



STEM CELL EPIGENETICS

REGULATORS OF GENE TRANSCRIPTION

ACTIVE  MOTIF®

STEM CELL EPIGENETICS

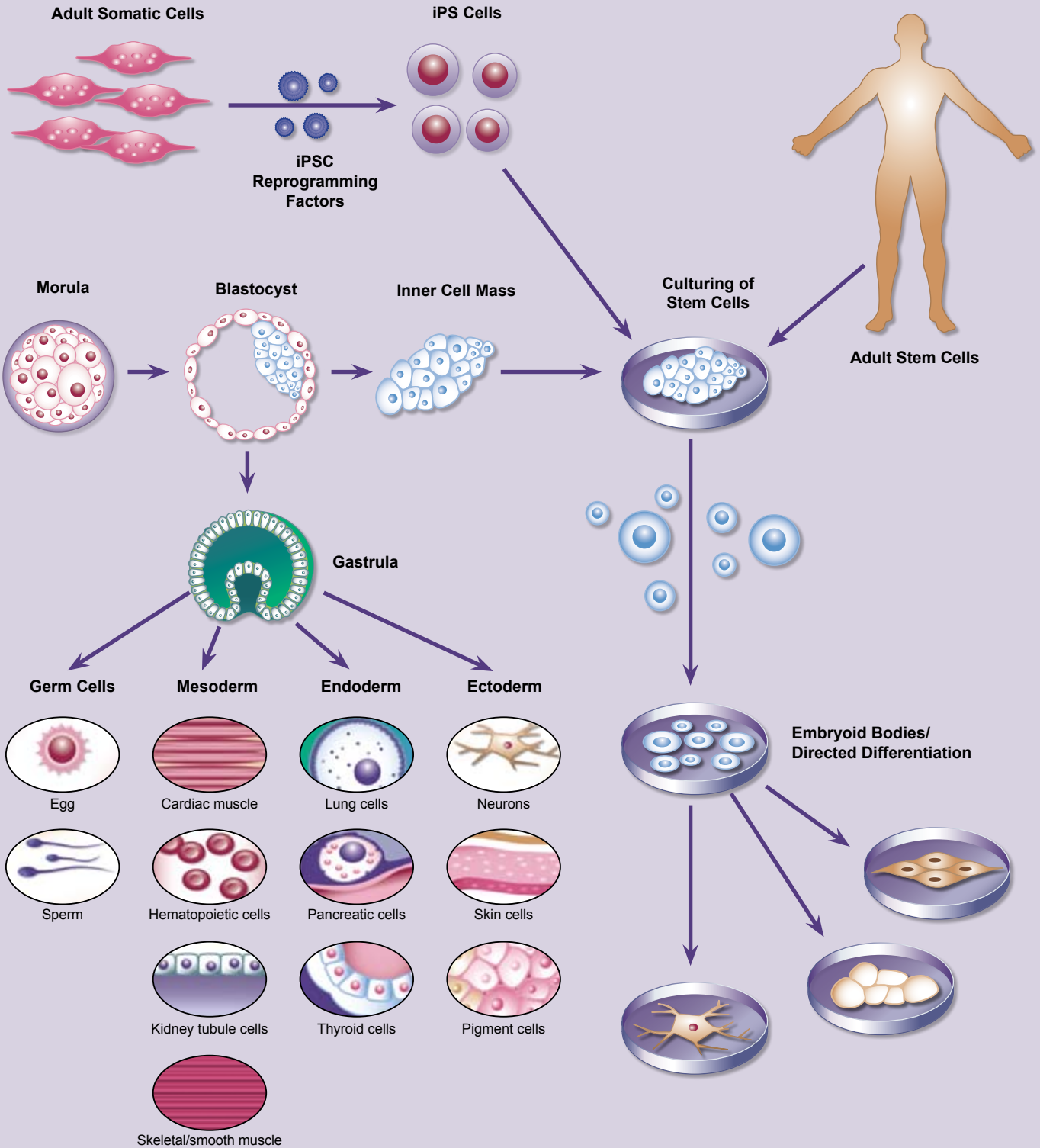
The two hallmark features of stem cells are pluripotency, the ability to differentiate into any mature cell type, and self-renewal, the capacity to undergo indefinite replicative cycles without losing stem cell identity. The general types of stem cells include embryonic stem cells (ESCs), adult stem cells and induced pluripotent stem cells (iPSCs). ESCs are derived from a population of cells in the blastocyst of a pre-implantation embryo called the inner cell mass that can differentiate into any cell type derivative of the three germ layers (endoderm, mesoderm and ectoderm). Adult stem cells can be found throughout the post-embryonic and adult organism, and function primarily in the maintenance, repair and regeneration of tissue and organs. iPSCs are pluripotent stem cells that are artificially derived when adult somatic cells are genetically reprogrammed to an embryonic stem cell phenotype.

