

<Plasmid DNA 導入文献>

導入細胞名	論文名
HEK293 (ヒト胎児腎臓)	M. Ohno, <i>et al.</i> , "Interactions of Thyroid Hormone Receptor with Ku Proteins and Interleukin Enhancer Binding Factor 3 Modulate the Promoter Activity of Thyroid-Stimulating Hormone Alpha", <i>Biol. Pharm. Bull.</i> , 2012 , 35(3), 380.
	S. Hino, <i>et al.</i> , "Discharge of solubilized and Dectin-1-reactive β -glucan from macrophage cells phagocytizing insoluble β -glucan particles: Involvement of reactive oxygen species (ROS)-driven degradation", <i>Biochem. Biophys. Res. Commun.</i> , 2012 , 421(2), 329.
	G. -A. Song, <i>et al.</i> , "Molecular Consequences of the ACVR1R206H Mutation of Fibrodysplasia Ossificans Progressiva", <i>J. Biol. Chem.</i> , 2010 , 285(29), 22542.
	S. Yamamoto, <i>et al.</i> , "Identification and Functional Characterization of the First Nucleobase Transporter in Mammals", <i>J. Biol. Chem.</i> , 2010 , 285(9), 6522.
	T. Shimizu, <i>et al.</i> , "Zinc finger genes Fezf1 and Fezf2 control neuronal differentiation by repressing Hes5 expression in the forebrain", <i>Development</i> , 2010 , 137(11), 1875.
	H. Nojima, <i>et al.</i> , "Syntabulin, a motor protein linker, controls dorsal determination.", <i>Development</i> , 2010 , 137(6), 923.
	S. Yamamoto, <i>et al.</i> , "Identification and functional characterization of rat riboflavin transporter 2", <i>J. Biol. Chem.</i> , 2009 , 145(4), 437.
HEK293T (ヒト胎児腎臓)	K. E. Lee, <i>et al.</i> , "HIV-1 Capsid-Targeting Domain of Cleavage and Polyadenylation Specificity Factor 6", <i>J. Virol.</i> , 2012 , 86(7), 3851.
	S. Ohata, <i>et al.</i> , "Dual roles of Notch in regulation of apically restricted mitosis and apicobasal polarity of neuroepithelial cells.", <i>Neuron</i> , 2011 , 69(2), 215.
	A. Shimahara, <i>et al.</i> , "Acetylation of Lysine 564 Adjacent to the C-terminal Binding Protein-binding Motif in EVI1 Is Crucial for Transcriptional Activation of GATA2", <i>J. Biol. Chem.</i> , 2010 , 285(22), 16967.
	V. Futak, <i>et al.</i> , "Perturbation of the P-body component Mov10 inhibits HIV-1 infectivity.", <i>PLoS One</i> , 2010 , 5(2), e9081.
	M. Ohno, <i>et al.</i> , "Repression of the Promoter Activity Mediated by Liver Receptor Homolog-1 through Interaction with Ku Proteins", <i>Biol. Pharm. Bull.</i> , 2010 , 33(5), 784.
	O.-J. Park, <i>et al.</i> , "FGF2-activated ERK Mitogen-activated Protein Kinase Enhances Runx2 Acetylation and Stabilization", <i>J. Biol. Chem.</i> , 2010 , 285(6), 3568.
	K. Koiwai, <i>et al.</i> , "BPOZ-2 directly binds to eEF1A1 to promote eEF1A1 ubiquitylation and degradation and prevent translation", <i>Genes to Cells</i> , 2008 , 13(6), 593.
T. Kohno, <i>et al.</i> , "Protein-tyrosine kinase CAKb/PYK2 is activated by binding Ca ²⁺ /calmodulin to FERM F2 α 2 helix and thus forming its dimer", <i>Biochem. J.</i> , 2008 , 410, 513.	
HeLa (ヒト子宮頸癌)	H. Kagawa, <i>et al.</i> , "A novel signaling pathway mediated by the nuclear targeting of C-terminal fragments of mammalian Patched 1.", <i>PLoS One</i> , 2011 , 6(4), e18638.
	K. Shiga, <i>et al.</i> , "Development of an intracellularly acting inhibitory peptide selective for PKN", <i>Biochem. J.</i> , 2010 , 425, 445.
