# **Epigenetics in Cancer**

**Adapted from Edition 3 of the Tocris Cancer Product Guide** To request a copy of the Tocris Cancer Product Guide, or to view the PDF, please visit www.tocris.com

Epigenetics can be defined as acquired changes in chromatin structure that arise independently of a change in the underlying DNA nucleotide sequence. Epigenetic modifications such as acetylation, methylation and ubiquitination can alter the accessibility of DNA to transcription machinery and therefore influence gene expression. The dysregulation of these epigenetic modifications has been shown to result in oncogenesis and cancer progression; the cell cycle and metastasis can be regulated by histone modification, DNA methylation and chromatin remodeling. Unlike genetic mutations, epigenetic alterations are considered to be reversible and thus make a promising therapeutic target.

# **Epigenetic Mechanism**

The fundamental unit of chromatin is the nucleosome, which consists of an octamer of the histone proteins H2A, H2B, H3 and H4 (two of each) tightly bound by DNA. Alterations in chromatin structure by post-translational modifications can regulate gene expression through the formation of heterochromatin or euchromatin, which usually repress or activate gene transcription, respectively.

Post-translational modifications include DNA methylation and the covalent methylation (Me) and acetylation (Ac) of histone tails. DNA methylation represses transcription by blocking the binding of transcription complexes to the gene promoter. The acetylation of histone tails usually loosens the DNA from around the nucleosomes, increasing the accessibility of gene promoters to transcription complexes, therefore promoting transcription. Histone tail methylation can repress or promote gene expression, depending on the site and extent of methylation, as well as the presence of other histone modifications in the vicinity. The pattern of these post-translational modifications on a nucleosome determines the transcriptional profile of nearby genes.

The functions of histone ubiquitination are less well understood. However increasing evidence points to an important role for this epigenetic modification in the DNA damage response.

# **Types of Epigenetic Modifiers**

Proteins that carry out these epigenetic modifications can be thought of as being either "writers", "readers" or "erasers".

• **Epigenetic writers** catalyze the addition of epigenetic marks onto either histone tails or the DNA itself.

• **Epigenetic reader domains** are effector proteins that recognize and are recruited to specific epigenetic marks. "Writer" and "eraser" enzymes may also contain such reader domains, leading to the coordination of "read-write" or "read-erase" mechanisms.

• Epigenetic erasers remove epigenetic marks to alter gene expression.

## **BRD** Inhibition Suppresses Tumor Growth and Metastasis

Bromodomains (BRDs) are epigenetic "readers" that selectively recognize acetylated lysine residues on histone protein tails. Of particular interest is the BET (bromodomain and extra-terminal) bromodomain family, which comprises the ubiquitously expressed proteins BRD2, BRD3, BRD4; and the testis-specific protein, BRDT. BET proteins are epigenome readers that play a key role at the interface between chromatin remodeling and transcriptional regulation, and are integral in the regulation of transcriptional elongation and the cell cycle. BRD4 influences mitotic progression and is a critical mediator of transcriptional elongation because it binds to transcriptional sites of genes expressed during the M/G<sub>1</sub> cell cycle transition. BRD4 increases expression of genes that promote growth by recruiting p-TEFb to mitotic chromosomes. Furthermore, it has been observed that BRD4 is significantly upregulated in both primary and metastatic melanomas. In vivo studies have shown that inhibition of BRD4 impairs tumor growth and metastasis.

Key BRD4 inhibitors include the potent, high affinity and selective, archetypical BET bromodomain inhibitor (+)-JQ1, which induces squamous cell differentiation and arrests tumor growth in BRD4-dependent carcinomas, including tumor growth in midline carcinoma cell xenograft models. I-BET 151 potently blocks recruitment of BRD3/4 to chromatin, inducing apoptosis and cell cycle arrest in MLL-fusion leukemia cell lines. This compound has also been shown to improve survival in rodent models of MLL-fusion leukemia.





Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals

www.tocris.com

### Products available from Tocris

14-3-3 Proteins

# Difopein

Aurora Kinases Anacardic acid, CCT 137690, Hesperadin, PF 03814735, SNS 314, TC-A 2317, TC-S 7010, ZM 447439 Bromodomains (BRD)

BAZ2-ICR, CP 203, GSK 2801, GSK 5959, I-BET 151, I-BRD9, (+)-JQ1, MS 436, NI 57, OF 1 Deubiquitinating Enzymes

P 22077, P005091, Spautin 1

DNA Methyltransferases (DNMT) 5-Azacytidine, Decitabine, EGCG, RG 108, SGI 1027, 6-Thioguanine, Zebularine

Histone Acetyltransferases (HAT) C 646, Garcinol, L002, NU 9056

Histone Deacetylases (HDAC) FK 228, MC 1568, MI 192, PCI 34051, SAHA, Trichostatin A, UF 010, Valproic acid

Histone Demethylases (KDM) GSK J1, GSK J2, GSK J4, IOX 1, JIB 04, RN 1, TC-E 5002

Lysine Methyltransferases (KMT) A 366, EPZ 004777, (*R*)-PFI 2, SGC 0946, UNC 0642, UNC 1999

MBT Domains UNC 1215, UNC 926

**Poly (ADP-ribose) Polymerase (PARP)** BYK 204165, EB 47, JW 55, MN 64, Nicotinamide, NU 1025, PJ 34, TC-E 5001, WIKI4, XAV 939

Protein Arginine Methyltransferases (PRMTs)

C 21, TC-E 5003

**Protein Ser/Thr Phosphatase** Ascomycin, GSK 2830371, Okadaic acid, Sanguinarine, Tautomycetin

Protein Tyrosine Phosphatase Alexidine dihydrochloride,

Sodium orthovanadate
RNA/DNA Polymerase

Mithramycin A, Thiolutin

Ubiquitin E3 Ligases SZL P1-41, SKPin C1, SMER 3

#### References:

Copeland *et al* (2009) Protein methyltransferases as a target class for drug discovery. *Nat. Rev. Drug Discov.* **8** 724. Dawson and Kouzarides (2012) Cancer

epigenetics: from mechanism to therapy. *Cell* **150** 12.

**Drouin** *et al* (2015) Structure enabled design of BAZ2-ICR, a chemical probe targeting the bromodomains of BAZ2A and BAZ2B. *J. Med* 

*Chem.* **58** 2553. **Muller** *et al* (2011) Bromodomains as therapeutic targets. *Expert Rev. Mol. Med.* **13** 

e29. Rodriguez-Paredes and Esteller (2011) Cancer epigenetics reaches mainstream oncology.

epigenetics reaches mainstream oncology. *Nat. Med.* **17** 330. **Suva** *et al* (2013) Epigenetic reprogramming in cancer. *Science* **339** 1567.

cancer. *Science* **339** 1567. **Virani** *et al* (2012) Cancer epigenetics: a brief review. *ILAR J.* **253** 359.

This poster is part of a series which compliments the Cancer Research Product Guide. For a copy of this poster or the Product Guide, please visit **www.tocris.com** © 2016 Tocris Cookson, Ltd. Tocris is a Bio-Techne brand