A stem cell's decision whether to maintain its stem cell identity or differentiate into a specific cell type is ultimately determined by the outcome of the complex crosstalk that occurs between extracellular signaling pathways, transcriptional regulatory networks, chromatin remodeling complexes and non-coding RNAs. At the transcriptional level, pluripotency is largely controlled by the 'master regulators' OCT4, SOX2 and NANOG^{1,2}. These transcription factors form the core of the ESC transcriptional network and are essential for induction and maintenance of the stem cell phenotype^{1,3}. The primary cell signaling pathways involved in maintenance of pluripotency and self-renewal of ESCs are the WNT, TGF β /ACTIVIN/NODAL and FGFR pathways^{4,5}. These signaling pathways regulate the activity of OCT4, SOX2 and NANOG as well as auxiliary transcription factors and cofactors to drive the expression of stem cell-specific genes. During differentiation, other signaling pathways, such as BMP and NOTCH, signal the activation of the downstream expression of lineage-specific genes that promote the loss of pluripotency and diminish the proliferative potential of the cell^{4,5,6}.

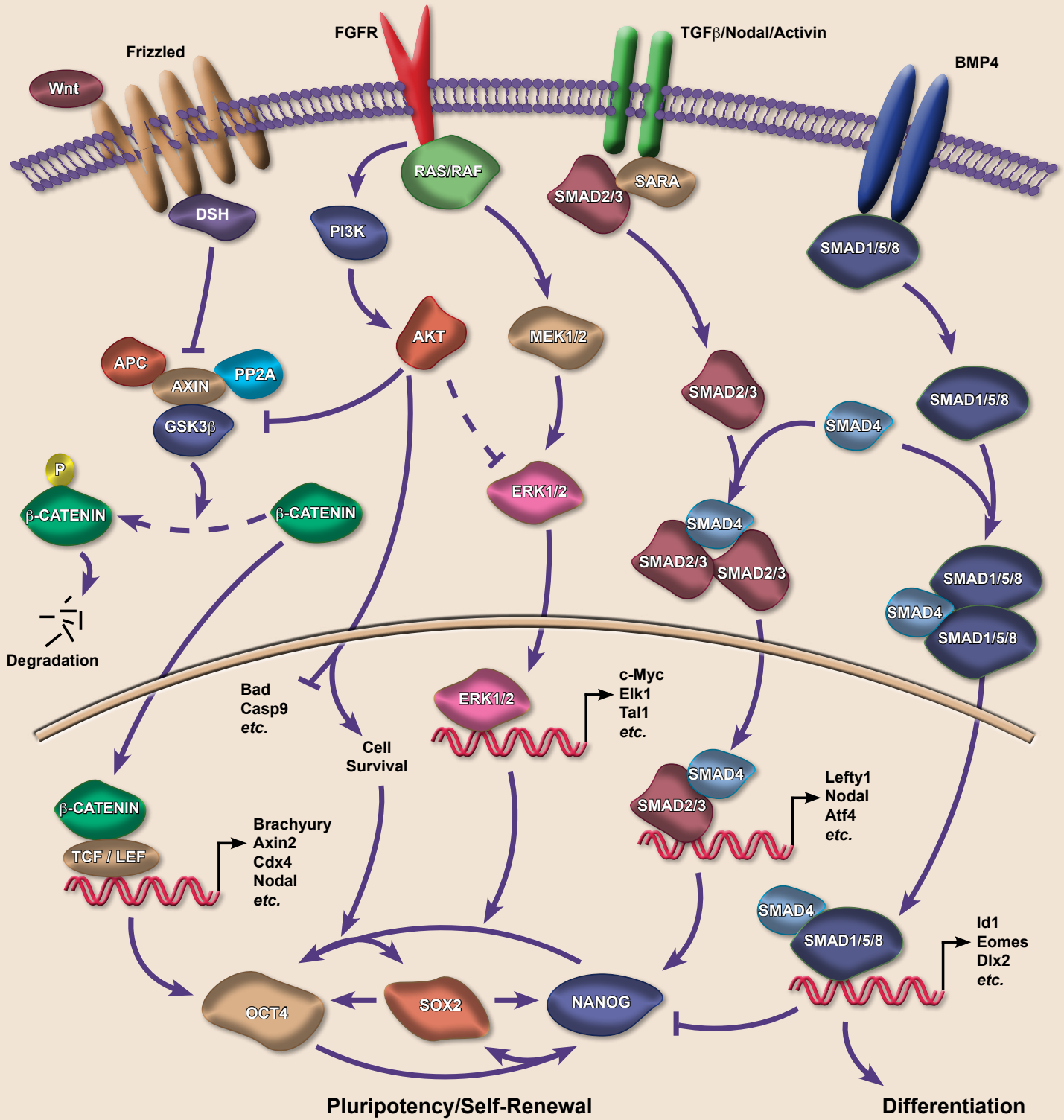
For a comprehensive understanding of the mechanisms that control self-renewal and pluripotency, it is essential to look beyond transcriptional networks towards the post-translational epigenetic events that modulate gene expression. These stochastic events set the epigenetic landscape within the cell by creating global changes that define regulatory networks, chromatin rearrangements, and the positioning of nuclear domains that determine the accessibility and transcriptional potential of underlying genes^{7,8,9}. Posttranslational epigenetic marks come in the form of acetylation, phosphorylation, methylation, citrullination and ubiquitination. By controlling the accessibility of DNA regulatory elements, these modifications modulate the interaction of transcription factor networks with other regulatory factors including transcriptional cofactors, chromatin remodeling proteins, histone modifiers, DNA methyltransferases and hydroxylases, and non-coding RNA regulators. This determines whether specific genes are actively transcribed, poised, or silenced at any given time^{7,8,10}. Having a multilayered transcriptional control mechanism serves as a system of checks and balances that allows fine-tuning and adaptability of the gene expression profile of a stem cell. This flexibility is the key to pluripotency, endowing stem cells the versatility to quickly modify gene expression in response to developmental and environmental cues, and to differentiate into essentially any cell type in the adult.