	Y. Enoki, <i>et al.</i> , "DNA-binding and transcriptional activities of human HSF4 containing mutations that associate with congenital and age-related cataracts", <i>Biochim. Biophys. Acta</i> , 2010 , 1802(9), 749.
	T. Murakami, <i>et al.</i> , "Ephedrae Herba, a Component of Japanese Herbal Medicine Mao-to, Efficiently Activates the Replication of Latent Human Immunodeficiency Virus Type 1 (HIV-1) in a Monocytic Cell Line", <i>Biol. Pharm. Bull.</i> , 2008 , 31(12), 2334.
	M. Soejima, <i>et al.</i> , "Distinct single nucleotide polymorphism pattern at the FUT2 promoter among human populations", <i>Ann. Hematol.</i> , 2008 , 87(1), 19.
	A. Ryo, <i>et al.</i> , "A Suppressive Role of the Prolyl Isomerase Pin1 in Cellular Apoptosis Mediated by the Death-associated Protein Daxx", <i>J. Biol. Chem.</i> , 2007 , 282(50), 36671.
MCF-7 (ヒト哺乳器)	Z. Zhang, <i>et al.</i> , "Fucoïdan extract induces apoptosis in MCF-7 cells via a mechanism involving the ROS-dependent JNK activation and mitochondria-mediated pathways", <i>PLoS One</i> , 2011 , 6(11), e27441.
	W. Keun, <i>et al.</i> , "Amurensin G, a Potent Natural SIRT1 Inhibitor, Rescues Doxorubicin Responsiveness via Down-Regulation of Multidrug Resistance 1", <i>Mol. Pharmacol.</i> , 2010 , 78(5), 855.
	M. R. Kim, <i>et al.</i> , "Involvement of Pin1 induction in epithelial-mesenchymal transition of tamoxifen-resistant breast cancer cells", <i>Cancer Science</i> , 2009 , 100(10), 1834.
	C. -Y. Han, <i>et al.</i> , "Role of FoxO1 activation in MDR1 expression in adriamycin-resistant breast cancer cells", <i>Carcinogenesis</i> , 2008 , 29(9), 1837.
A549 (ヒト肺上皮癌)	Y. Nishimoto, <i>et al.</i> , "Glycyrrhizin Attenuates Mucus Production by Inhibition of MUC5AC mRNA Expression In Vivo and In Vitro", <i>J. Pharmacol. Sci.</i> , 2010 , 113(1), 76.
	T. Kawai, <i>et al.</i> , "Associations between decay-accelerating factor polymorphisms and allergic respiratory diseases", <i>Clin. exp. Allergy</i> , 2009 , 39(10), 1508.
	H. Takei, <i>et al.</i> , "Glycyrrhizin Inhibits Interleukin-8 Production and Nuclear Factor- κ B Activity in Lung Epithelial Cells, but Not Through Glucocorticoid Receptors", <i>J. Pharmacol. Sci.</i> , 2008 , 106(3), 460.
HepG2 (ヒト肝癌)	M. Taura, <i>et al.</i> , "MEF/ELF4 transactivation by E2F1 is inhibited by p53", <i>Nucl. Acid Res.</i> , 2011 , 39(1), 76.
	Y. Kuwahara, <i>et al.</i> , "Enhancement of autophagy is a potential modality for tumors refractory to radiotherapy", <i>Cell Death Dis.</i> , 2011 , 2, e177.
	S. Morino, <i>et al.</i> , "Mild Electrical Stimulation with Heat Shock Ameliorates Insulin Resistance via Enhanced Insulin Signaling", <i>PLoS One</i> , 2008 , 3(12), e4068
BAEC (ウシ大動脈内皮細胞)	川崎 那緒人, <i>et al.</i> , "光転換型蛍光タンパク質を用いた血管内皮細胞の微小管構造の可視化", <i>生体医工学</i> , 2009 , 47(6), 595.
	A. Endo, <i>et al.</i> , "1,2-Naphthoquinone disrupts the function of cAMP response element-binding protein through covalent modification", <i>Biochem. Biophys. Res. Commun.</i> , 2007 , 361(1), 243.
C2C12 (マウス筋芽細胞)	S. Ohata, <i>et al.</i> , "Dual roles of Notch in regulation of apically restricted mitosis and apicobasal polarity of neuroepithelial cells.", <i>Neuron</i> , 2011 , 69(2), 215.
