

RECOMMENDED READING

Accell™ siRNA References

Accell™ siRNA reagents are specially modified for use in difficult-to-transfect cells without the need for transfection reagents, virus, or electroporation. The following selected peer-reviewed publications have cited their successful use in a variety of experimental systems.

*For more references that use our siRNA for in vivo applications, please see our [in vivo siRNA reading list](#).

Established cell lines

ARPE-19 (human retinal epithelial cells) – 35
C1 tumor derived cells – 47
CD4+ primary human T cells – 4
CD14+ primary monocytes – 20
DG-75 human B lymphocytes – 71
GH3 (rat somatotroph pituitary cell line) – 56
HBECs (human bronchial epithelial cells) – 84
HCT-116 (colorectal carcinoma) – 25
HUVEC – 26
JJN3(plasma cell leukemia) – 40
Human lymphoblastoid cell lines – 82
LUHMES (Lund human mesencephalic cells) – 62
MEC1 (human chronic B cell leukemia) – 13
MN-1 – 32
MS1 (mouse pancreatic islet endothelia cells) – 21
NOD CD4+CD25+splenic cells – 37
NOXA – 40
OVCA 420 (ovarian carcinoma) – 52
PGA-1 (lymphocytic leukemia B cell line) – 61
RAW264.7 macrophages – 49
SHSY5Y (neuroblastoma) – 11, 23, 60, 85
SKBR3 (ER-/PR-/HER2+ breast cancer cell line) – 65
SNB19 glioma cells – 12
T98 glioma cells – 12
THP-1 monocytes – 10, 24, 42, 46, 57, 83
U266 (peripheral blood B lymphocyte myeloma) – 40
U937 – 46

Primary cells and in vivo

β-islet cells – 14
Bone marrow cells – 9, 16
Bronchial smooth muscle cells (BSMC) – 27, 28
Cardiomyocytes – 5
Cerebellar granule neurons (CGN) – 63
Colon stem/progenitor cells – 69
Corneal endothelial cells (adult human CECs), and ex vivo human corneal endothelium – 68
Cortical neurons – 1, 8, 41, 53, 66
Endometrial cells – 15
Endothelial cells – 7, 33
Extravillous trophoblasts (EVT) – 29
Fibroblasts (primary) – 66
Hepatocytes – 36, 38, 45
Human fetal liver cells – 87
Immortalized B cells – 59
Keratinocytes – 51
Lung epithelial cells – 73
Lymphocytes – 43
Macrophages – 3, 34, 48, 88
Mantle cell lymphoma cells (MCL) – 44
Mast cells – 92
Monocytes – 18
Mouse embryonic fibroblasts (MEF) – 22
Naïve fetal T cells – 75
Natural killer (NK) cell line – 55
Neonatal mouse ovary – 74
Neurons (primary rat) – 19
Neurons derived from iPS cells – 67
Oligodendrocyte precursors – 54
Peripheral blood mononuclear cells (PBMC) – 6, 31, 78
Primary human T-cells – 64, 86
Primary mouse T-cells – 70
Regulatory T cells – 76
Stem cell-derived peripheral neurons – 72

Vascular smooth muscle cells (VSMC) – 2, 58
 In vivo skin delivery – 17
 In vivo rat periodontal model – 30
 In vivo mouse intradermal injection – 39
 In vivo mouse model – 50, 77
 In vivo mouse brain – 62
 In vivo rat brain – 79, 80, 81, 89
 In vivo mouse, injection into tibialis anterior (TA) muscle – 90