1. Boyer, L.A. *et al.* (2005) *Cell* **122**: 947-956.
2. Silva, J. and Smith, A. (2008) *Cell* **132**: 532-536.
3. Jaenisch, R. and Young, R. (2008) *Cell* **132**: 567-582.
4. Boiani, M. and Schöler, H.R. (2005) *Nat Rev Mol Cell Biol.* **6**: 872-884.
5. Okita, K. and Yamanaka, S. (2006) *Curr Stem Cell Res Ther.* **1**: 103-111.
6. Varga, A.C. and Wrana, J.L. (2005) *Oncogene* **24**: 5713-5721.
7. Young, R.A. (2011) *Cell* **144**: 940-954.
8. Han, J.W. and Yoon, Y.S. (2012) *Antioxid Redox Signal* **17**: 205-223.
9. Bilic, J. and Izpisua Belmonte, J.C. (2012) *Stem Cells* **30**: 33-41.
10. Chen, X. *et al.* (2008) *Cell* **133**: 1106-1117.

The Fate of Stem Cells



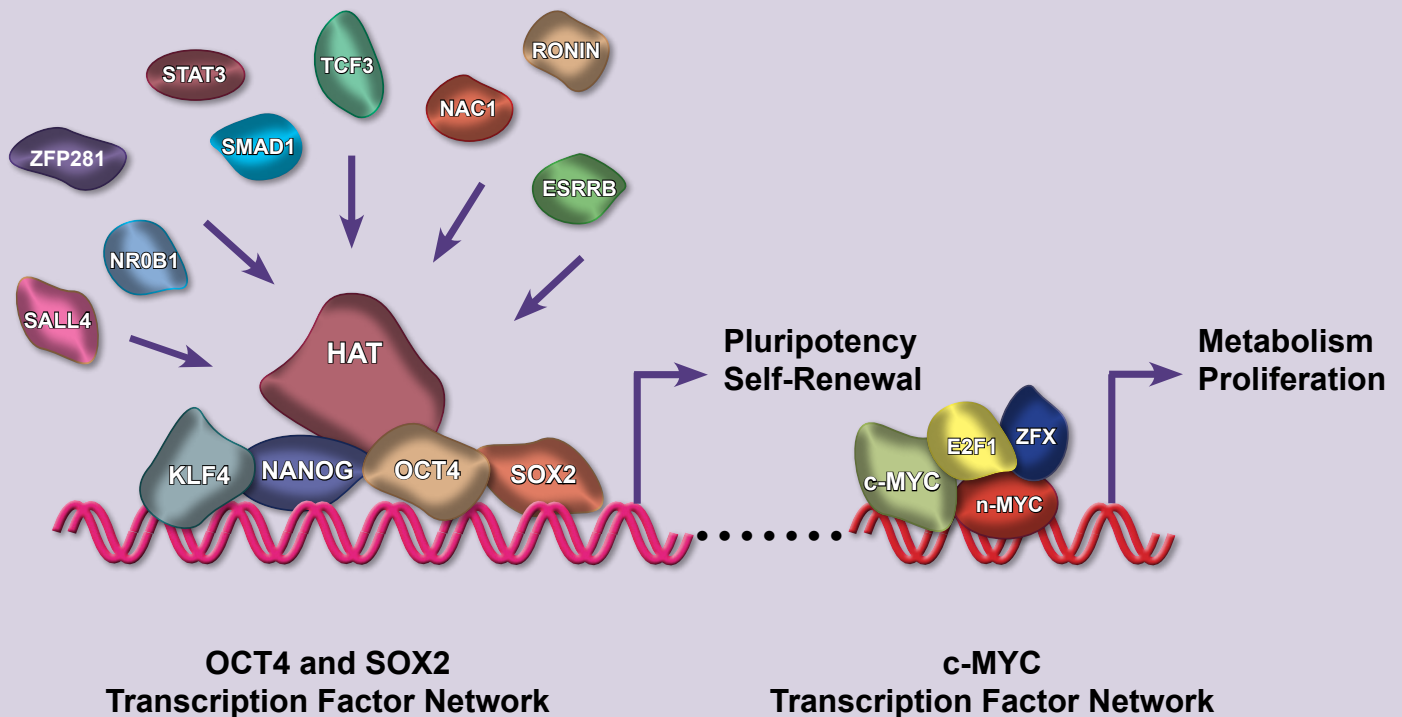
Cell Signaling Pathways Controlling Pluripotency and Self-Renewal