	M. Ogawa, <i>et al.</i> , "17 β -Estradiol Represses Myogenic Differentiation by Increasing Ubiquitin-specific Peptidase 19 through Estrogen Receptor α ", <i>J. Biol. Chem.</i> , 2011 , 286(48), 41455.

HaCaT (ヒト皮膚角化細胞)	K. Park, <i>et al.</i> , "Regulation of cathelicidin antimicrobial peptide expression by an endoplasmic reticulum (ER) stress signaling, vitamin D receptor-independent pathway", <i>J. Biol. Chem.</i> , 2011 , 286(39), 34121. K. Park, <i>et al.</i> , "Regulation of cathelicidin expression by an ER stress-signaling, vitamin D receptor-independent pathway", <i>J. Biol. Chem.</i> , 2011 , 286(39), 34121.
NIH3T3 (マウス胎仔線維芽細胞)	A. Takahashi, <i>et al.</i> , "A novel potent tumour promoter aberrantly overexpressed in most human cancers", <i>Sci. Rep.</i> , 2011 , 1, 15. R. Doi, <i>et al.</i> , "CLOCK Regulates Circadian Rhythms of Hepatic Glycogen Synthesis through Transcriptional Activation of Gys2", <i>J. Biol. Chem.</i> , 2010 , 285(29), 22114.
HCT116 p53 (ヒト結腸癌)	M. Taura, <i>et al.</i> , "MEF/ELF4 transactivation by E2F1 is inhibited by p53", <i>Nucl. Acid. Res.</i> , 2011 , 39(1), 76. M. Taura, <i>et al.</i> , "p53 Regulates Toll-Like Receptor 3 Expression and Function in Human Epithelial Cell Lines", <i>Mol. Cell. Biol.</i> , 2008 , 28(21), 6557.
adrenal chromaffin cells (ウシ副腎髄質細胞)	T. Memoto, <i>et al.</i> , "Homologous posttranscriptional regulation of insulin-like growth factor-I receptor level via glycogen synthase kinase-3 β and mammalian target of rapamycin in adrenal chromaffin cells: Effect on tau phosphorylation", <i>Neuropharmacology</i> , 2010 , 58(7), 1097.
B16 (マウスメラノーマ)	I. Shirasugi, <i>et al.</i> , "Novel Screening Method for Potential Skin-Whitening Compounds by a Luciferase Reporter Assay", <i>Biosci. Biotechnol. Biochem.</i> , 2010 , 74(11), 2253.
Caco2 (ヒト腸管上皮)	M. Saeki, <i>et al.</i> , "Functional analysis of genetic variations in the 5'-flanking region of the human MDR1 gene", <i>Mol. Genet. Metabol.</i> , 2011 , 102(1), 91.
CHO (チャイニーズハムスター卵巣)	I. Tanida, <i>et al.</i> , "Consideration about negative controls for LC3 and expression vectors for four colored fluorescent protein-LC3 negative controls", <i>Autophagy.</i> , 2008 , 4(1), 131.
COS7 (アフリカミドリサル腎臓)	A. Shimahara, <i>et al.</i> , "Acetylation of Lysine 564 Adjacent to the C-terminal Binding Protein-binding Motif in EVI1 Is Crucial for Transcriptional Activation of GATA2", <i>J. Biol. Chem.</i> , 2010 , 285(22), 16967.
DLD-1 (ヒト結腸腺癌)	M. Koyama, <i>et al.</i> , "Histone Deacetylase Inhibitors and 15-Deoxy- Δ 12,14-Prostaglandin J2 Synergistically Induce Apoptosis", <i>Clin. Cancer. Res.</i> , 2010 , 16(8), 2320.
GH3 (ラット下垂体癌)	S. Hasimoto, <i>et al.</i> , "Protein disulfide isomerase regulates the thyroid hormone receptor-mediated gene expression via redox factor-1 through thiol reduction-oxidation", <i>J. Biol. Chem.</i> , 2012 , 288, 1706.
H4IIE (ラット肝癌)	H. Hashizume, <i>et al.</i> , "Histone Methyltransferase PR-Set7 and Histone Variant H2A.Z, Induced during Hepatocarcinogenesis, Repress the Promoter Activity of the Tumor Marker Gene and the Ras-Induced Colony Formation Activity", <i>J. Health Sci.</i> , 2011 , 57(3), 264.
