

Active Motif Epigenetic Services Publications

The papers below contain data generated by Active Motif's Epigenetic Services team. To learn more about our services, please give us a call or visit us at www.activemotif.com/services.

Technique	Target	Journal	Year	Reference
ChIP-Seq and ChIP-qPCR	CBP, p300, H3K27Ac	Cell Reports	2019	Denise de Almeida Nagata <i>et al.</i> Regulation of Tumor-Associated Myeloid Cell Activity by CBP/EP300 Bromodomain Modulation of H3K27 Acetylation. <i>Cell Reports</i> . 27:269
ChIP-qPCR	ChREBP	Nature Metabolism	2019	Pauline Morigny <i>et al.</i> Interaction between hormone-sensitive lipase and ChREBP in fat cells controls insulin sensitivity. <i>Nature Metabolism</i> . 1: 133
ChIP-Seq	H3K4me1, H3K4me3, H3K27Ac	Cell Metabolism	2019	Liming Du <i>et al.</i> IGF-2 Preprograms Maturing Macrophages to Acquire Oxidative Phosphorylation-Dependent Antiinflammatory Properties. <i>Cell Metabolism</i> . 29: 1
ChIP-qPCR	SRF and CEBP	Nature Communications	2019	Shilpita Sarcar <i>et al.</i> Next-generation muscle-directed gene therapy by in silico vector design. <i>Nat Commun</i> . https://doi.org/10.1038/s41467-018-08283-7
ChIP-Seq	H3K27ac, H3K4me3, CTCF, and NR4A3	Nature Communications	2019	Florian Haller <i>et al.</i> Enhancer hijacking activates oncogenic transcription factor NR4A3 in acinic cell carcinomas of the salivary glands. <i>Nat Commun</i> . 10: 336
ChIP-Seq	Gfi1, Lsd1	Nature Communications	2019	Catherine Lee <i>et al.</i> Lsd1 as a therapeutic target in Gfi1-activated medulloblastoma. <i>Nat Commun</i> . 10: 332
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K9me3, H3K36me3, EZH2, SMARCB1, SUZ12, SMARCA4, REST	Cancer Cell	2019	Serap Erkek <i>et al.</i> Comprehensive Analysis of Chromatin States in Atypical Teratoid/Rhabdoid Tumors Identified Diverging Roles for SWI/SNF and Polycomb in Gene Regulation. <i>Cancer Cell</i> . 35: 95
ChIP-Seq	BRG1	The Journal of Clinical Investigation	2019	Yufeng Ding <i>et al.</i> Chromatin remodeling ATPase BRG1 and PTEN are synthetic lethal in prostate cancer. <i>J Clin Invest</i> . https://doi.org/10.1172/JCI123557
ChIP-Seq, RNA-Seq	GR	Journal of the American Heart Association	2019	Elena Severinova <i>et al.</i> Glucocorticoid Receptor Binding and Transcriptome Signature in Cardiomyocytes. <i>J Am Heart Assoc</i> . 8:e011484. DOI: 10.1161/JAHA.118.011484.
ChIP-Seq	H4K20me1	Journal of Autoimmunity	2019	Keqi Fan <i>et al.</i> CRL4DCAF2 is required for mature T-cell expansion via Aurora B-regulated proteasome activity. <i>Journal of Autoimmunity</i> . 96:74
MeDIP-Seq	—	Epigenomes	2019	Liliana Ferreira <i>et al.</i> Uncovering Differentially Methylated Regions (DMRs) in a Salt-Tolerant Rice Variety under Stress: One Step towards New Regulatory Regions for Enhanced Salt Tolerance. <i>Epigenomes</i> . 3:4
ChIP-Seq	RNA Pol II, H3K27Ac	Journal of Neuro-Oncology	2019	Isabel Tegeder <i>et al.</i> Functional relevance of genes predicted to be affected by epigenetic alterations in atypical teratoid/rhabdoid tumors. <i>J. Neurooncol</i> . https://doi.org/10.1007/s11060-018-03018-6
ChIP-Seq	EGRI, H3K27Ac	Journal of Steroid Biochemistry & Molecular Biology	2019	Maria Szwarc <i>et al.</i> Early growth response 1 transcriptionally primes the human endometrial stromal cell for decidualization. <i>J. Steroid Biochem. Mol. Biol</i> . https://doi.org/10.1016/j.jsbmb.2019.01.021

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3, H3K4me3	Acta Neuropathologica	2019	Andre Silveira <i>et al.</i> H3.3 K27M depletion increases differentiation and extends latency of diffuse intrinsic pontine glioma growth in vivo. <i>Acta Neuropathol.</i> https://doi.org/10.1007/s00401-019-01975-4
ChIP-Seq	RNA Pol II	Agronomy	2019	Sonja Klemme <i>et al.</i> Selection of Salicylic Acid Tolerant Epilines in <i>Brassica napus</i> . <i>Agronomy</i> . 9: 92
ChIP-Seq	LSD1, H3K4me2, H3K27Ac	Science Signaling	2019	Arnaud Augert <i>et al.</i> Targeting NOTCH activation in small cell lung cancer through LSD1 inhibition. <i>Sci.Signal.</i> 12, eaau2922
ChIP-qPCR	PPAR gamma	International Journal of Obesity	2019	Laura Butruille <i>et al.</i> Maternal high-fat diet during suckling programs visceral adiposity and epigenetic regulation of adipose tissue stearoyl-CoA desaturase-1 in offspring. <i>Int. J. Obes.</i> https://doi.org/10.1038/s41366-018-0310-z
ChIP-Seq	HA-tagged JP2NT, TBP, MEF2C	Science	2018	Ang Guo <i>et al.</i> E-C coupling structural protein junctophilin-2 encodes a stress-adaptive transcription regulator. <i>Science</i> . 362: 6421
ChIP-qPCR	p65/RelA	Human Molecular Genetics	2018	Jeffrey R. Gehlhausen <i>et al.</i> A proteasome-resistant fragment of NIK mediates oncogenic NF- κ B signaling in schwannomas. <i>Clin Cancer Res.</i> 25:1601. doi: 10.1093/hmg/ddy361
ChIP-Seq	Nrf2	Journal of Biological Chemistry	2018	Junsheng Fu <i>et al.</i> Hyperactivity of the transcription factor Nrf2 causes metabolic reprogramming in mouse esophagus. <i>J. Biol. Chem.</i> 294: 327
ChIP-Seq	H3K36me2	Cell reports	2018	Jingjing Chen <i>et al.</i> Methyltransferase Nsd2 Ensures Germinal Center Selection by Promoting Adhesive Interactions between B Cells and Follicular Dendritic Cells. <i>Cell Reports</i> . 25: 3393
ChIP-Seq	H3K27Ac	Clinical Cancer Research	2018	Shori Saito <i>et al.</i> Eradication of central nervous system leukemia of T-cell origin with a brain-permeable LSD1 inhibitor. <i>Clin Cancer Res.</i> 25:1601
ChIP-Seq	RNA Pol II, Brd4, H3K9me3	Cell Reports	2018	Lu Want <i>et al.</i> Retinal Cell Type DNA Methylation and Histone Modifications Predict Reprogramming Efficiency and Retinogenesis in 3D Organoid Cultures. <i>Cell Reports</i> . 22: 2601
ATAC-Seq	NA	Cell Reports	2018	Xiaolong Zhang <i>et al.</i> OX40 Costimulation inhibits Foxp3 expression and Treg induction via BATF3-dependent and independent mechanisms. <i>Cell Reports</i> . 24: 607.
Mod Spec	NA	Acta Neuropathologica	2018	Leah Katz <i>et al.</i> Loss of histone H3K27me3 identifies a subset of meningiomas with increased risk of recurrence. <i>Acta Neuropathol.</i> 135: 6.
RIME	TRPS1	Cell Reports	2018	Robert M. Witwicki <i>et al.</i> TRPS1 is a lineage-specific transcriptional dependency in breast cancer. <i>Cell</i> . 175: 1.
ChIP-Seq	H3K27Ac	Cell	2018	Xun Huang <i>et al.</i> Targeting epigenetic crosstalk as a therapeutic strategy for EZH2-aberrant solid tumors. <i>Cell</i> . 175: 1.
ChIP-Seq	BRD4, RNA Pol II	Cancer Cell	2018	Elizabeth Stewart <i>et al.</i> Identification of therapeutic targets in rhabdomyosarcoma through integrated genomic, epigenomic, and proteomic analyses. <i>Cancer Cell</i> . 34: 1.
ChIP-Seq	Sox17	Nature Communications	2018	Xiaoqi Wang <i>et al.</i> SOX17 regulates uterine epithelial-stromal cross-talk acting via a distal enhancer upstream of <i>Ihh</i> . <i>Nat Commun.</i> 9: 4421
ChIP-Seq, RNA-Seq	NRF2, RNA Pol II	Nature Communications	2018	David Olganier <i>et al.</i> Nrf2 negatively regulates STING indicating a link between antiviral sensing and metabolic reprogramming. <i>Nat Commun.</i> 9:3506