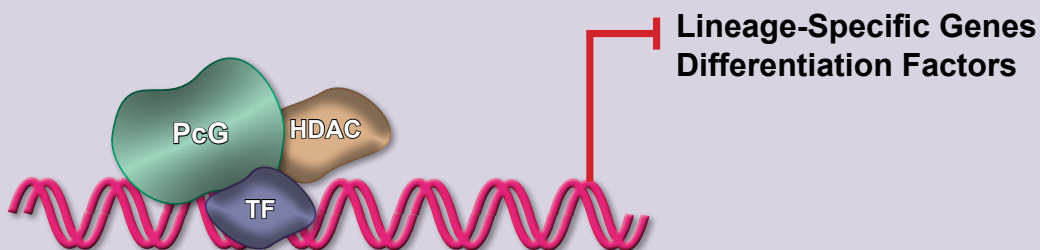
This diagram depicts the interaction between key intracellular signaling pathways regulating stem cell pluripotency and self-renewal.

Transcriptional Networks and Stem Cell Identity

ACTIVATION



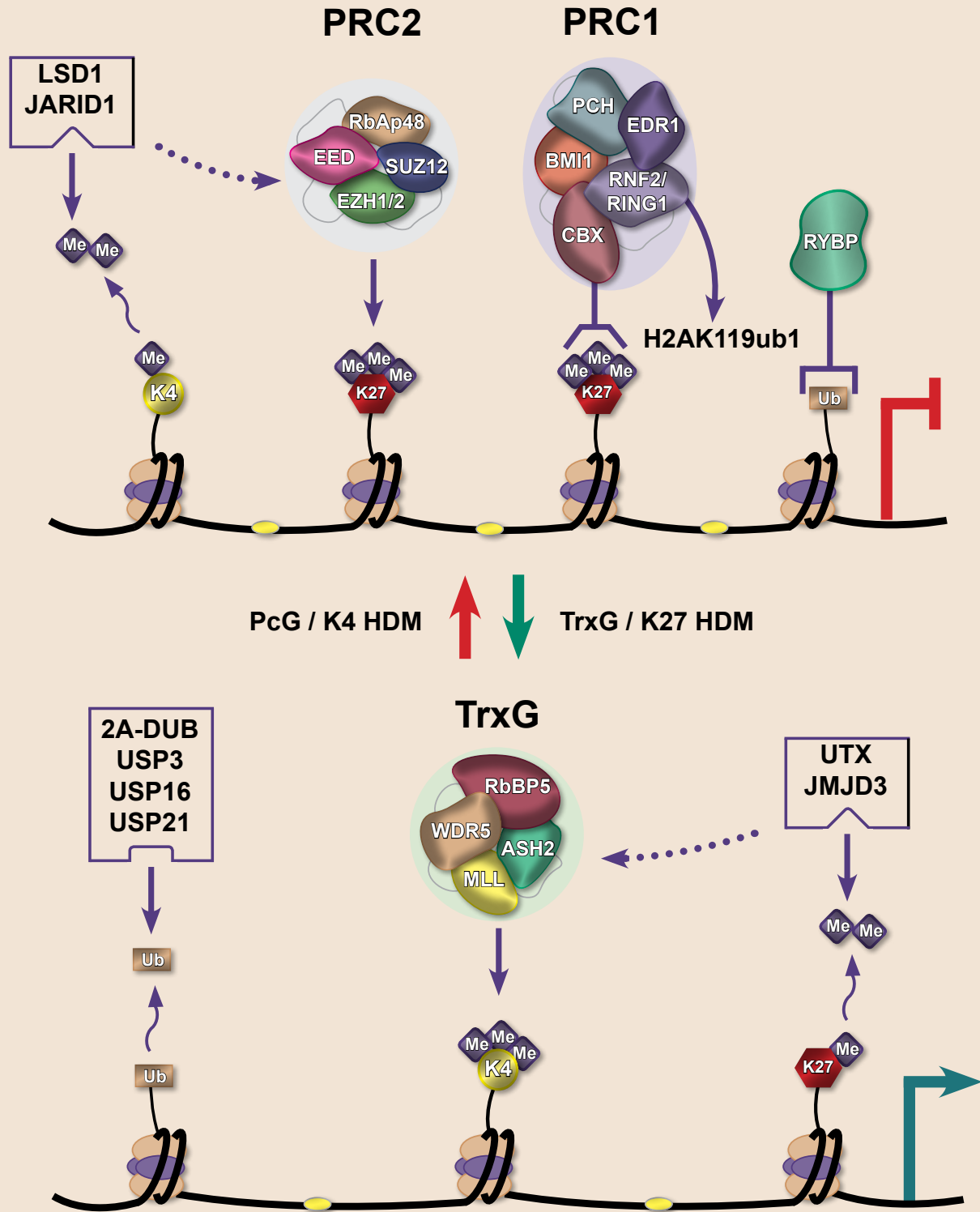
REPRESSION



A transcriptional network built around the master regulators OCT4, SOX2 and NANOG co-occupies promoter regions of genes regulating pluripotency and self-renewal. Recruitment of coactivators, such as histone acetyltransferases (HATs), signals transcriptional *activation*. In contrast, c-MYC functions to regulate the transcriptional efficiency of POLII, as