References

1. A. H. Mortiboys, J. Aasly, et al. [Ursodeoxycholic acid rescues mitochondrial function in common forms of familial Parkinson's disease](#). *Brain* **136**, 3038–3050 (2013). [Mouse primary cortical neurons]
2. D. Gomez, K. Kessler, et al. [Modifications of Chromatin Dynamics Control Smad2 Pathway Activation in Aneurysmal Smooth Muscle Cells](#). *Circ. Res.* **113**, 881–890 (2013). [Human vascular smooth muscle cells (VSMC)]
3. S. Jiang, D. Part, et al. [Mitochondria and AMP-activated Protein Kinase-dependent Mechanism of Efferocytosis](#). *J. Biol. Chem.* **288**, 26013–26026 (2013). [mouse primary perinatal macrophages]
4. S. Sumitomo, K. Fujio, et al. [Transcription Factor Early Growth Response 3 Is Associated with the TGF-β1 Expression and the Regulatory Activity of CD4+Positive T Cells In vivo](#). *J. Immunol.* **191**, 2351–2359 (2013). [CD4+ primary human T cells]
5. F. Beig, J. Schmeckpeper, et al. [C3orf58, a Novel Paracrine Protein, Stimulates Cardiomyocyte Cell-Cycle Progression Through the PI3K-AKT-CDK7 Pathway](#). *Circ. Research* **113**, 372–380 (2013). [neonatal cardiomyocytes]
6. T. J. Mulrooney, P. E. Posch, et al. [DAP12 impacts trafficking and surface stability of killer immunoglobulin-like receptors on natural killer cells](#). *J. Leukoc. Biology* **94**, 301–313 (2013). [PMBCs]
7. R. Arita, S. Nakao, et al. [A Key Role for ROCK in TNF-α-Mediated Diabetic Microvascular Damage](#). *Invest. Ophthalmol. Vis. Science* **54**, 2373–2383 (2013). [cultured microvascular endothelial cells]
8. S. Hannila, M. Siddiq, et al. [Secretory Leukocyte Protease Inhibitor Reverses Inhibition by CNS Myelin, Promotes Regeneration in the Optic Nerve, and Suppresses Expression of the Transforming Growth Factor-β Signaling Protein Smad2](#). *J. Neuroscience* **33**, 5138–5151 (2013). [P6 CGNs and P1 rat cortical neurons]
9. S. Yasunaga, M. Ohtsubo et al. [Scmh1 Has E3 Ubiquitin Ligase Activity for Geminin and Histone H2A and Regulates Geminin Stability Directly or Indirectly via Transcriptional Repression of Hoxa9 and Hoxb4](#). *Mol. Cell. Biology* **33**, 644–660 (2013). [mouse BM cells]
10. G. Lopez-Castejon, N. M. Luheshi, et al. [Deubiquitinases Regulate the Activity of Caspase-1 and Interleukin-1β Secretion via Assembly of the Inflamasome](#). *J. Biol. Chem.* **288**, 2721–2733 (2013). [THP-1]
11. G. R. Tundo, D. Sbardella, et al. [Insulin-degrading Enzyme \(IDE\): A NOVEL HEAT SHOCK-LIKE PROTEIN](#). *J. Biol. Chemistry* **288**, 2281–2289 (2013). [SHSY5Y cells]
12. H. J. Seol, J. H. Chang, et al. [Overexpression of CD99 Increases the Migration and Invasiveness of Human Malignant Glioma Cells](#). *Genes & Cancer* **3**, 535–549 (2012). [T98 and SNB19 glioma cells]
13. A. G. Ramsay, A. J. Clear, et al. [Multiple inhibitory ligands induce impaired T-cell immunologic synapse function in chronic lymphocytic leukemia that can be blocked with lenalidomide: establishing a reversible immune evasion mechanism in human cancer](#). *Blood* **120**, 1412–1421 (2012). [MEC1 cells – Human chronic B cell leukemia]
14. M.R. Metukuri, P. Zhang, et al. [ChREBP Mediates Glucose-Stimulated Pancreatic β-Cell Proliferation](#). *Diabetes* **61**, 2004–2015 (2012). [human and rat islet cells]
15. E. M. De La Garza, P. A. Binkley, et al. [Raf-1, a Potential Therapeutic Target, Mediates Early Steps in Endometriosis Lesion Development by Endometrial Epithelial and Stromal Cells](#). *Endocrinology* **153**, 3911–3921 (2012). [primary endometrial cells]