Technique	Target	Journal	Year	Reference
ChIP-Seq, RIME	p300, CBP, CDK9, BRD4, H3K27Ac, H3K4me1, H3K4me3, H3K18Ac	Cell Reports	2018	Ryan Raisner <i>et al.</i> Enhancer activity requires CBP/P300 bromodomain-dependent histone H3K27 acetylation. <i>Cell Reports</i> . 24: 1722.
ChIP-Seq, RNA-Seq	H3K27Ac, Med1, RNA Pol II	The Journal of Clinical Investigation	2018	Maria Florencia Martinez <i>et al.</i> Super-Enhancers maintain renin-expressing cell identity and memory to preserve multi-system homeostasis. <i>J. Clin Invest.</i> Advance Online Publication doi: 10.1172/JCI121361.
ChIP-Seq,	BRD4, H3K27Ac	JCI Insight	2018	Anisley Valenciana <i>et al.</i> Transcriptional targeting of oncogene addiction in medullary thyroid cancer. <i>JCI Insight</i> . 3: e122225.
RNA-Seq	NA	PAIN	2018	Pradipta Ray <i>et al.</i> Comparative transcriptome profiling of the human and mouse dorsal root ganglia: an RNA-Seq based resource for pain and sensory neuroscience research. <i>Pain</i> . 159: 7.
ChIP-Seq	ELL2	The Journal Of Immunology	2018	Ashley M. Nelson <i>et al.</i> RNA splicing in the transition from B cells to antibody-secreting cells: The influences of ELL2, small nuclear RNA, and endoplasmic reticulum stress. <i>J Immunol.</i> Advance online publication doi: 10.4049/jimmunol
ChIP-Seq	MTA1	Cancer Medicine	2018	Nasir A. Butt <i>et al.</i> Targeting MTA1/HIF-1 α signaling by pterostilbene in combination with histone deacetylase inhibitor attenuates prostate cancer progression. <i>Cancer Med</i> . 6: 2673
ChIP-Seq, Next-Gen Bisulfite-Seq	H3K4me3, RNA Pol II	Agronomy	2018	Martin Schmidt <i>et al.</i> Methylome and epialleles in rice epilines selected for energy use efficiency. <i>Agronomy</i> . 8: 163
ChIP-Seq	GATA2	The American Journal of Human Genetics	2018	Katelyn M. Mika <i>et al.</i> An ancient fecundability-associated polymorphism creates a GATA2 binding site in a distal enhancer of HLA-F. <i>Am J Hum Genet</i> . 103: 509
ChIP-Seq	FOXL2	Human Molecular Genetics	2018	Barbara Nicol <i>et al.</i> Genome-wide identification of FOXL2 binding and characterization of FOXL2 feminizing action in the fetal gonads. <i>Hum Mol Genet.</i> Advance online publication doi: 10.1093/hmg/ddy312
MeDIP-Seq	5-Methylcytosine	Scientific Reports	2018	Jose P. Silva <i>et al.</i> Analysis of diet-induced differential methylation, expression, and interactions of lncRNA and protein-coding genes in mouse liver. <i>Sci Rep</i> . 8: 11537
ChIP-Seq	ZFP24	Cell Reports	2018	Benayahu Elbaz <i>et al.</i> Phosphorylation state of ZFP24 controls oligodendrocyte differentiation. <i>Cell Reports</i> . 23: 2254
ChIP-Seq, Histone PTM Quantitation	H3K9me3, H3K27me3	Cell Reports	2018	Jessica Camacho <i>et al.</i> The memory of environmental chemical exposure in <i>C. elegans</i> is dependent on the jumonji demethylases <i>jmjd-2</i> and <i>jmjd-3/utx-1</i> . <i>Cell Reports</i> . 23: 2392
ChIP-Seq	TET1, DNMT1	Nature Genetics	2018	Nipun Verma <i>et al.</i> TET proteins safeguard bivalent promoters from de novo methylation in human embryonic stem cells. <i>Nat Genet</i> . 50: 83
ChIP-Seq	EZH2	Journal of Experimental Medicine	2018	Xingli Zhang <i>et al.</i> Macrophage/microglial Ezh2 facilitates autoimmune inflammation through inhibition of <i>Socs3</i> . <i>J Exp Med</i> . 215: 1365
ChIP-Seq	STAT3, STAT5, ROCK2	Scientific Reports	2018	Wei Chen <i>et al.</i> ROCK2, but not ROCK1 interacts with phosphorylated STAT3 and co-occupies TH17/TFH gene promoters in TH17-activated human T cells. <i>Sci Rep</i> . 8: 16636
ChIP-Seq	Bhlhe40	PLoS One	2018	Kelly A. Hamilton <i>et al.</i> Mice lacking the transcriptional regulator Bhlhe40 have enhanced neuronal excitability and impaired synaptic plasticity in the hippocampus. <i>PLoS One</i> . 13: e0196223

Technique	Target	Journal	Year	Reference
ChIP-Seq	RNA Pol II	PLoS One	2018	Liana Basova <i>et al.</i> Dopamine and its receptors play a role in the modulation of CCR5 expression in innate immune cells following exposure to Methamphetamine: Implications to HIV infection. <i>PLoS One</i> . 13: e0199861
ChIP-qPCR	Glucocorticoid Receptor, TFI	Arthritis & Rheumatology	2018	Yanhua Hu <i>et al.</i> Development of a molecular signature to monitor pharmacodynamic response mediated by in vivo administration of glucocorticoids. <i>Arthritis Rheumatol</i> . 70: 1331
ChIP-Seq	H3K27me3	Acta Neuropathologica Communications	2018	David Castel <i>et al.</i> Transcriptomic and epigenetic profiling of 'diffuse midline gliomas, H3 K27M-mutant' discriminate two subgroups based on the type of histone H3 mutated and not supratentorial or infratentorial location. <i>Acta Neuropathol Commun</i> . 6: 17
ChIP-Seq	NRF2	Molecular Pharmacology	2018	Rance Nault <i>et al.</i> Comparison of hepatic NRF2 and AHR binding in 2,3,7,8- tetrachlorodibenzo-p-dioxin (TCDD) treated mice demonstrates NRF2- independent PKM2 induction. <i>Mol Pharmacol</i> . 94: 876
ChIP-Seq	HDAC2, H3K27Ac, H3K27me3	Cancer Immunology, Immunotherapy	2018	David Briere <i>et al.</i> The class I/IV HDAC inhibitor mocetinostat increases tumor antigen presentation, decreases immune suppressive cell types and augments checkpoint inhibitor therapy. <i>Cancer Immunol Immunother</i> . 67: 381
ChIP-Seq, ChIP-qPCR	TRRAP	Journal of Cell Biology	2018	Zhao Wang <i>et al.</i> TRRAP is a central regulator of human multiciliated cell formation. <i>J Cell Biol</i> . 217: 1941
ChIP-Seq	BRD4	Blood	2018	Rebecca Kohnken <i>et al.</i> Diminished microRNA-29b level is associated with BRD4 mediated activation of oncogenes in cutaneous T-cell lymphoma. <i>Blood</i> . 131: 771
RNA-Seq	NA	Cell Communication and Signaling	2018	Keri Callegari <i>et al.</i> Pharmacological inhibition of LSD1 activity blocks REST-dependent medulloblastoma cell migration. <i>Cell Commun Signal</i> . 16: 60
ChIP-Seq	MTA1	Molecular Oncology	2018	Avinash Kumar <i>et al.</i> MTA drives malignant progression and bone metastasis in prostate cancer
ChIP-Seq	RNA Pol II, H2A.Z, CDK9, ANP32e	Biochimica et Biophysica Acta	2018	Hyewon Shin <i>et al.</i> Transcriptional regulation mediated by H2A.Z via ANP32e-dependent inhibition of protein phosphatase 2A. <i>Biochim Biophys Acta</i> . 1861: 481
ChIP-Seq	H3K4me3, H3K9me3, H3K27me3	Proceedings of the Royal Society B	2018	Theresa K. Kelly <i>et al.</i> Epigenetic regulation of transcriptional plasticity associated with developmental song learning. <i>Proc Biol Sci</i> . 285: 20180160
ChIP-Seq	SIRT1	Journal of Neuroimmune Pharmacology	2018	Nikki Bortell <i>et al.</i> Sirtuin 1-chromatin-binding dynamics points to a common mechanism regulating inflammatory targets in SIV infection and in the aging brain. <i>J Neuroimmune Pharmacol</i> . 13: 163
ChIP-Seq	H3K27me3	G3 - Genes, Genomics Genetics	2018	James Ferguson <i>et al.</i> PRC2 is dispensable in vivo for b-catenin-mediated repression of chondrogenesis in the mouse embryonic cranial mesenchyme. <i>G3</i> . 8: 491
MeDIP-Seq	5-Methylcytosine	Nature	2017	Rama S. Akondy <i>et al.</i> Origin and differentiation of human memory CD8 T cells after vaccination. <i>Nature</i> . 552: 362
ChIP-Seq	10 different Histone mods plus CTCF, RNA Pol II, & BRD4	Nature	2017	Elizabeth Stewart <i>et al.</i> Orthotopic patient-derived xenografts of paediatric solid tumours. <i>Nature</i> . 549: 96