well as to co-occupy and regulate genes involved in metabolism and proliferation. The subset of transcription factors (TF) co-occupying promoters varies in

response to epigenetic and intracellular signals. Low-level occupancy of promoters by reprogramming factors signals transcriptional *repression* and is accompanied by recruitment of corepressors, including the histone deacetylase (HDACs) complexes NuRD, NCoR/SMRT, SIN3A and REST, as well as Polycomb Group (PcG) proteins. In stem cells, this leads to repression of the somatic cell program.

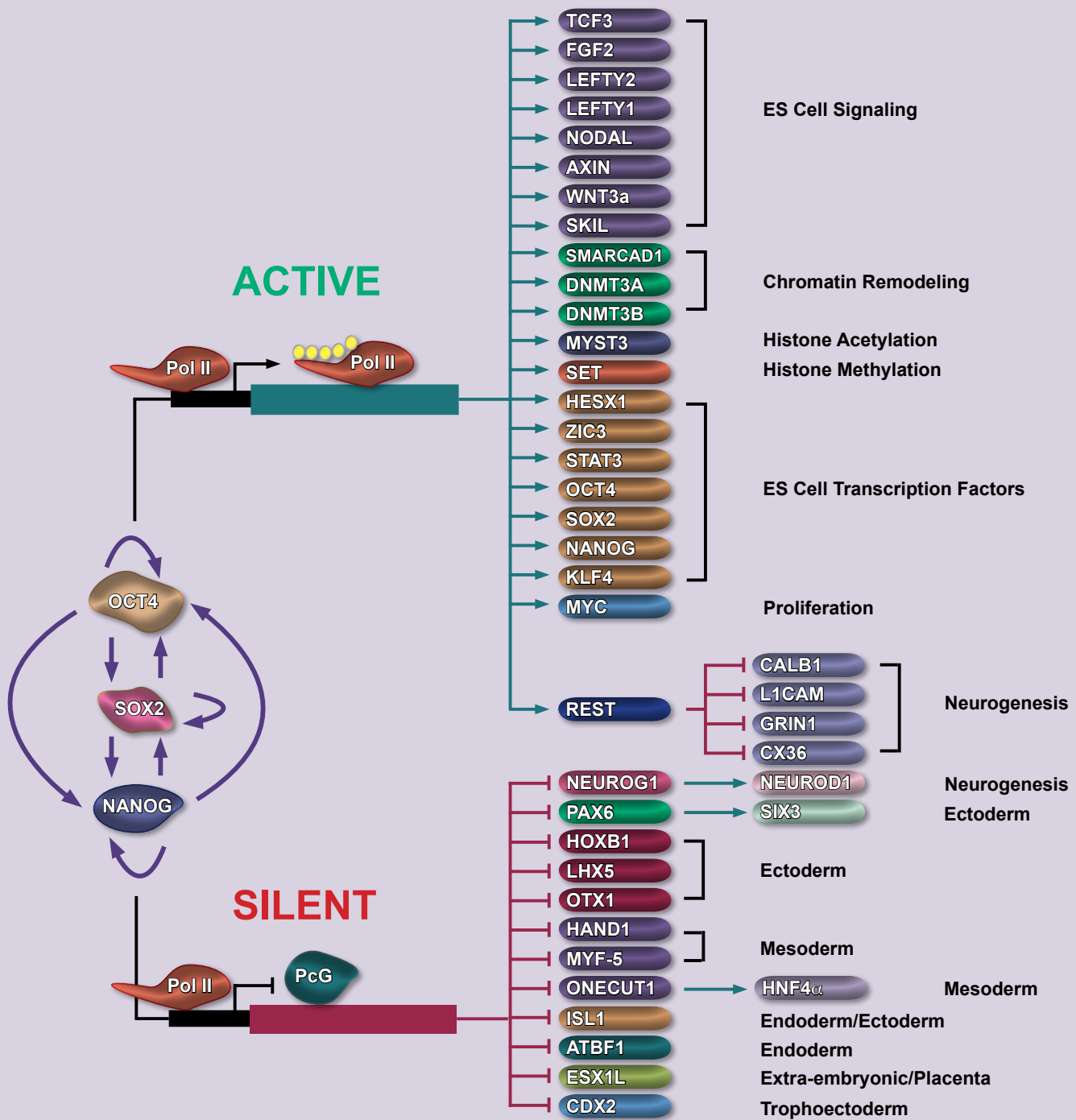
Transcriptional Regulation by Polycomb and Trithorax Groups



