i> promoter during STAT5-induced transcription, <i>Nucleic Acids Res</i> 31, 6882–6890 (2003)
CHO-K1	Hemmings-Mieszczak M et al., Independent combinatorial effect of antisense oligonucleotides and RNAi-mediated specific inhibition of the recombinant rat P2X <sub>3</sub> receptor, <i>Nucleic Acids Res</i> 31, 2117–2126 (2003)
CML (chronic myeloid leukemia)	Aichberger K et al., Identification of <i>mcl-1</i> as a BCR/ABL-dependent target in chronic myeloid leukemia (CML): evidence for cooperative antileukemic effects of imatinib and <i>mcl-1</i> antisense oligonucleotides, <i>Blood</i> 105, 3303–3311 (2005)
CWR22-Rv1	Zha S et al., α-Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
DU145	Zha S et al., α-Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
E10 (immature mouse thymocyte)	McManus MT et al., Small interfering RNA-mediated gene silencing in T lymphocytes, <i>J Immunol</i> 169, 5754–5760 (2002)
E10 (immature mouse thymocyte)	McManus MT et al., Gene silencing using micro-RNA designed hairpins, <i>RNA</i> 8, 842–850 (2002)
FRTL-5 (thyroid cells)	Kim H et al., Thyrotropin-mediated repression of class II trans-activator expression in thyroid cells: involvement of STAT3 and suppressor of cytokine signaling, <i>J Immunol</i> 171, 616–627 (2003)
HaCaT	Seo M et al., Cdc42-dependent mediation of UV-induced p38 activation by G protein βγ subunits, <i>J Biol Chem</i> 279, 17366–17375 (2004)
HeLa	Quensel C et al., In vivo analysis of importin α proteins reveals cellular proliferation inhibition and substrate specificity, <i>Mol Cell Biol</i> 24, 10246–10255 (2004)
HeLa S3	Zha S et al., α-Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
HeLa SS6	Schaffert N et al., RNAi knockdown of hPrp31 leads to an accumulation of U4/U6 di-snRNPs in Cajal bodies, <i>EMBO J</i> 23, 3000–3009 (2004)
HL-60/Vinc (human acute myelogenous leukemia cells)	Zhong X and Safa AR, RNA helicase A in the MEF1 transcription factor complex up-regulates the MDR1 gene in multidrug-resistant cancer cells, <i>J Biol Chem</i> 279, 17134–17141 (2004)
Human precursor B cell (acute lymphoblastic leukemia)	Abrams MT et al., Inhibition of glucocorticoid-induced apoptosis by targeting the major splice variants of BIM mRNA with small interfering RNA and short hairpin RNA, <i>J Biol Chem</i> 279, 55809–55817 (2004)
Jurkat	Coburn GA and Cullen BR, Potent and specific inhibition of human immunodeficiency virus type 1 replication by RNA interference, <i>J Virol</i> 76, 9225–9231 (2002)
K562	Lunghi P et al., Treatment with arsenic trioxide (ATO) and MEK1 inhibitor activates the p73-p53AIP1 apoptotic pathway in leukemia cells, <i>Blood</i> 104, 519–525 (2004)

(continued)

Cell Line	Reference
K562	Liao B et al., Targeted knockdown of the RNA-binding protein CRD-BP promotes cell proliferation via an insulin-like growth factor II-dependent pathway in human K562 leukemia cells, <i>J Biol Chem</i> 279, 48716–48724 (2004)
Karpas 299	Bacchiocchi R et al., Activation of $\alpha$ -diacylglycerol kinase is critical for the mitogenic properties of anaplastic lymphoma kinase, <i>Blood</i> 106, 2175–2182 (2005)
L428 (Hodgkin/Reed-Sternberg)	Mathas S et al., c-FLIP mediates resistance of Hodgkin/Reed-Sternberg cells to death receptor-induced apoptosis, <i>J Exp Med</i> 199, 1041–1052 (2004)
L540Cy	Mathas S et al., c-FLIP mediates resistance of Hodgkin/Reed-Sternberg cells to death receptor-induced apoptosis, <i>J Exp Med</i> 199, 1041–1052 (2004)
LAPC-4	Zha S et al., $\alpha$ -Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
LNCaP	Zha S et al., $\alpha$ -Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
MCF-7	Yeh S et al., Abnormal mammary gland development and growth retardation in female mice and MCF7 breast cancer cells lacking androgen receptor, <i>J Exp Med</i> 198, 1899–1908 (2003)
MDCK (Madin-Darby canine kidney)	Ge Q et al., RNA interference of influenza virus production by directly targeting mRNA for degradation and indirectly inhibiting all viral RNA transcription, <i>Proc Natl Acad Sci USA</i> 100, 2718–2723 (2003)
NB4	Lunghi P et al., Treatment with arsenic trioxide (ATO) and MEK1 inhibitor activates the p73-p53AIP1 apoptotic pathway in leukemia cells, <i>Blood</i> 104, 519–525 (2004)
NB4	Chou WC et al., Role of NADPH oxidase in arsenic-induced reactive oxygen species formation and cytotoxicity in myeloid leukemia cells, <i>Proc Natl Acad Sci USA</i> 101, 4578–4583 (2004)
PC3	Zha S et al., $\alpha$ -Methylacyl-CoA racemase as an androgen-independent growth modifier in prostate cancer, <i>Cancer Res</i> 63, 7365–7376 (2003)
TF-1	Wen L et al., TL1A-induced NF- $\kappa$ B activation and c-IAP2 production prevent DR3-mediated apoptosis in TF-1 cells, <i>J Biol Chem</i> 278, 39251–39258 (2003)

For direct link to abstracts, go to [www.bio-rad.com/siRNAElectroporation/](http://www.bio-rad.com/siRNAElectroporation/)



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD **Australia** 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 4029300/5013478 **Israel** 03 963 6050  
**Italy** 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 55 5200 05 20 **The Netherlands** 0318 540666 **New Zealand** 64 9415 2280  
**Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 095 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 0861 246 723  
**Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189/2578 7241 **United Kingdom** 020 8328 2000