Technique	Target	Journal	Year	Reference
ChIP-Seq	BACH2	Nature Communications	2017	Nicolas Hipp <i>et al.</i> IL-2 imprints human naive B cell fate towards plasma cell through ERK/ELK1-mediated BACH2 repression. <i>Nat Commun.</i> 8: 1443
ChIP-qPCR	H3K9Ac, RNA Pol II, ChREBP, PPARα	Cell Reports	2017	Alison Iroz <i>et al.</i> A specific ChREBP and PPARα cross-talk is required for the glucose-mediated FGF21 response. <i>Cell Reports.</i> 21: 403
ChIP-Seq	H3K79me2	Molecular Cancer Therapy	2017	Carly T. Campbell <i>et al.</i> Mechanisms of Pinometostat (EPZ-5676) treatment-emergent resistance in MLL-rearranged leukemia. <i>Mol Cancer Ther.</i> 16: 1669
Next-Gen Bisulfite-Seq, ChIP-Seq	KLF4	Leukemia	2017	Y Shen <i>et al.</i> Inactivation of KLF4 promotes T-cell acute lymphoblastic leukemia and activates the MAP2K7 pathway. <i>Leukemia.</i> 31: 1314
ChIP-Seq	H3K27ac, H3K4me1	Epigenomics	2017	John P. Thomson <i>et al.</i> Defining baseline epigenetic landscapes in the rat liver. <i>Epigenomics.</i> 9: 1503
ChIP-Seq	H3K36me3	The Journal of Clinical Investigation	2017	Huairui Yuan <i>et al.</i> Histone methyltransferase SETD2 modulates alternative splicing to inhibit intestinal tumorigenesis. <i>J Clin Invest.</i> 127: 3375
RIME	Estrogen Receptor, Progesterone Receptor	Cancer Research	2017	Jessica Finlay-Schultz <i>et al.</i> Breast cancer suppression by progesterone receptors is mediated by their modulation of estrogen receptors and RNA polymerase III. <i>Cancer Res.</i> 77: 4934
ChIP-Seq	EZH2, H3K27me3, H3K4me3	Molecular Cancer Therapy	2017	Dorothy Brach <i>et al.</i> EZH2 inhibition by tazemetostat results in altered dependency on B-cell activation signaling in DLBCL. <i>Mol Cancer Ther.</i> 16:2586
ChIP-Seq Spike-in	p300, H3K27Ac, Androgen Receptor	Cancer Research	2017	Lingyan Jin <i>et al.</i> Therapeutic targeting of the CBP/p300 bromodomain blocks the growth of castration-resistant prostate cancer. <i>Cancer Res.</i> 77: 5564
ChIP-qPCR	IRF5, NFκB (p65)	Journal of Biological Chemistry	2017	Leah Cushing <i>et al.</i> IRAK4 kinase activity controls Toll-like receptor induced inflammation through the transcription factor IRF5 in primary human monocytes. <i>J Biol Chem.</i> 292: 18689
ChIP-Seq	RNA Pol II	Journal of Biological Chemistry	2017	Leah A. Gates <i>et al.</i> Acetylation on histone H3 lysine 9 mediates a switch from transcription initiation to elongation. <i>J Biol Chem.</i> 292: 14456
ChIP-Seq ChIP-qPCR	Estrogen Receptor	Endocrinology	2017	Sylvia C. Hewitt <i>et al.</i> Role of ERα in mediating female uterine transcriptional Responses to IGF1. <i>Endocrinology.</i> 158: 2427
ATAC-Seq ChIP-Seq	H3K4me3, H3K9me3	Cancer Cell	2017	Gulfem Dilek Guler <i>et al.</i> Repression of stress-induced LINE-1 expression protects cancer cell subpopulations from lethal drug exposure. <i>Cancer Cell.</i> 32: 1
ChIP-Seq	BRD4, H3K9me3, RNA Pol II	Neuron	2017	Issam Aldiri <i>et al.</i> The dynamic epigenetic landscape of the retina during development, reprogramming and tumorigenesis. <i>Neuron.</i> 94: 550
ChIP-Seq	Lmx1b	Development	2017	Endika Haro <i>et al.</i> Lmx1b-targeted cis-regulatory modules involved in limb dorsalization. <i>Dev.</i> 11: 2009
ChIP-Seq	EZH2	Proceedings of the Natl. Academy of Sciences	2017	Yongfeng Liu <i>et al.</i> Epithelial EZH2 serves as an epigenetic determinant in experimental colitis by inhibiting TNFα-mediated inflammation and apoptosis. <i>Proc Natl Acad Sci.</i> 114: E3796
ChIP-Seq	Estrogen Receptor	Endocrinology	2017	Laurel A. Coons <i>et al.</i> DNA sequence constraints define functionally active steroid nuclear receptor binding sites in chromatin. <i>Endocrinology.</i> 10: 3212

Technique	Target	Journal	Year	Reference
ChIP-Seq	ETV5	Proceedings of the Natl. Academy of Sciences	2017	Zhen Zhang <i>et al.</i> Transcription factor Etv5 is essential for the maintenance of alveolar type II cells. <i>Proc Natl Acad Sci.</i> 114: E3903
ChIP-Seq	Androgen Receptor	Molecular and Cellular Endocrinology	2017	Claire Nash <i>et al.</i> Genome-wide analysis of AR binding and comparison with transcript expression in primary human fetal prostate fibroblasts and cancer associated fibroblasts. <i>Mol Cell Endocrinol.</i> 471: 1
ChIP-Seq Analysis	H3K27me3	Scientific Reports	2017	Brid O'Leary <i>et al.</i> Long non-coding RNA PARTICLE bridges histone and DNA methylation. <i>Sci Rep.</i> 7: 1790
ChIP-Seq	H3K36me2	The Journal of Clinical Investigation	2017	Ni Li <i>et al.</i> AKT-mediated stabilization of histone methyltransferase WHSC1 promotes prostate cancer metastasis. <i>J Clin Invest.</i> 127: 1284
ChIP-Seq ChIP-qPCR	EZH2, H3K27me3, H3K27ac	Cancer Cell	2017	Eric E. Gardner <i>et al.</i> Chemosensitive relapse in small cell lung cancer proceeds through an EZH2-SLFN11 axis. <i>Cancer Cell.</i> 31: 286
ChIP-Seq	TRIM28	Cell Reports	2017	Per Ludvik Brattas <i>et al.</i> TRIM28 controls a gene regulatory network based on endogenous retroviruses in human neural progenitor cells. <i>Cell Reports</i> 18: 1
ChIP-Seq	Cbx3/HPI	Scientific Reports	2017	Michael Sun <i>et al.</i> Cbx3/HPI γ deficiency confers enhanced tumor-killing capacity on CD8 ⁺ T cells. <i>Sci Rep.</i> 7: 42888
ChIP-Seq	H3K27Ac	Scientific Reports	2017	Yao Shen <i>et al.</i> Epigenetic and genetic dissections of UV-induced global gene dysregulation in skin cells through multi-omics analyses. <i>Sci Rep.</i> 7: 42646
ChIP-qPCR	SIRT1	Scientific Reports	2017	Jung-Yoon Yoo <i>et al.</i> KRAS activation and over-expression of SIRT1/BCL6 contributes to the pathogenesis of endometriosis and progesterone resistance. <i>Sci Rep.</i> 7: 6765
MeDIP-Seq	5-methylcytosine	Cerebral Cortex	2017	Daniela Grassi <i>et al.</i> Neuronal activity, TGF-Signaling and unpredictable chronic stress modulate transcription of Gadd45 family members and DNA methylation in the hippocampus. <i>Cereb Cortex.</i> 27: 4166
ChIP-Seq	SRC-2	PLoS Genetics	2017	Shruthy Suresh <i>et al.</i> SRC-2-mediated coactivation of anti-tumorigenic target genes suppresses MYC-induced liver cancer. <i>PLoS Genet.</i> 13: e1006650
ChIP-Seq	SLY1	Cell Death and Differentiation	2017	Charlotte Moretti <i>et al.</i> SLY regulates genes involved in chromatin remodeling and interacts with TBLIXR1 during sperm differentiation. <i>Cell Death Differ.</i> 24: 1029
ChIP-Seq	G34R	Acta Neuropathologica Communications	2017	Farhana Haque <i>et al.</i> Evaluation of a novel antibody to define histone 3.3 G34R mutant brain tumours. <i>Acta Neuropathol Commun.</i> 5: 45
ChIP-Seq	SALL4	Journal of Hematology & Oncology	2017	Lina Yang <i>et al.</i> The stem cell factor SALL4 is an essential transcriptional regulator in mixed lineage leukemia-rearranged leukemogenesis. <i>J Hematol Oncol.</i> 10: 159
ChIP-qPCR	EZH2	Mechanisms of Ageing and Development	2017	Khyobeni Mozhui <i>et al.</i> Conserved effect of aging on DNA methylation and association with EZH2 polycomb protein in mice and humans. <i>Mech Ageing Dev.</i> 162: 27
ChIP-Seq Spike-in	H3K27me3, H3K4me3, H3K9me3	PLoS One	2016	Brian Egan <i>et al.</i> An alternative approach to ChIP-Seq normalization enables detection of genome-wide changes in histone H3 lysine 27 trimethylation upon EZH2 inhibition. <i>PLoS One.</i> 11: e0166438