16. X. Qu, G. Zhuang, et al. [Induction of Bv8 Expression by Granulocyte Colony-stimulating Factor in CD11b+Gr1+ Cells](#). *J. Biol. Chemistry* **287**, 19574–19584 (2012). [primary mouse bone marrow cells]
17. R. P. Hickerson, M. A. Flores, et al. [Use of self-delivery siRNAs to inhibit gene expression in an organotypic pachyonychia congenita model](#). *J Invest Dermatol* **131**(5), 1037–44 (2011) [In vivo skin delivery]
18. A. Nijnik, J. Pistolic, et al. [The role of the Src family kinase Lyn in the immunomodulatory activities of cathelicidin peptide LL-37 on monocytic cells](#). *J. Leukoc. Biology* **91**, 599–607 (2012) [monocytes]
19. A. Vagnoni, M. S. Perkinton, et al. [Calsyntenin-1 mediates axonal transport of the amyloid precursor protein and regulates Aβ production](#). *Human Molecular Genetics* **21**(13), 2845–54 (2012). [Rat E18 primary neurons]
20. P. Kuo, M. Huang, et al. [Lung Cancer-derived Galectin-1 Enhances Tumorigenic Potentiation of Tumor-associated Dendritic Cells by Expressing Heparin-binding EGF-like Growth Factor](#). *J. Biol. Chem.* **287**, 9753–9764 (2012). [CD14+ primary monocytes]
21. M. Franco, P. Roswall, et al. [Pericytes promote endothelial cell survival through induction of autocrine VEGF-A signaling and Bcl-w expression](#). *Blood* **118**, 2906–2917 (2011). [MS1cells; Mouse pancreatic islet endothelia]
22. C. X. George and C. E. Samuel, [Host Response to Polyomavirus Infection Is Modulated by RNA Adenosine Deaminase ADAR1 but Not by ADAR2](#). *Journal of Virology* **85**, 8338–8347 (2011). [MEF cells]
23. M. Meguro-Horiike, D. H. Yasui, et al. [Neuron-specific impairment of inter-chromosomal pairing and transcription in a novel model of human 15q-duplication syndrome](#). *Human Molecular Genetics* **20**(19), 3798–810 (2011). [SH-SY5Y]
24. Emily Turner-Brannen, Ka-Yee Grace Choi, et al. [Inflammatory Cytokines IL-32 and IL-17 Have Common Signaling Intermediates despite Differential Dependence on TNF-Receptor 1](#). *J. Immunology* **186**, 7127–7135 (2011). [THP-1 monocytes]
25. R. J. Boohaker, G. Zhang, et al. [BAX supports the mitochondrial network, promoting bioenergetics in nonapoptotic cells](#). *Am J Physiol Cell Physiology* **300**, C146–C1478 (2011). [HCT-116]
26. K. He, G. Sui, et al. [Feedback Regulation of Endothelial Cell Surface Plasmin Generation by PKC-dependent Phosphorylation of Annexin A2](#). *J. Biol. Chemistry* **286**, 15428–15439 (2011). [HUVEC cells]
27. P. L. Kuo, M. S. Huang, et al. [Signaling pathway of isophorone diisocyanate-responsive interleukin-8 in airway smooth muscle cells](#). *Eur. Respir. J* **37**, 1226–1236 (2011). [BSMCs]
28. P. Kuo, Y. Hsu, et al. [Bronchial Epithelium-Derived IL-8 and RANTES Increased Bronchial Smooth Muscle Cell Migration and Proliferation by Krüppel-like Factor 5 in Areca Nut-Mediated Airway Remodeling](#). *Toxicol. Sci.* **121**, 177–190 (2011). [BSMCs]
29. K. Biadasiewicz, S. Sonderegger, et al. [Transcription Factor AP-2α Promotes EGF-Dependent Invasion of Human Trophoblast](#). *Endocrinology* **152**, 1458–1469 (2011). [primary EVTs]
30. Q. Li, H. Yu, et al. [Silencing Mitogen-Activated Protein Kinase-Activated Protein Kinase-2 Arrests Inflammatory Bone Loss](#). *J. Pharmacol. Exp. Ther* **336**, 633–642 (2011). [ex vivo and In vivo rat periodontal model]
31. K. Billot, J. Soeur, et al. [Deregulation of Aiolos expression in chronic lymphocytic leukemia is associated with epigenetic modifications](#). *Blood* **117**, 1917–1927 (2011). [PBMCs]