This diagram depicts the antagonistic relationship between the repressive Polycomb Group (PcG) and activating Trithorax Group (TrxG) proteins in regulating the transcriptional dynamics of embryonic stem cells. PcG gene silencing results from PRC2-mediated trimethylation of H3K27 by the histone methyltransferase (HMT) EZH1/2. H3K27me3 is recognized by the chromobox (CBX) subunit of PRC1, which leads to recruitment of PRC1 to the chromatin and the subsequent ubiquitination of H2AK119 via

RNF2/RING1 ubiquitin ligase. The RYBP repressor protein recognizes H2A mono-ubiquitination, contributing to transcriptional repression. Opposing this activity, TrxG recruitment of HMTs, such as MLL, mediate trimethylation of H3K4, leading to transcriptional activation and inhibition of PcG binding. The respective histone demethylases (HDMs) and deubiquitinating enzymes (DUB) are also shown.

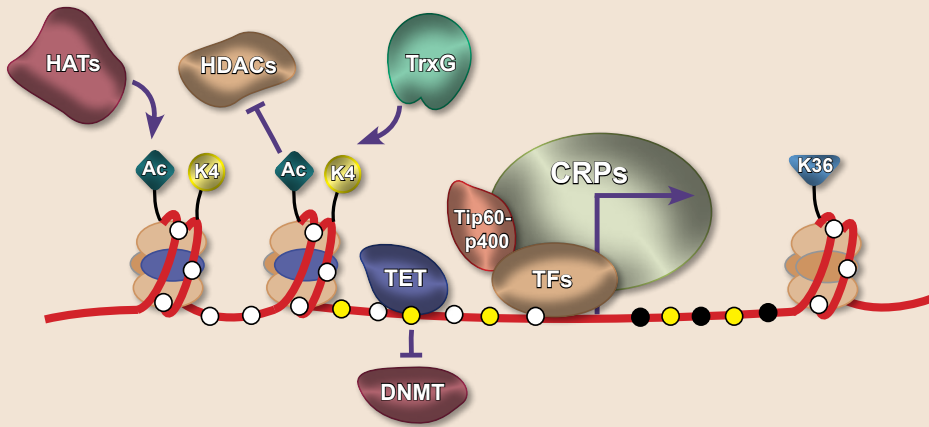
The Transcriptional Regulatory Circuitry and Nuclear Reprogramming



This transcriptional wiring diagram represents the core transcriptional regulatory circuitry in human embryonic stem cells based on expression data for OCT4, SOX2 and NANOG target genes. The core transcription factor interconnected autoregulatory loop is depicted on the left and the

activation and silencing of specific gene promoters and corresponding gene products (middle) are also shown. (Permission for use of this image was kindly provided by Dr. Rudolf Jaenisch, Professor of Biology at the Whitehead Institute for Biomedical Research at MIT).

Epigenetic Control of Chromatin Remodeling In Stem Cells

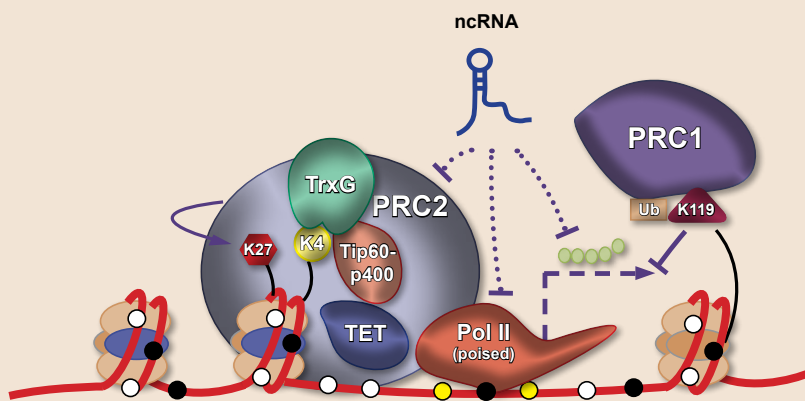


Active Chromatin (Euchromatin)

Signaling and stem cell maintenance proteins
FGF8, FGFR3, Lefty1, Inhba, Ezh1, etc.

Transcription factors
Stat3, Tcf3, Sall4, Esrrb, etc.