Technique	Target	Journal	Year	Reference
Low Cell ChIP-Seq	H3K27me3, CTCF	Nature Genetics	2016	Joachim Weischenfeldt <i>et al.</i> Pan-cancer analysis of somatic copy-number alterations implicates IRS4 and IGF2 in enhancer hijacking. <i>Nat Genet.</i> 49: 65
ChIP-Seq	OTX2, MITF, BRD4, H3K27Ac	Cancer Cell	2016	Pascal D. Johann <i>et al.</i> Atypical teratoid/rhabdoid tumors are comprised of three epigenetic subgroups with distinct enhancer landscapes. <i>Cancer Cell.</i> 29: 379
ChIP-Seq	GATA2, Progesterone Receptor	Cell Reports	2016	Cory A. Rubel <i>et al.</i> A Gata2-dependent transcription network regulates uterine progesterone responsiveness and endometrial function. <i>Cell Reports.</i> 17: 1414
ChIP-Seq	Androgen Receptor	Molecular Cell	2016	Boyu Zhang <i>et al.</i> Non-Cell-Autonomous Regulation of Prostate Epithelial Homeostasis by Androgen Receptor. <i>Mol Cell.</i> 63: 976
ChIP-Seq Spike-in	BRD4	The Journal of Clinical Investigation	2016	Mark L. McClelland <i>et al.</i> CCAT1 is an enhancer-templated RNA that predicts BET sensitivity in colorectal cancer. <i>J Clin Invest.</i> 126: 639
hMeDIP-Seq	5-Hydroxymethylcytosine	Nature Communications	2016	Carolina M. Greco <i>et al.</i> DNA hydroxymethylation controls cardiomyocyte gene expression in development and hypertrophy. <i>Nat Commun.</i> 7: 12418
ChIP-qPCR	Androgen Receptor	Cell Reports	2016	Karyn Schmidt <i>et al.</i> The lncRNA SLNCR1 mediates melanoma invasion through a conserved SRA1-like region. <i>Cell Rep.</i> 15: 2025
Low Cell ChIP-Seq	H3K4me3, H3K27Ac, H3K27me3	Biology of Reproduction	2016	Kazadi Mutoji <i>et al.</i> TSPAN8 expression distinguishes spermatogonial stem cells in the prepubertal mouse testis. <i>Biol Reprod.</i> 95: 17
ChIP-Seq	HDAC1, H3K27Ac	Cancer Discovery	2016	Anjali Mishra <i>et al.</i> Mechanism, consequences, and therapeutic targeting of abnormal IL15 signaling in cutaneous T-cell lymphoma. <i>Cancer Discov.</i> 6: 986
ChIP-Seq	V5-tagged DUX4	Human Molecular Genetics	2016	Jocelyn O. Eidahl <i>et al.</i> Mouse Dux is myotoxic and shares partial functional homology with its human paralog DUX4. <i>Hum Mol Genet.</i> 25: 4577
ChIP-Seq	H3K27me3	Molecular and Cellular Biology	2016	Kyung Hyun Yoo <i>et al.</i> Histone demethylase KDM6A controls the mammary luminal lineage through enzyme-independent mechanisms. <i>Mol Cell Biol.</i> 36: 2108
ChIP-Seq	H3K27Ac	Nature Genetics	2016	Pratiti Bandopadhyay <i>et al.</i> MYB-QKI rearrangements in angiocentric glioma drive tumorigenicity through a tripartite mechanism. <i>Nat Genet.</i> 48: 273
ChIP-Seq Spike-in	KDM5C	Cell Reports	2016	Shigeki Iwase <i>et al.</i> A mouse model of X-linked intellectual disability associated with impaired removal of histone methylation. <i>Cell Reports.</i> 14: 1
ChIP-Seq ChIP-qPCR	Wiz	eLIFE	2016	Luke Isbel <i>et al.</i> Wiz binds active promoters and CTCF-binding sites and is required for normal behaviour in the mouse. <i>Elife.</i> 5: e15082
ChIP-Seq	BRD4, H3K27Ac, H3K4me1, H3K27me3, HLX, LHX2, LMX1A	Nature	2016	Charles Y. Lin <i>et al.</i> Active medulloblastoma enhancers reveal subgroup specific cellular origins. <i>Nature.</i> 530: 57
ChIP-qPCR	H3K9me3 H3 pan-acetyl	Nucleic Acids Research	2016	Xiaoyu Chen <i>et al.</i> Probing the impact of chromatin conformation on genome editing tools. <i>Nucleic Acids Res.</i> 44: 6482
ChIP-Seq ChIP-qPCR	H3K9Ac	Stem Cells Translational Medicine	2016	Dalia Ali <i>et al.</i> Epigenetic library screen identifies abexinostat as novel regulator of adipocytic and osteoblastic differentiation of human skeletal (mesenchymal) stem cells. <i>Stem Cells Transl Med.</i> 5: 1036

Technique	Target	Journal	Year	Reference
ChIP-Seq	H3K27me3	Cancer Research	2016	John P. Thomson <i>et al.</i> Loss of Tet1 associated 5-hydroxymethylcytosine is concomitant with aberrant promoter hypermethylation in liver cancer. <i>Cancer Res.</i> 76: 3097
ChIP-Seq, ChIP-qPCR	H3K36me3	Oncogene	2016	Thai H. Ho <i>et al.</i> High-resolution profiling of histone h3 lysine 36 trimethylation in metastatic renal cell carcinoma. <i>Oncogene.</i> 35: 1565
ChIP-Seq	HDAC1, HDAC2, GATA2	PLoS One	2016	Jeffrey R. Shearstone <i>et al.</i> Chemical Inhibition of histone deacetylases 1 and 2 Induces fetal hemoglobin through activation of GATA2. <i>PLoS One.</i> 11: e0153767
ChIP-Seq	LL-37 (antimicrobial peptide)	Journal of Cancer	2016	Mindy Munoz <i>et al.</i> Antimicrobial peptide LL-37 participates in the transcriptional regulation of melanoma cells. <i>J Cancer.</i> 26: 2341
ChIP-Seq ChIP-qPCR	p53	Molecular Oncology	2016	Cheryl Chan <i>et al.</i> Global re-wiring of p53 transcription regulation by the hepatitis B virus X protein. <i>Mol Oncol.</i> 10: 1183
ChIP-Seq	PLZF, SALL4	Development	2016	Dawn L. Lovelace <i>et al.</i> The regulatory repertoire of PLZF and SALL4 in undifferentiated spermatogonia. <i>Development.</i> 143: 1893
ChIP-Seq	PLZF	PLoS Genetics	2016	Ramakrishna Kommagani <i>et al.</i> The Promyelocytic leukemia zinc finger transcription factor is critical for human endometrial stromal cell decidualization. <i>PLoS Genet.</i> 12: e1005937
ChIP libraries, sequencing and analysis	H3K27me3	Developmental Dynamics	2016	Oyvind Dahle <i>et al.</i> Inhibiting smad2/3 signaling in pluripotent mouse embryonic stem cells enhances endoderm formation by increasing transcriptional priming of lineage-specifying target genes. <i>Dev Dyn.</i> 245: 807.
ChIP-Seq	NR4A1	PLoS One	2016	Ryan P. Duren <i>et al.</i> Genome wide mapping of NR4A binding reveals cooperativity with ETS factors to promote epigenetic activation of distal enhancers in acute myeloid leukemia cells. <i>PLoS One.</i> 11: e0150450
ChIP-Seq	MTA1	Oncotarget	2016	Swati Dhar <i>et al.</i> Dietary pterostilbene is a novel MTA1-targeted chemopreventive and therapeutic agent in prostate cancer. <i>Oncotarget.</i> 7: 18469
ChIP-Seq	Myc-tagged Pet-1	The Journal of Neuroscience	2016	Steven C. Wyler <i>et al.</i> Pet-1 switches transcriptional targets postnatally to regulate maturation of serotonin neuron excitability. <i>J Neurosci.</i> 36:1758
ChIP-Seq	EGR2	Nature Communications	2015	Tomohisa Okamura <i>et al.</i> TGF- β 3-expressing CD4 ⁺ CD25 ⁻ LAG3 ⁺ regulatory T cells control humoral immune responses. <i>Nat Commun.</i> 6: 6329
ChIP-Seq	HA-tagged Shox2	Development	2015	Wenduo Ye <i>et al.</i> A common Shox-2-Nkx2-5 antagonistic mechanism primes the pacemaker cell fate in the pulmonary vein myocardium and sinoatrial node. <i>Development.</i> 142: 2521
ChIP-Seq	BA180/PBRM1	Molecular Cell	2015	Bokai Zhu <i>et al.</i> Coactivator-dependent oscillation of chromatin accessibility dictates circadian gene amplitude via REV-ERB loading. <i>Mol Cell.</i> 60: 769
ChIP-qPCR	EZH2, SUZ12, H3K-27me3, RNA Pol II	The EMBO Journal	2015	Stephen G. Dann <i>et al.</i> Reciprocal regulation of amino acid import and epigenetic state through Lat1 and EZH2. <i>EMBO J.</i> 34: 1773
ChIP-Seq	CCND1 (Cyclin D1)	Oncotarget	2015	Mathew C. Casimiro <i>et al.</i> Kinase-independent role of cyclin D1 in chromosomal instability and mammary tumorigenesis. <i>Oncotarget.</i> 6: 8524
ChIP-Seq	TRIM33	PLoS Genetics	2015	Luke Isbel <i>et al.</i> Trim33 binds and silences a class of young endogenous retroviruses in the mouse testis; a novel component of the arms race between retrotransposons and the host genome. <i>PLoS Genetics.</i> 11: e1005693