32. L. Hubers, H. Valderrama-Carvajal, et al. [HuD interacts with survival motor neuron protein and can rescue spinal muscular atrophy-like neuronal defects](#). *Hum. Mol. Genetics* **20**, 553–579 (2011). [MN-1 cells]
33. S. F. Leicht, T. M. Schwarz, et al. [Adiponectin Pretreatment Counteracts the Detrimental Effect of a Diabetic Environment on Endothelial Progenitors](#). *Diabetes* **60**, 652–661 (2011). [endothelial colony-forming cells (ECFC)]
34. D. N. Petrusca, Y. Gu, et al. [Sphingolipid-mediated Inhibition of Apoptotic Cell Clearance by Alveolar Macrophages](#). *Journal of Biological Chemistry* **285**, 40322–40332 (2010). [Rat alveolar macrophages (AM)]
35. A. Giddabasappa, M. Bauler, et al. [17-β Estradiol Protects ARPE-19 Cells from Oxidative Stress through Estrogen Receptor-β](#). *Invest. Ophthalmol. Vis. Sci.* **51**, 5278–5287 (2010). [Human retinal epithelial cells ARPE-19 cells]
36. E. Murakami et. al. [Mechanism of Activation of PSI-7851 and its Diastereoisomer PSI-7977](#). *Journal of Biological Chemistry* **285**, 34337–34347 (2010). [primary human hepatocytes]
37. J. Zhang et al. [MEKK3 Overexpression Contributes to the Hyperresponsiveness of IL-12-Overproducing Cells and CD4+ T Conventional Cells in Nonobese Diabetic Mice](#). *Journal of Immunology* **185**, 3554–3563 (2010). [NOD CD4+CD25+ splenic cells]
38. M. Liao et al. [Inhibition of Hepatic Organic Anion-transporting Polypeptide by RNA Interference in Sandwich-cultured Human Hepatocytes: An in vitro Model to Assess Transporter-mediated Drug-drug Interactions](#). *Drug Metabolism and Disposition* **38**, 9 1612–1622 (2010). [freshly isolated human hepatocytes]
39. E. Gonzalez-Gonzalez et al. [Silencing of Reporter Gene Expression in Skin Using siRNAs and Expression of Plasmid DNA Delivered by a Soluble Protrusion Array Device \(PAD\)](#). *Molecular Therapy* **18**(9), 1667–74 (2010). [mouse intradermal injection]
40. B. Tunquist et al. [Mcl-1 Stability Determines Mitotic Cell Fate of Human Multiple Myeloma Tumor Cells Treated with the Kinesin Spindle Protein Inhibitor ARRY-520](#). *Molecular Cancer Therapeutics* **9**, 2046–2056 (2010). [multiple myeloma cell lines: JJN3 NOXA in RPMI 8226 and U266]
41. S. Suzuki et al. [Differential Roles of Epac in Regulating Cell Death in Neuronal and Myocardial Cells](#). *Journal of Biological Chemistry* **285**, 24248–24259 (2010). [primary mouse cortical neurons (E15–17)]
42. S. Winning et al. [Acute hypoxia induces HIF-independent monocyte adhesion to endothelial cells through increased intercellular adhesion molecule-1 expression: the role of hypoxic inhibition of prolyl hydroxylase activity for the induction of NF-κappa B](#). *Journal of Immunology* **185**, 1786–1793 (2010). [THP-1 monocytes]
43. M. Chetane et al. [Interleukin-7 mediates glucose utilization in lymphocytes through transcriptional regulation of the hexokinase II gene](#). *Am J Physiol Cell Physiol* **298**, C1560–C1571 (2010). [lymphocytes]
44. S. Desai et al. [PRDM1 Is Required for Mantle Cell Lymphoma Response to Bortezomib](#). *Molecular Cancer Research* **8**, 907–918 (2010). [mantle cell lymphoma (MCL) cells]
45. B. Mukhopadhyay et al. [Transcriptional Regulation of Cannabinoid Receptor-1 Expression in the Liver by Retinoic Acid Acting via Retinoic Acid Receptor](#). *Journal of Biological Chemistry* **285**, 19002–19011 (2010). [cultured mouse hepatocytes]
46. V. Saini et al. [CXC Chemokine Receptor 4 Is a Cell Surface Receptor for Extracellular Ubiquitin](#). *Journal of Biological Chemistry* **285**, 15566–15576 (2010). [THP-1 monocytes & U937]
47. I. Barbieri et al. [Constitutively Active Stat3 Enhances Neu-Mediated Migration and Metastasis in Mammary Tumors via Upregulation of Ctn](#). *Cancer Research* **70**, 2558–2567 (2010). [C1 tumor derived cells]
48. J. W. Perry et al. [Endocytosis of Murine Norovirus 1 into Murine Macrophages Is Dependent on Dynamin II and Cholesterol](#). *Journal of Virology* **84**, 6163–6176 (2010). [murine macrophages]