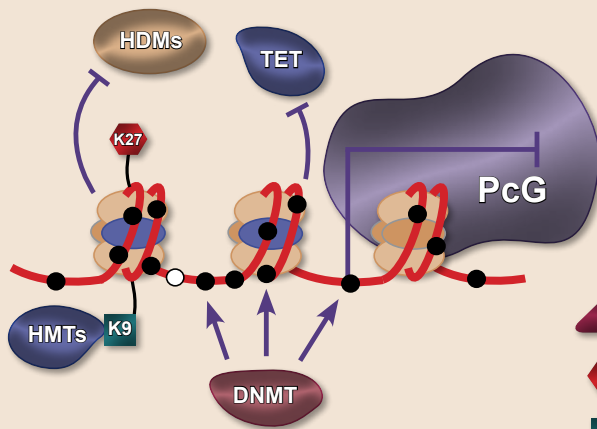
Regulators of proliferation/homeostasis
c-Myc, Tbx3, p53, GAPDH, Evi1, miR302-367, etc.



Poised Chromatin (Permissive/Repressed)

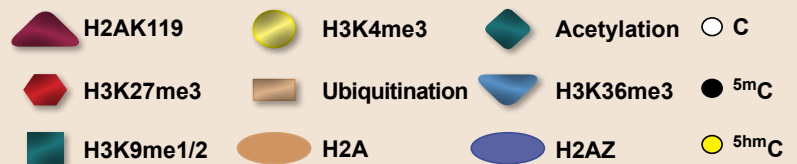
Core Transcription factors
Oct4, Sox2, Nanog, etc.

Early developmental regulators
Myf5, MyoD, Brachyury, Irx3, Pax6, etc.



Silent Chromatin (Heterochromatin)

Lineage specification factors
hCG, Adcy4, Anxa1, Actn1, Ebf1, etc.



In embryonic stem cells, H3K4 & H3K36 methylation and H3 & H4 acetylation are characteristic, active marks exclusively found within euchromatin. In addition, DNA regulatory elements of active genes are characterized by DNA hypomethylation. In combination, these modifications function to neutralize histone charges and recruit chromatin remodeling proteins (CRPs) that lead to unraveling of the chromatin structure, allowing access to the basal transcriptional machinery. In contrast, silenced genes are associated with condensed heterochromatin and are characteristically marked by H3K27 & H3K9 methylation and DNA hypermethylation.

'Bivalent domains,' where both activating H3K4 and repressive H3K27 methylation are present, mark genomic loci of early developmental regulators and HOX genes. These opposing marks silence genes while keeping them 'poised' for activation. Together, the cofactors and regulatory proteins effecting these epigenetic modifications define the chromatin landscape that dictates the expression profile of the cell.

TrxG, Trithorax group; PcG, Polycomb group; HATs, Histone acetyltransferases; HDACs, Histone deacetylases; TFs, Transcription factors; HDMs, Histone demethylases; DNMT, DNA methyltransferase; TET, Ten-eleven translocation enzymes; HMTs, Histone methyltransferases; PRC, Polycomb Repressive Complex.

Stem Cell Epigenetics Antibodies

Description	Applications	Cat. No.	Description	Applications	Cat. No.
MASTER REGULATORS					
c-Myc pAb	WB	39012	MyoD mAb	IF, WB	39991
KLF4 pAb	WB	39745	N-Myc pAb	WB	61185
LIN28A pAb	WB	61191	NKX2.5 pAb	WB	61267
Oct-4 pAb	WB	39811	Notch1 mAb	WB	61147
Sox2 pAb	Ch, IF, IHC, IP, WB	39823	Notch3 mAb	WB	61149
			NR2C2 pAb	WB	61279
			p53 pAb	Ch, EMSA	39334
			PAX7 mAb	IF, WB	39803
			PBX1b mAb	Ch, IHC, IP, WB	61165
			PDX1 pAb	WB	61289
			PLZF mAb	Ch, IF, WB	39987
			PP2A pAb	Ch, IP, WB	39192
			RNA pol II mAb	Ch IF, IP, WB	39097
			SALL4 pAb	WB	39957
			SIP1 mAb	IF, IHC, IP, WB	61095
			SMAD3 pAb	WB	61249
			Sox11 pAb	WB	61181
			Sp1 pAb	Ch, WB	39058
			STAT3 phospho Ser727 pAb	DB, WB	39613
			STAT3 phospho Tyr705 pAb	DB, WB	39595
			TAL-1 mAb	EMSA, IHC, WB	61259
			TAZ / WWTR1 pAb	WB	61265
			TCF7L1 / TCF3 pAb	WB	61125
			UTF1 pAb	WB	61253
			YY1 pAb	Ch, WB	39071
POLYCOMB GROUP			EPIGENETICS & CHROMATIN REMODELING		
BMI-1 mAb	Ch, IP	39993	5-Carboxylcytosine (5-caC) pAb	DB	61229
CBX8 pAb	WB	61237	5-Formylcytosine (5-fC) pAb	DB, IF	61223
EED mAb	IHC, WB	61203	5-Hydroxymethylcytosine (5-hmC) mAb	DB, MeDIP	39999
EZH2 mAb	Ch, IF, IP	39875	5-Hydroxymethylcytosine (5-hmC) pAb	DB, IF, IHC, MeDIP	39769
EZH2 phospho Thr345 pAb	DB, WB	61241	5-Methylcytosine (5-mC) mAb	DB, FACS, IHC, IP, MeDIP	39649
GCN5 mAb	ELISA, IF, WB	39975	Ago1/2/3 mAb	Ch, IF, IHC, IP, WB	39937
PCL2 mAb	WB	61153	BRG-1 mAb	IF, WB	39807
Phc1 mAb	IF, IP, WB	39723	BRM mAb	Ch, IF, WB	39805
Phc2 mAb	Ch, IF, IP	39661	CGBP pAb	WB	39203
Ring1B mAb	Ch, IF, IP, WB	39663	CHD1 pAb	Ch, WB	39729
Suz12 pAb	Ch, WB	39357	CoREST pAb	WB	39955
			Dicer mAb	WB	39817
			DNMT1 mAb	Ch, IHC, IP, WB	39204
TRITHORAX GROUP					
ASH2 pAb	IF, IP, WB	39099			
MLL pAb	Ch, WB	61295			
MLL1/HRX mAb	Ch, IP, WB	39829			
TRANSCRIPTION & REPROGRAMMING					
AKT1 phospho Ser473 mAb	WB	40902			
DAX1 / NR0B1 mAb	ICC, IF, IHC, IP, WB	39983			
FOXO1/FKHR pAb	WB	39629			
GATA-1 pAb	Ch, WB	39025			
GATA-4 pAb	WB	39893			
GATA-6 pAb	WB	61063			
GLI1 pAb	WB	61215			
Goosecoid pAb	WB	61121			
HNF-3β/ FOXA2 pAb	IHC, WB	39827			
HNF4A pAb	WB	61189			
HOXA9 pAb	WB	39825			
KLF5 pAb	WB	61099			
KLF6 mAb	IHC, WB	61297			
Myf-5 mAb	IF, WB	39801			