Technique	Target	Journal	Year	Reference
ChIP-Seq	Progesterone Receptor	Cell Reports	2015	Vincent J. Lynch <i>et al.</i> Ancient transposable elements transformed the uterine regulatory landscape and transcriptome during the evolution of mammalian pregnancy. <i>Cell Rep.</i> 10: 551
ChIP-Seq	EGR1	PLoS One	2015	Anthony A. Portale <i>et al.</i> Characterization of FGF23-dependent Egr-1 cistrome in the mouse renal proximal tubule. <i>PLoS One.</i> 10: e0142924
ChIP-Seq	FOXM1	Breast Cancer Research and Treatment	2015	Christina Yau <i>et al.</i> FOXM1 cistrome predicts breast cancer metastatic outcome better than FOXM1 expression levels or tumor proliferation index. <i>Breast Cancer Res Treat.</i> 154: 23
ChIP-Seq	BRD4	Molecular Cancer Therapy	2015	Ryan Lenhart <i>et al.</i> Sensitivity of small cell lung cancer to BET inhibition is mediated by regulation of ASCL1 gene expression. <i>Mol Cancer Ther.</i> 14: 2167
Next-Gen Bisulfite Seq	NA	The Journal of Clinical Investigation	2015	Coralie Hoareau-Aveilla <i>et al.</i> Reversal of microRNA-150 silencing disadvantages crizotinib-resistant NPM-ALK(+) cell growth. <i>J Clin Invest.</i> 125:3505
ChIP-Seq ChIP-qPCR	β -Arrestin-1	Cancer Research	2015	Smitha Pillai <i>et al.</i> β -arrestin-1 mediates nicotine-induced metastasis through E2F1 target genes that modulate epithelial-mesenchymal transition. <i>Cancer Res.</i> 75: 1009
ChIP-Seq	H3K9ac, RNA Pol II, TFIIB	Circulation: Heart Failure	2015	Danish Sayed <i>et al.</i> Acute targeting of general transcription factor IIB restricts cardiac hypertrophy via selective inhibition of gene transcription. <i>Circ Heart Fail.</i> 8: 138
ChIP-Seq	H3K4me3	Journal of the National Cancer Institute	2015	Cristian Taccioli <i>et al.</i> Repression of esophageal neoplasia and inflammatory signaling by anti-miR-31 delivery in vivo. <i>J Natl Cancer Inst.</i> 107: djv220
ChIP-Seq	FLAG-tagged Twist2	American Journal of Human Genetics	2015	Shannon Marchegiani <i>et al.</i> Recurrent mutations in the basic domain of TWIST2 cause ablepharon macrostomia and barber-say syndromes. <i>Am J Hum Genet.</i> 97: 99
ChIP-Seq	LHX6	Human Molecular Genetics	2015	Jeffry M. Cesario <i>et al.</i> Lhx6 and Lhx8 promote palate development through negative regulation of a cell cycle inhibitor gene, p57 ^{Kip2} . <i>Hum Mol Genet.</i> 24: 5024
ChIP-Seq	SOX9	Nucleic Acids Research	2015	Zhongcheng Shi <i>et al.</i> Context-specific role of SOX9 in NF-Y mediated gene regulation in colorectal cancer cells. <i>Nucleic Acids Res.</i> 43: 6257
ChIP-Seq	H3K27me3	Nucleic Acids Research	2015	Kyung Hyun Yoo <i>et al.</i> Loss of EZH2 results in precocious mammary gland development and activation of STAT5-dependent genes. <i>Nucleic Acids Res.</i> 43: 8774
ChIP-Seq	H3K4me1	BMC Biology	2015	Sara K. Harten <i>et al.</i> The recently identified modifier of murine metastable epialleles, Rearranged L-Myc Fusion, is involved in maintaining epigenetic marks at CpG island shores and enhancers. <i>BMC Biol.</i> 13: 21
ChIP-Seq	ZNF384	Molecular Endocrinology	2015	Paul Childress <i>et al.</i> Genome-wide mapping and interrogation of the Nmp4 anti-anabolic bone axis. <i>Mol Endocrinol.</i> 29: 1269
ChIP-Seq	FOXO1, RNA Pol II	Molecular Endocrinology	2015	Yasmin M. Vasquez <i>et al.</i> FOXO1 is Required for Binding of PR on IRF4, Novel Transcriptional Regulator of Endometrial Stromal Decidualization. <i>Mol Endocrinol.</i> 29: 421
ChIP-Seq	Progesterone Receptor, FOSL2	Endocrinology	2015	Erik C. Mazur <i>et al.</i> Progesterone receptor transcriptome and cistrome in decidualized human endometrial stromal cells. <i>Endocrinology.</i> 156: 2239
ChIP-Seq	H3K27me3, H3K4me1, H3K4me3, H3K27Ac, p300, RAR α	Immunity	2015	Chrysothemis C. Brown <i>et al.</i> Retinoic acid is essential for Th1 cell lineage stability and prevents transition to a Th17 cell program. <i>Immunity.</i> 42: 1