49. M. Steenport et al., [Matrix Metalloproteinase \(MMP\)-1 and MMP-3 Induce Macrophage MMP-9: Evidence for the Role of TNF-α and Cocl oxygenase-2](#). *Journal of Immunology* **183**(12), 8119–27 (2009). [RAW264.7 macrophages]
50. A. DiFeo et al. [KLF6-SV1 Is a Novel Antiapoptotic Protein That Targets the BH3-Only Protein NOXA for Degradation and Whose Inhibition Extends Survival in an Ovarian Cancer Model](#). *Cancer Research* **69**, 4733–41 (2009). [*In vivo* mouse model]
51. A. Kovalenko et al. [Caspase-8 Deficiency in Epidermal Keratinocytes Triggers an Inflammatory Skin Disease](#). *Journal of Experimental Medicine* **206**: 2161–2177 (2009). [Keratinocytes]
52. C. Bartholomeusz et al. [PEA-15 Induces Autophagy in Human Ovarian Cancer Cells and is Associated with Prolonged Overall Survival](#). *Cancer Research* **68**, 9302–9310 (2008). [OVCA 420; ovarian carcinoma]
53. A.M. Dolga et al. [TNF-alpha-mediates neuroprotection against glutamate-induced excitotoxicity via NF-κappaB-dependent up-regulation of K2.2 channels](#). *Journal of Neurochemistry* **107**, 1158–1167 (2008). [mouse primary cortical neurons]
54. F. Mir and G.C. Le Breton. [A Novel Nuclear Signaling Pathway for Thromboxane A2 Receptors in Oligodendrocytes: Evidence for Signaling Compartmentalization during Differentiation](#). *Mol. Cell. Biology* **28**, 6329–6341 (2008). [oligodendrocyte precursors]
55. C.B Lai, Y.Zhang, et al. [Creation of the two isoforms of rodent NKG2D was driven by a B1 retrotransposon insertion](#). *Nucleic Acids Research* **37**(9), 3032–43 (2009). [mouse NK cell line]
56. G. A. Peters et al. [The double-stranded RNA-binding protein, PACT, is required for postnatal anterior pituitary proliferation](#). *PNAS* **106**(26), 10696–10701 (2009). [GH3; rat somatotrophs (pituitary cell line) and LβT2 gonadotrophs]
57. N. Mookherjee et al. [Intracellular Receptor for Human Host Defense Peptide LL-37 in Monocytes](#). *J. Immunol.* 2009; 183. 2688–2696 [THP-1; human monocytes]
58. B. Zheng et. al. [Krüppel-like Factor 4 Inhibits Proliferation by Platelet-derived Growth Factor Receptor β-mediated, Not by Retinoic Acid Receptor α-mediated, Phosphatidylinositol 3-Kinase and ERK Signaling in Vascular Smooth Muscle Cells](#). *Journal of Biological Chemistry* **284** (34), 22773–22785 (2009). [VUMC; primary rat vascular smooth muscle cells]
59. D. Smirnov et. al. [Genetic Analysis of Radiation-induced Changes in Human Gene Expression](#). *Nature* **459**, 587–591 (2009). [immortalized B cells]
60. U. Dreses-Werringloer et. al. [A Polymorphism in CALHM1 Influences Ca²⁺ Homeostasis, Ab Levels, and Alzheimer's Disease Risk](#). *Cell* **133**, 1149–1161 (2008). [SHSY-5Y; human neuroblastoma]
61. A.M. McElligott et. al. [The Novel Tubulin-Targeting Agent Pyrrolo-1,5-Benzoxazepine-15 Induces Apoptosis in Poor Prognostic Subgroups of Chronic Lymphocytic Leukemia](#). *Cancer* **69**(21), 8366–75 (2009). [PGA-1; EBV-transformed chronic lymphocytic leukemia (CLL) B cell line]
62. H. Xu et. al. [Tau Silencing by siRNA in the P301S Mouse Model of Tauopathy](#). *Current Gene Therapy* **14**, 343–351 (2014) [*In vivo* mouse brain injection]
63. A. Collado-Alsina et. al. [The Regulation of Synaptic Vesicle Recycling by cGMP-Dependent Protein Kinase Type II in Cerebellar Granule Cells under Strong and Sustained Stimulation](#). *J. Neuroscience*. 2014; 34:8788–8799. [cerebellar granule cells]
64. M. Freeley et. al. [RNAi Screening with Self-Delivering Synthetic siRNAs for Identification of Genes That Regulate Primary Human T Cell Migration](#). *J. Biomolecular Screening*. 2015; DOI: 10.1177/1087057115588288 [Primary human T-cells]
65. S. Das, G. Sondarva, et. al. [Human Epidermal Growth Factor Receptor 2 \(HER2\) Impedes MLK3 Kinase Activity to Support Breast Cancer Cell Survival](#). *J. Biol. Chemistry* **290**, 35, 21705–21712 (2015). [SKBR3 (ER-/PR-/HER2+ breast cancer cell line)]