HT29 (ヒト結腸癌)	M. Soejima, <i>et al.</i> , "Distinct single nucleotide polymorphism pattern at the FUT2 promoter among human populations", <i>Ann., Hematol.</i> , 2008 , 87(1), 19.
HUVEC (ヒト臍帯静脈内皮細胞)	X. Zhu, <i>et al.</i> , "A high-throughput fluorescence resonance energy transfer (FRET)-based endothelial cell apoptosis assay and its application for screening vascular disrupting agents", <i>Biochem. Biophys. Res. Commun.</i> , 2012 , 418(4), 641.
MCAS (ヒト卵巣癌)	M. Soejima, <i>et al.</i> , "Distinct single nucleotide polymorphism pattern at the FUT2 promoter among human populations", <i>Ann. Hematol.</i> , 2008 , 87(1), 19.
REF (ラット初代線維芽細胞)	T. Kuroshima, <i>et al.</i> , "Viral-mediated stabilization of AU-rich element containing mRNA contributes to cell transformation", <i>Oncogene</i> , 2011 , 30(26), 2912.
ROS 17/2.8 (ラット骨芽細胞様細胞)	Y. D. Cho, <i>et al.</i> , "Molecular Regulation of Matrix Extracellular Phosphoglycoprotein Expression by Bone Morphogenetic Protein-2", <i>J. Biol. Chem.</i> , 2009 , 284(37), 25230.
Tn5 (昆虫細胞)	A. I. Guce, <i>et al.</i> , "Catalytic Mechanism of Human α -Galactosidase", <i>J. Biol. Chem.</i> , 2010 , 285(6), 3625.

<siRNA 導入文献>

導入細胞名	論文名
bovine fibroblast cell (ウシ線維芽)	K. Yamanaka, <i>et al.</i> , "Gene Silencing of DNA Methyltransferases by RNA Interference in Bovine Fibroblast Cells", <i>J. Reprod. Dev.</i> , 2010 , 56(1), 60.
EF/CT-1	T. Sakurai, <i>et al.</i> , "Function of a Transcription Factor CDX2 Beyond Its Trophoblast Lineage Specification", <i>Endocrinology</i> , 2010 , 151(12), 5873.
HaCaT (ヒト皮膚角化細胞)	H. Takeuchi, <i>et al.</i> , "Characterization of PDK as a Protein Involved in Epidermal Growth Factor Receptor Trafficking", <i>Mol. Cell. Biol.</i> , 2010 , 30(7), 1689.
HEK293T (ヒト胎児腎臓)	K. Koiwai, <i>et al.</i> , "TdfI2 is a nucleolar protein that promotes rRNA gene promoter activity", <i>Genes. Cells</i> , 2011 , 16(7), 748.
HeLa (ヒト子宮頸癌)	Y. Kim, <i>et al.</i> , "Redox regulation of the tumor suppressor PTEN by glutathione", <i>FEBS Letters</i> , 2010 , 584(16), 3550.
HSG (ヒト顎下腺細胞)	Y. Onishi, "HSG cells, a model in the submandibular clock", <i>Biosci. Rep.</i> , 2011 , 31(1), 57.
NHLF (ヒト正常肺線維芽)	S. Y. Kim, <i>et al.</i> , "Cigarette smoke induces Akt protein degradation by the ubiquitin-proteasome system", <i>J. Biol. Chem.</i> , 2011 , 286(37), 31932.
NIH3T3 (マウス胎仔線維芽細胞)	Y. Onishi, <i>et al.</i> , "Rhythmic binding of Topoisomerase I impacts on the transcription of Bmal1 and circadian period", <i>Nucl. Acids Res.</i> , 2012 , 40(19), 9482.
THP-1 (ヒト単球性白血病)	S. -H. Lee, <i>et al.</i> , "The major outer membrane protein of a periodontopathogen induces IFN- β and IFN-stimulated genes in monocytes via lipid raft and TANK-binding kinase 1/IFN regulatory factor-3.", <i>J. Immunol.</i> , 2009 , 182(9), 5823.

○ お問い合わせ先 ○
株式会社同仁化学研究所
カスタマーリレーション部
E-mail: info@dojindo.co.jp Tel: 0120-489-548