Technique	Target	Journal	Year	Reference
ChIP-Seq	SIX1	Cancer Cell	2015	Jenny Wegert <i>et al.</i> Mutations in the SIX1/2 pathway and the DROSHA/DGCR8 miRNA microprocessor complex underlie high-risk blastemal type Wilms tumors. <i>Cancer Cell.</i> 27: 298
ChIP-Seq	H3K4me3, H3K9me3	Epigenetics	2015	Nioka C. Chisholm <i>et al.</i> Histone methylation patterns in astrocytes are influenced by age following ischemia. <i>Epigenetics.</i> 10: 142
ChIP-Seq Data Analysis	KDM3A	PLoS Biology	2015	Mo-bin Cheng <i>et al.</i> Specific phosphorylation of histone demethylase KDM3A determines target gene expression in response to heat shock. <i>PLOS Biol.</i> 12: e1002026
ChIP-Seq	SRC1	Molecular Endocrinology	2014	Mounia Tannour-Louet <i>et al.</i> Hepatic SRC-1 activity orchestrates transcriptional circuitries of amino acid pathways with potential relevance for human metabolic pathogenesis. <i>Mol Endocrinol.</i> 28: 1707
ChIP-Seq	Fish H3K4me3	Molecular Ecology Resources	2014	Claudius F. Kratochwil <i>et al.</i> Mapping active promoters by ChIP-seq profiling of H3K4me3 in cichlid fish – a first step to uncover cis-regulatory elements in ecological model teleosts. <i>Mol Ecol Resour.</i> 15: 761
ChIP-Seq	GATA2, FOXA1, p300, CBP, SRC1, SRC2, SRC3, Androgen Receptor	Proceedings of the National Academy of Sciences	2014	Bin He <i>et al.</i> GATA2 facilitates steroid receptor coactivator recruitment to the androgen receptor complex. <i>Proc Natl Acad Sci.</i> 111: 18261
Next-Gen Bisulfite Seq	NA	Cell Reports	2014	Deepak Kumar <i>et al.</i> Fibroblast growth factor maintains chondrogenic potential of limb bud mesenchymal cells by modulating DNMT3A recruitment. <i>Cell Rep.</i> 8: 1419
ChIP-Seq	PPAR γ	Biology of Reproduction	2014	Kelsey E. Brooks <i>et al.</i> Peroxisome proliferator activator receptor gamma (PPAR γ) regulates conceptus elongation in sheep. <i>Biol Reprod.</i> 92: 42.
ChIP-Seq, LightSwitch™	PXR, p300, H3K4me1, H3K27ac	PLoS Genetics	2014	Robin P. Smith <i>et al.</i> Genome-wide discovery of drug-dependent human liver regulatory elements. <i>PLOS Genet.</i> 10: e1004648
ChIP-Seq	H3K9me3, H3K36me3, H3K27me3	Genome Research	2014	Goran Kungulovski <i>et al.</i> Application of histone modification-specific interaction domains as an alternative to antibodies. <i>Genome Res.</i> 24: 1842
ChIP-Seq ChIP-qPCR	SRF	Physiological Genomics	2014	Sharolyn V. Kawakami-Schulz. Serum response factor: positive and negative regulation of an epithelial gene expression network in the destrin mutant cornea. <i>Physiol Genomics.</i> 46: 277
ChIP-Seq ChIP-qPCR	SRC3	The Journal of Clinical Investigation	2014	Jun Qin <i>et al.</i> Androgen deprivation–induced NCoA2 promotes metastatic and castration-resistant prostate cancer. <i>J Clin Invest.</i> 124: 5013
ChIP-Seq	H3K27Ac	Journal of Biological Chemistry	2014	Andre Landin Malt <i>et al.</i> Identification of a face enhancer reveals direct regulation of LIM homeobox 8 (Lhx8) by wntless-int (WNT)/ β -catenin Signaling. <i>J Biol Chem.</i> 289: 30289
ChIP-Seq ChIP-qPCR	ZBTB20	Journal of Medical Genetics	2014	Malene B. Rasmussen <i>et al.</i> Neurodevelopmental disorders associated with dosage imbalance of ZBTB20 correlate with the morbidity spectrum of ZBTB20 candidate target genes. <i>J Med Genet.</i> 51: 605
ChIP-Seq	H3K9Ac, H3K27Ac	Nature	2014	Paul A. Northcott <i>et al.</i> Enhancer hijacking activates GFII family oncogenes in medulloblastoma. <i>Nature.</i> 511: 428
ChIP-Seq	H3K4me3, H3K27me3	Nature	2014	Volker Hovestadt <i>et al.</i> Decoding the regulatory landscape of medulloblastoma using DNA methylation sequencing. <i>Nature.</i> 510: 537
ChIP-qPCR	EZH2, Suz12, H3K-27me3	Molecular Cancer Therapy	2014	Sarah K. Knutson <i>et al.</i> Selective Inhibition of EZH2 by EPZ-6438 Leads to Potent Antitumor Activity in EZH2 Mutant Non-Hodgkin Lymphoma. <i>Mol Cancer Ther.</i> 13: 842

Technique	Target	Journal	Year	Reference
ChIP-Seq	RAR- α , RAR- β , RXR- α	American Journal of Physiology – Gastrointestinal and Liver Physiology	2014	Yuqi He <i>et al.</i> Biological functional annotation of retinoic acid alpha and beta in mouse liver based on genome-wide binding. <i>Am J Physiol Gastrointest Liver Physiol.</i> 307: G205
ChIP-Seq	H3K4me3, H3K27me3	Nature Communications	2014	Brian C. Belyea <i>et al.</i> Identification of renin progenitors in the mouse bone marrow that give rise to B-cell leukaemia. <i>Nat Commun.</i> 5: 3273
ChIP-Seq	Estrogen Receptor	Molecular Endocrinology	2014	Sylvia C. Hewitt <i>et al.</i> Novel DNA motif binding activity observed In vivo with an estrogen receptor a mutant mouse. <i>Mol Endocrinol.</i> 26: 899
ChIP-Seq	H3K27me3	Bioinformatics	2014	Yanxiao Zhang <i>et al.</i> PePr: A peak-calling prioritization pipeline to identify consistent or differential peaks from replicated ChIP-Seq data. <i>Bioinformatics.</i> 30: 2568
ChIP-Seq	H3K36me3	Genome Research	2014	Jeremy M. Simon <i>et al.</i> Variation in chromatin accessibility in human kidney cancer links H3K36 methyltransferase loss with widespread RNA processing defects. <i>Genome Res.</i> 24: 241
ChIP-Seq	SRC-2	Cell Reports	2014	Erin Stashi <i>et al.</i> SRC-2 Is an Essential Coactivator for Orchestrating Metabolism and Circadian Rhythm. <i>Cell Rep.</i> 6: 663
Next-Gen Bisulfite Seq	—	Nature Neuroscience	2014	Brian G. Dias <i>et al.</i> Parental olfactory experience influences behavior and neural structure in subsequent generations. <i>Nat Neurosci.</i> 17: 89
ChIP-Seq ChIP-qPCR	Zbtb20	Cerebral Cortex	2014	Jakob V. Nielsen <i>et al.</i> Zbtb20 Defines a Hippocampal Neuronal Identity Through Direct Repression of Genes That Control Projection Neuron Development in the Isocortex. <i>Cereb Cortex.</i> 24: 1216.
ChIP-qPCR	H3K9me3, H3	Nature Immunology	2013	Patrick M. Gubser <i>et al.</i> Rapid effector function of memory CD8+ T cells requires an immediate-early glycolytic switch. <i>Nat Immunol.</i> 14: 1064
ChIP-Seq	RNA Pol II	PLoS One	2013	Jonathan P. Riley <i>et al.</i> PARP-14 Binds Specific DNA Sequences to Promote Th2 Cell Gene Expression. <i>PLoS One.</i> 8: e83127
Next-Gen Bisulfite Seq	—	Clinical Cancer Research	2013	David S. Shames <i>et al.</i> Loss of NAPRT1 Expression by Tumor-Specific Promoter Methylation Provides a Novel Predictive Biomarker for NAMPT Inhibitors. <i>Clin Cancer Res.</i> 19: 6912
ChIP-Seq	Cyclin D1	Cancer Research	2013	Xiaoming Ju <i>et al.</i> Identification of a cyclin D1 network in prostate cancer that antagonizes epithelial-mesenchymal restraint. <i>Cancer Res.</i> 74: 508
ChIP-Seq	H3K27me3	Cancer Cell	2013	Sebastian Bender <i>et al.</i> Reduced H3K27me3 and DNA Hypomethylation Are Major Drivers of Gene Expression in K27M Mutant Pediatric High-Grade Gliomas. <i>Cancer Cell.</i> 24: 660
ChIP-Seq	ASXL1	Journal of Experimental Medicine	2013	Omar Abdel-Wahab <i>et al.</i> Deletion of Asxl1 results in myelodysplasia and severe developmental defects in vivo. <i>J Exp Med.</i> 210: 2641
ChIP-Seq	COUP-TFII (NR2F2)	Molecular Endocrinology	2013	Xilong Li <i>et al.</i> COUP-TFII Regulates Human Endometrial Stromal Genes Involved in Inflammation. <i>Mol Endocrinol.</i> 27: 2041
ChIP-Seq	Gfi1	Nature Immunology	2013	Chauncey J. Spooner <i>et al.</i> Specification of type 2 innate lymphocytes by the transcriptional determinant Gfi1. <i>Nat Immunol.</i> 14: 1229
ChIP-Seq	RNA Pol II	Journal of Allergy and Clinical Immunology	2013	Purvi Mehrotra <i>et al.</i> PARP-14 and its enzyme activity regulates Th2 differentiation and allergic airway disease. <i>J Allergy Clin Immunol.</i> 131: 52
ChIP-Seq	Progesterone Receptor	Molecular Endocrinology	2013	Ashlee R. Lain <i>et al.</i> Research Resource: Progesterone Receptor Targetome Underlying Mammary Gland Branching Morphogenesis. <i>Mol Endocrinol.</i> 27: 1743