66. J. Bruban, G. Voloudakis, et. al. [Presenilin 1 is necessary for neuronal, but not glial, EGFR expression and neuroprotection via gamma-secretase-independent transcriptional mechanisms](#). *FASEB J.* **29**, 3702–3712 (2015) [murine primary cortical neurons (E15.5) and primary fibroblasts]
67. M. Komatsu, H. E. Wheeler, et. al. [Pharmacogenetic Discovery in Paclitaxel-Induced Sensory Peripheral Neuropathy](#). *Clin. Cancer Res.* **21**, 4337–4346 (2015) [Neurons derived from iPS cells]
68. J. G. Lee and M. Heur. [WNT10B Enhances Proliferation through β-Catenin and RAC1 GTPase in Human Corneal Endothelial Cells](#). *J. Biol. Chem.* **290**, 26752–26764 (2015) [Adult human corneal endothelial cells (CECs) and ex vivo human corneal endothelium]
69. I. Nasri, D. Bonnet, et. al. [PAR2-dependent activation of GSK3β regulates the survival of colon stem/progenitor cells](#). *Am J Physiol Gastrointest Liver Physiol* **311**, G221–G236 (2016) [colon stem/progenitor cells]
70. H. Lee, J.W. Lee, et. al. [Inflammation and Inflammatory Mediators in Kidney Disease: Cln 3-requiring 9 is a negative regulator of Th17 pathway-driven inflammation in anti-glomerular basement membrane glomerulonephritis](#). *Am J Physiol Renal Physiol.* **311**, F505–F519 (2016). [Primary mouse T-cells]
71. T. Tagawa, M. Albanese, et. al. [Epstein-Barr viral miRNAs inhibit antiviral CD4+ T cell responses targeting IL-12 and peptide processing](#). *J Exp Med.* **213**(10), 2065–2080 (2016) [human B cells, DG-75]
72. D.L. Hertz, K. Owzar, et. al. [Pharmacogenetic Discovery in CALGB \(Alliance\) 90401 and Mechanistic Validation of a VAC14 Polymorphism that Increases Risk of Docetaxel-Induced Neuropathy](#). *Clin Cancer Res.* **22**, 4890–4900 (2016). [stem cell-derived peripheral neuronal cells]
73. S. Das, M. Miller, et. al. [GSDMB induces an asthma phenotype characterized by increased airway responsiveness and remodeling without lung inflammation](#). *PNAS.* **113** (46) 13132–13137 (2016) [human lung epithelial cells]
74. I.B. Sharum, S. Granados-Aparici, et. al. [Serine threonine kinase receptor associated protein regulates early follicle development in the mouse ovary](#). *Reproduction*, Dec 2016; **153**: 221–231. [neonatal mouse ovary]
75. T.D. Burt, J.M. McCune et. al. [Lin28b Regulates Fetal Regulatory T Cell Differentiation through Modulation of TGF-β Signaling](#). *J. Immunol.*, Dec 2016; **197**: 4344–4350. [naïve fetal T cells]
76. K. Morita, T. Okamura, et. al. [Egr2 and Egr3 in regulatory T cells cooperatively control systemic autoimmunity through Ltpb3-mediated TGF-β3 production](#). *PNAS*, Nov 2016; **113**(10): 1611286114. [regulatory T cells]
77. N.K. Banda, S. Acharya, et. al. [Mannan-Binding Lectin–Associated Serine Protease 1/3 Cleavage of Pro–Factor D into Factor D In vivo and Attenuation of Collagen Antibody-Induced Arthritis through Their Targeted Inhibition by RNA Interference–Mediated Gene Silencing](#). *J. Immunol.*, Nov 2016; **197**: 3680–3694. [*In vivo* mouse injection; liver and adipose tissue knockdown]
78. Y. Deng, J. Zhao, et. al. [Decreased SMG7 expression associates with lupus-risk variants and elevated antinuclear antibody production](#). *Ann Rheum Dis*, Nov 2016; **75**: 2007–2013. [PBMCs]
79. K.A. Mitchnick, S. Creighton, et. al. [Differential contributions of de novo and maintenance DNA methyltransferases to object memory processing in the rat hippocampus and perirhinal cortex—a double dissociation](#). *European Journal of Neuroscience*, 1–14 (2014) [infusion via intracranial cannula (rat)]