Description	Applications	Cat. No.	Description	Applications	Cat. No.
EPIGENETICS & CHROMATIN REMODELING, cont					
DNMT2 pAb	WB	39205	Histone H3 acetyl Lys27 pAb	Ch, ChC, ChS, DB, IF, WB	39133
DNMT3A mAb	Ch, IF, IHC, WB	39206	Histone H3 di/trimethyl Lys27 mAb	Ch, ChC, ChS, WB	39535
DNMT3B mAb	Ch, IF, IP, WB	39207	Histone H3 trimethyl Lys27 pAb	Ch, DB, ELISA, WB	39156
Drosha pAb	WB	39783	Histone H3 trimethyl Lys36 pAb	Ch, ChC, ChS, DB, WB	61101
HDAC1 pAb	Ch, ChC, ChS, WB	40967	HMGAI pAb	IF, WB	39615
Histone H2A pAb	WB	39209	HPI α mAb	Ch, ELISA, ICC, IF, IHC	39977
Histone H2AX pAb	IF, WB	39689	LSD1 pAb	Ch, ChC, ChS, IP, WB	39186
Histone H2AX phospho Ser139 pAb	DB, IF, WB	39117	MBD2 pAb	WB	39547
Histone H2A.Z pAb	Ch, WB	39113	MBD3 mAb	WB	39216
Histone H3, C-terminal pAb	Ch, WB	61277	MeCP2 mAb	Ch, IF, IHC, IP, WB	61291
Histone H3 mAb	Ch, IF, WB	39763	Mili / PiwiL2 mAb	IF, IHC, IP, WB	61143
Histone H3 monomethyl Lys4 pAb	Ch, ChC, ChS, DB, IF, WB	39297	MMSET / WHSC1 mAb	Ch, IF, IP, WB	39879
Histone H3 dimethyl Lys4 pAb	Ch, ChC, ChS, DB, WB	39141	PHF8 pAb	WB	39711
Histone H3 trimethyl Lys4 pAb	Ch, ChC, ChS, DB, IF, WB	39159	PRMT5 pAb	WB	61001
Histone H3 acetyl Lys9 pAb	Ch, ChC, ChS, DB, IF, WB	39917	PRMT6 pAb	WB	61003
Histone H3 dimethyl Lys9 pAb	Ch, DB, IF, WB	39239	SATB1 pAb	WB	39839
Histone H3 pan-methyl Lys9 pAb	DB, IF, WB	39241	SIN3A pAb	Ch, WB	39865
Histone H3 trimethyl Lys9 mAb	Ch, DB, IF, IP, WB	61013	SIRT1 mAb	IF, IP, WB	39353
Histone H3 trimethyl Lys9 pAb	Ch, ChC, ChS, DB, IF, WB	39161	SMRT / NCoR2 mAb	WB	61105
			SUV39H1 mAb	Ch, IP, WB	39785

Applications Key

Ch	Chromatin immunoprecipitation	IF	Immunofluorescence
ChC	ChIP-chip	IHC	Immunohistochemistry
ChS	ChIP-Seq	IP	Immunoprecipitation
DB	Dot blot	WB	Western blot
ICC	Immunocytochemistry		

For an up-to-date list of available stem cell antibodies, please visit www.activemotif.com/stemcellabs.