Technique	Target	Journal	Year	Reference
ChIP-Seq	PPAR α	Chemico-Biological Interactions	2013	Patrick D. McMullen <i>et al.</i> A map of the PPAR α transcription regulatory network for primary human hepatocytes. <i>Chem Biol Interact.</i> 209: 104
ChIP-Seq	H3K36me3	Cancer Discovery	2013	Lynn Bjerke <i>et al.</i> Histone H3.3 Mutations Drive Pediatric Glioblastoma through Upregulation of MYCN. <i>Cancer Discov.</i> 3: 512
ChIP-Seq	p53	Physiological Genomics	2013	Yuwen Li <i>et al.</i> Genome-wide analysis of the p53 gene regulatory network in the developing mouse kidney. <i>Physiol Genomics.</i> 45: 948
ChIP-Seq	FOXA2	The FASEB Journal	2013	Justyna Filant <i>et al.</i> Integrated chromatin immunoprecipitation sequencing and microarray analysis identifies FOXA2 target genes in the glands of the mouse uterus. <i>FASEB J.</i> 28: 230
ChIP-Seq	p300	PLoS Genetics	2013	Aaron M. Wenger <i>et al.</i> The Enhancer Landscape during Early Neocortical Development Reveals Patterns of Dense Regulation and Co-option. <i>PLoS Genet.</i> 9: e1003728
ChIP-Seq	MYRF	PLoS Biology	2013	Helena Bujalka <i>et al.</i> MYRF is a membrane-associated transcription factor that autoproteolytically cleaves to directly activate myelin genes. <i>PLoS Biol.</i> 11: e1001625
ChIP-Seq ChIP-qPCR	RXR- α , RNA Pol II	PLoS One	2013	Astrid Kosters <i>et al.</i> Sexually Dimorphic Genome-Wide Binding of Retinoid X Receptor alpha (RXR α) Determines Male-Female Differences in the Expression of Hepatic Lipid Processing Genes in Mice. <i>PLoS One.</i> 8: e71538
ChIP-qPCR	NF κ B (p52 & p65)	Journal of Biological Chemistry	2013	Sarah L. Doyle <i>et al.</i> Nuclear factor κ B2 p52 protein has a role in antiviral immunity through I κ B kinase epsilon-dependent induction of Sp1 protein and interleukin 15. <i>J Biol Chem.</i> 288: 25066
ChIP-chip ChIP-qPCR	C/EBP β , RNA Pol II	Journal of Biological Chemistry	2013	Hana Vakili <i>et al.</i> CCAAT-enhancer-binding protein β (C/EBP β) and downstream human placental growth hormone genes are targets for dysregulation in pregnancies complicated by maternal obesity. <i>J Biol Chem.</i> 288: 22849
ChIP-Seq	KDM2B	The Journal of Clinical Investigation	2013	Alexandros Tzatsos <i>et al.</i> KDM2B promotes pancreatic cancer via Polycomb-dependent and -independent transcriptional programs. <i>J Clin Invest.</i> 123: 727
ChIP-Seq	Vitamin D Receptor	BMC Medicine	2013	Adam E. Handel <i>et al.</i> Vitamin D receptor ChIP-seq in primary CD4+ cells: relationship to serum 25-hydroxyvitamin D levels and autoimmune disease. <i>BMC Med.</i> 11: 163
ChIP-Seq	COUP-TFII (NR2F2)	Developmental Cell	2013	San-pin Wu <i>et al.</i> Atrial Identity Is Determined by a COUP-TFII Regulatory Network. <i>Dev Cell.</i> 25: 417
ChIP-Seq	H3K4me3	PLoS One	2013	Alex Gutteridge <i>et al.</i> Novel Pancreatic Endocrine Maturation Pathways Identified by Genomic Profiling and Causal Reasoning. <i>PLoS One.</i> 8: e56024.
ChIP-Seq	Glucocorticoid Receptor	PLoS One	2012	Marie-José C. van Lierop <i>et al.</i> Org 214007-0: A Novel Non-Steroidal Selective Glucocorticoid Receptor Modulator with Full Anti- Inflammatory Properties and Improved Therapeutic Index. <i>PLoS One.</i> 7: e48385
ChIP-Seq	KDM1A (LSD1), H3K4me2	Molecular and Cellular Biology	2012	Venugopalan D. Nair <i>et al.</i> Involvement of Histone Demethylase LSD1 in Short-Time-Scale Gene Expression Changes during Cell Cycle Progression in Embryonic Stem Cells. <i>Mol Cell Biol.</i> 32: 4861
ChIP-Seq	EZH2	Nature	2012	Michael T. McCabe <i>et al.</i> EZH2 inhibition as a therapeutic strategy for lymphoma with EZH2-activating mutations. <i>Nature.</i> 492: 108

Technique	Target	Journal	Year	Reference
Data Analysis	pTyr37 H2B	Nature Structural and Molecular Biology	2012	Kiran Mahajan <i>et al.</i> H2B Tyr37 phosphorylation suppresses expression of replication-dependent core histone genes. <i>Nat Struct Mol Biol.</i> 19: 930
ChIP-chip	MYBL2 (B-MYB)	PLoS One	2012	Ming Zhan <i>et al.</i> The B-MYB Transcriptional Network Guides Cell Cycle Progression and Fate Decisions to Sustain Self-Renewal and the Identity of Pluripotent Stem Cells. <i>PLoS One.</i> 7: e42350
ChIP-Seq	FLAG-BAP1, OGT, HCF1	Science	2012	Anwasha Dey <i>et al.</i> Loss of the Tumor Suppressor BAP1 Causes Myeloid Transformation. <i>Science.</i> 337: 1541
ChIP-Seq	RNA Pol II	Journal of Biological Chemistry	2012	Danish Sayed <i>et al.</i> Transcriptional regulation patterns revealed by high-resolution chromatin immunoprecipitation during cardiac hypertrophy. <i>J Biol Chem.</i> 288: 2546
ChIP-Seq	Androgen Receptor	BMC Genomics	2012	Zhou Zhu <i>et al.</i> Dose-dependent effects of small-molecule antagonists on the genomic landscape of androgen receptor binding. <i>BMC Genomics.</i> 13: 355
ChIP-Seq	Progesterone Receptor	Molecular Endocrinology	2012	Cory A. Rubel <i>et al.</i> Genome-Wide Profiling of Progesterone Receptor Binding in the Mouse Uterus. <i>Mol Endocrinol.</i> 26: 1428
ChIP-Seq	Estrogen Receptor, RNA Pol II	Molecular Endocrinology	2012	Sylvia C. Hewitt <i>et al.</i> Whole-Genome Estrogen Receptor α Binding in Mouse Uterine Tissue Revealed by ChIP-Seq. <i>Mol Endocrinol.</i> 26: 887
ChIP-qPCR	PDX1, TCF3	Chemistry & Biology	2012	Alice Kiselyuk <i>et al.</i> HNF4 α antagonists discovered by a high-throughput screen for modulators of the human insulin promoter. <i>Chem Biol.</i> 19: 806
ChIP-chip	ROR α	PLoS One	2012	Yongjun Wang <i>et al.</i> Regulation of p53 Stability and Apoptosis by a ROR Agonist. <i>PLoS One.</i> 7: e34921
ChIP-Seq	Nkx3.1	The Journal of Clinical Investigation	2012	Philip D. Anderson <i>et al.</i> Nkx3.1 and Myc crossregulate shared target genes in mouse and human prostate tumorigenesis. <i>J Clin Invest.</i> 122: 1907
ChIP-Seq	RNA Pol II	Cardiovascular Research	2012	Mingyue Han <i>et al.</i> GATA4 expression is primarily regulated via a miR-26b-dependent post-transcriptional mechanism during cardiac hypertrophy. <i>Cardiovasc Res.</i> 93: 645
ChIP-Seq	FXR (Farnesoid X Receptor)	The American Journal of Physiology: Gastrointestinal and Liver Physiology	2012	Julia Yue Cui <i>et al.</i> Bile acids via FXR initiate the expression of major transporters involved in the enterohepatic circulation of bile acids in newborn mice. <i>Am J Physiol Gastrointest Liver Physiol.</i> 302: G979
ChIP-Seq	Progesterone Receptor	PLoS One	2012	Ping Yin <i>et al.</i> Genome-wide progesterone receptor binding: cell type-specific and shared mechanisms in T47D breast cancer cells and primary leiomyoma cells. <i>PLoS One.</i> 7: e29021
ChIP-Seq	FLAG-CCND1 (Cyclin D1)	The Journal of Clinical Investigation	2012	Mathew C. Casimiro <i>et al.</i> ChIP sequencing of cyclin D1 reveals a transcriptional role in chromosomal instability in mice. <i>J Clin Invest.</i> 122: 833
ChIP-Seq	RBPJ	Stem Cells	2012	Yaochen Li <i>et al.</i> Genome-wide Analysis of N1ICD/RBPJ Targets In Vivo Reveals Direct Transcriptional Regulation of Wnt, SHH, and Hippo Pathway Effectors by Notch1. <i>Stem Cells.</i> 30: 741
ChIP-chip	Progesterone Receptor	The FASEB Journal	2011	Heather L. Franco <i>et al.</i> Epithelial progesterone receptor exhibits pleiotropic roles in uterine development and function. <i>FASEB J.</i> 26: 1218
ChIP-Seq	Msgn1	Nature Communication	2011	Ravindra B. Chalamalasetty <i>et al.</i> The Wnt3a/ β -catenin target gene Mesogenin1 controls the segmentation clock by activating a Notch signalling program. <i>Nat Commun.</i> 2: 390