80. L. Huang et al. [Phosphoinositide 3-Kinase Gamma Contributes to Neuroinflammation in a Rat model of Surgical Brain Injury](#). *The Journal of Neuroscience*, **35**(29): 10390–10401 (2015) [rat brain, intracerebroventricular administration]
81. H. Nakajima et al. [Nuclear-translocated Glyceraldehyde-3-phosphate Dehydrogenase Promotes Poly\(ADP-ribose\) Polymerase-1 Activation during Oxidative/Nitrosative Stress in Stroke](#). *Journal of Biological Chemistry*, Vol. 290 No. 23 (14493-14503)(2015) [rat brain-intracerebroventricular injection]
82. S. L. Jaslow et al. [Salmonella Activation of STAT3 Signaling by SarA Effector Promotes Intracellular Replication and Production of IL-10](#). *Cell Reports* **19**;23(12):3525-3536 (2018) [LCLs, human lymphoblastoid cell lines]
83. M. Hemshekhar et al. [Host Defense Peptide LL-37-Mediated Chemoattractant Properties, but Not Anti-Inflammatory Cytokine IL-1RA Production, Is Selectively Controlled by Cdc42 Rho GTPase via G Protein-Coupled Receptors and JNK Mitogen-Activated Protein Kinase](#). *Frontiers in Immunology* **9**:1871 (2018). [Human monocytic THP-1]
84. X. Zhou et al. [Modulation of Mutant KrasG12D-Driven Lung Tumorigenesis In Vivo by Gain or Loss of PCDH7 Function](#). *Molecular Cancer Research* **17**(2):594-603 (2019). [HBECs, human bronchial epithelial cells]
85. M. B. Maina et al. [The involvement of tau in nucleolar transcription and the stress response](#). *Acta Neuropathologica Communications* **6**, Article number: 70. [SHSY5Y, human neuroblastoma cells]
86. G. H. Wabnitz et al. [Protein Phosphatase 1 and Cofilin Regulate Nuclear Translocation of NF-κB and Promote Expression of the Anti-Inflammatory Cytokine Interleukin-10 by T Cells](#). *Molecular and Cellular Biology* **38**(22) (2018). [Primary human T-cells]
87. J.C. Gerlach et al. [Effects of Delta-Like Noncanonical Notch Ligand 1 Expression of Human Fetal Liver Hepatoblasts on Hematopoietic Progenitors](#). *Stem Cells International* **7916275** (2019). [Human fetal liver cells]
88. J. Schulthess et al. [The Short Chain Fatty Acid Butyrate Imprints an Antimicrobial Program in Macrophages](#). *Immunity* **50**(2) (2019). [Mouse, macrophages]
89. T.J. Jarome et al. [EZH2 Methyltransferase Activity Controls Pten Expression and mTOR Signaling during Fear Memory Reconsolidation](#). *The Journal of Neuroscience* **38**(35):7635-7648 (2018). [*In vivo* rat brain]
90. S. Schwörer et al. [Epigenetic stress responses induce muscle stem-cell ageing by Hoxa9 developmental signals](#). *Nature* **540**(7633):428-432 (2016). [*In vivo* mouse tibialis anterior (TA) muscle]
91. S. Sin Min Dong et al. [Transfection of hard-to-transfect primary human macrophages with Bax siRNA to reverse Resveratrol-induced apoptosis](#). *RNA Biology* **17**(6):755-764 (2020). [Primary human macrophages]
92. T. Hazzan et al. [An efficient method for gene knock-down by RNA interference in human skin mast cells](#). *Experimental Dermatology* **26**(11):1136-1139 (2017). [Human skin mast cells]

For more information

To find the contact information in your country for your technology of interest, please visit us at [horizondiscovery.com/contact-us](#)

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

©2021 The Horizon logo and other trademarks are the property of Horizon Discovery Limited, unless otherwise stated. DHARMACON and EDIT-R are trademarks of Dharmacon Inc.

horizon