Technique	Target	Journal	Year	Reference
ChIP-qPCR	IRF4, Myc	British Journal of Haematology	2011	Antonia Lopez-Girona <i>et al.</i> Lenalidomide downregulates the cell survival factor, interferon regulatory factor-4, providing a potential mechanistic link for predicting response. <i>Br J Haematol.</i> 154: 325
ChIP-chip	IRF4	Allergy	2011	Barrenas S. Bruhn <i>et al.</i> Increased expression of IRF4 and ETS1 in CD4+ cells from patients with intermittent allergic rhinitis. <i>Allergy.</i> 67: 33
ChIP-Seq	COBRA1 (Nelf-b)	Journal of Biological Chemistry	2011	Jianlong Sun <i>et al.</i> Genetic and Genomic Analyses of RNA Polymerase II-pausing Factor in Regulation of Mammalian Transcription and Cell Growth. <i>J Biol Chem.</i> 286: 36248
ChIP-Seq	Foxp3	Nucleic Acids Research	2011	Fabian Birzele <i>et al.</i> Next-generation insights into regulatory T cells: expression profiling and FoxP3 occupancy in Human. <i>Nucleic Acids Res.</i> 39: 7946
ChIP-qPCR	BRD4	Proceedings of the National Academy of Sciences	2011	Jennifer A. Mertz <i>et al.</i> Targeting MYC dependence in cancer by inhibiting BET bromodomains. <i>Proc Natl Acad Sci.</i> 108: 16669
ChIP-chip	C/EBP β	Journal of Molecular Endocrinology	2011	Aristides Lytras <i>et al.</i> Identification of functional CCAAT/enhancer-binding protein and Ets protein binding sites in the human chorionic somatomammotropin enhancer sequences. <i>J Mol Endocrinol.</i> 47: 179
ChIP-qPCR	Progesterone Receptor	Biochemical Pharmacology	2011	Matthew R. Yudt <i>et al.</i> Discovery of a novel mechanism of steroid receptor antagonism: WAY-255348 modulates progesterone receptor cellular localization and promoter interactions. <i>Biochem Pharmacol.</i> 82: 1709
ChIP-chip	RNA Pol II, H3K9Ac	Journal of Biological Chemistry	2011	Hong Hao <i>et al.</i> The Transcription Factor Neural Retina Leucine Zipper (NRL) Controls Photoreceptor-specific Expression of Myocyte Enhancer Factor Mef2c from an Alternative Promoter. <i>J Biol Chem.</i> 286: 34893
ChIP-chip	DAF-12	PLoS Genetics	2011	Daniel Hochbaum <i>et al.</i> DAF-12 Regulates a Connected Network of Genes to Ensure Robust Developmental Decisions. <i>PLoS Genet.</i> 7: e1002179
ChIP-qPCR	CTCF	DNA and Cell Biology	2011	Yan Jin <i>et al.</i> Enhancer-Blocking Activity Is Associated with Hypersensitive Site V Sequences in the Human Growth Hormone Locus Control Region. <i>DNA Cell Biol.</i> 30: 995
Bisulfite-Seq	Bisulfite Sequencing	PLoS One	2011	Michela Deleidi <i>et al.</i> Oct4-Induced Reprogramming Is Required for Adult Brain Neural Stem Cell Differentiation into Midbrain Dopaminergic Neurons. <i>PLoS ONE.</i> 6: e19926
MeDIP-chip	5-methylcytosine (5-mC)	BMC Biology	2011	Genevieve Lavoie <i>et al.</i> PKC isoforms interact with and phosphorylate DNMT1. <i>BMC Biol.</i> 9: 31
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor)	Toxicology and Applied Pharmacology	2011	K. Nadira De Abrew <i>et al.</i> Regulation of Bach2 by the aryl hydrocarbon receptor as a mechanism for suppression of B-cell differentiation by 2,3,7,8-tetrachlorodibenzo-p-dioxin. <i>Toxicol Appl Pharmacol.</i> 252: 150
ChIP-qPCR	PPAR γ	Investigative Ophthalmology and Visual Science	2011	Gerard A. Rodrigues <i>et al.</i> Differential Effects of PPAR γ Ligands on Oxidative Stress-Induced Death of Retinal Pigmented Epithelial Cells. <i>Invest Ophthalmol Vis Sci.</i> 52: 890
ChIP-chip	p73	Proceedings of the National Academy of Sciences	2011	Jennifer M. Rosenbluth <i>et al.</i> Differential regulation of the p73 cistrome by mammalian target of rapamycin reveals transcriptional programs of mesenchymal differentiation and tumorigenesis. <i>Proc Natl Acad Sci.</i> 108: 2076

Technique	Target	Journal	Year	Reference
ChIP-chip	LXR, RXR	Journal of Biological Chemistry	2011	Qi Shen <i>et al.</i> Liver X Receptor-Retinoic X Receptor (LXR-RXR) Heterodimer Cistrome Reveals Coordination of LXR and AP1 Signaling in Keratinocytes. <i>J Biol Chem.</i> 286: 14554
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor), MAF	Nature Immunology	2010	Lionel Apetoh <i>et al.</i> The aryl hydrocarbon receptor interacts with c-Maf to promote the differentiation of type 1 regulatory T cells induced by IL-27. <i>Nat Immunol.</i> 11: 854
ChIP-qPCR	p73	Molecular Cancer	2010	Kathryn G. Eby <i>et al.</i> ISG20L1 is a p53 family target gene that modulates genotoxic stress-induced autophagy. <i>Mol Cancer.</i> 9: 95
ChIP-Seq	Vitamin D Receptor	Genome Research	2010	Sreeram V. Ramagopalan <i>et al.</i> A ChIP-seq defined genome-wide map of vitamin D receptor binding: associations with disease and evolution. <i>Genome Res.</i> 20: 1352
ChIP-chip	O-GlcNAc, RNA Pol II	Proceedings of the National Academy of Sciences	2010	Dona C. Love <i>et al.</i> Dynamic O-GlcNAc cycling at promoters of <i>Caenorhabditis elegans</i> genes regulating longevity, stress, and immunity. <i>Proc Natl Acad Sci.</i> 107: 7413
ChIP-Seq	PXR (Pregnane X Receptor)	Toxicological Sciences	2010	Julia Yue Cui <i>et al.</i> Genetic and Epigenetic Regulation and Expression Signatures of Glutathione S-Transferases in Developing Mouse Liver. <i>Toxicol Sci.</i> 116: 32
ChIP-Seq	SRC3, RNA Pol II	Molecular Endocrinology	2010	Rainer B. Lanz <i>et al.</i> Global Characterization of Transcriptional Impact of the SRC-3 Coregulator. <i>Mol Endocrinol.</i> 24: 859
ChIP-qPCR	Maf1, Rpc39, mTOR, Brf1, Raptor	Journal of Biological Chemistry	2010	Boris Shor <i>et al.</i> Requirement of the mTOR Kinase for the Regulation of Maf1 Phosphorylation and Control of RNA Polymerase III-dependent Transcription in Cancer Cells. <i>J Biol Chem.</i> 285: 15380
ChIP-chip	AHR (Aryl Hydrocarbon Receptor)	Toxicological Sciences	2010	K. Nadira De Abrew <i>et al.</i> An Integrated Genomic Analysis of Aryl Hydrocarbon Receptor-Mediated Inhibition of B-Cell Differentiation. <i>Toxicol Sci.</i> 118: 454
ChIP-qPCR	Myc-tagged-Pet-1	Nature Neuroscience	2010	Chen Liu <i>et al.</i> Pet-1 is required across different stages of life to regulate serotonergic function. <i>Nat Neurosci.</i> 13: 1190
ChIP-chip	Androgen Receptor	Molecular Endocrinology	2010	Anastasia Wyce <i>et al.</i> The Androgen Receptor Modulates Expression of Genes with Critical Roles in Muscle Development and Function. <i>Mol Endocrinol.</i> 24: 1665
ChIP-qPCR	NFκB (p50 and p65)	Cardiovascular Research	2010	Alina G. Sofronescu <i>et al.</i> FGF-16 is a target for adrenergic stimulation through NF-κB activation in postnatal cardiac cells and adult mouse heart. <i>Cardiovasc Res.</i> 87: 102
ChIP-Seq	PXR (Pregnane X Receptor)	Nucleic Acids Research	2010	Julia Yue Cui <i>et al.</i> ChIPing the cistrome of PXR in mouse liver. <i>Nucleic Acids Res.</i> 38: 7943
ChIP-chip	RNA Pol II	Molecular Vision	2010	Padmaja Tummala <i>et al.</i> Temporal ChIP-on-Chip of RNA-Polymerase-II to detect novel gene activation events during photoreceptor maturation. <i>Mol Vis.</i> 16: 252
ChIP-chip	STAT4	The Journal of Immunology	2009	Seth R. Good <i>et al.</i> Temporal Induction Pattern of STAT4 Target Genes Defines Potential for Th1 Lineage-Specific Programming. <i>J Immunol.</i> 183: 3839
ChIP-Seq	p73	Molecular and Cellular Biology	2008	Jennifer M. Rosenbluth <i>et al.</i> A Gene Signature-Based Approach Identifies mTOR as a Regulator of p73. <i>Mol Cell Biol.</i> 28: 5951
ChIP-qPCR	AHR (Aryl Hydrocarbon Receptor)	Nature	2008	Francisco J. Quintana <i>et al.</i> Control of Treg and TH17 cell differentiation by the aryl hydrocarbon receptor. <i>Nature.</i> 453: 65