



Helping advance science,
one protein at a time.

NOV
2017

Mitochondrial Acetylation: Emerging Concepts and Therapeutic Potential

Related Publications

Research Tools

Meetings

Society for Neuroscience
2017

Booth # 2901

November 11-15,
Washington, D.C.

American Society for Cell
Biology 2017

Booth # 818

December 2-6,
Philadelphia, PA

Cytoskeleton Products

Actin Proteins
Activation Assays
Antibodies
ECM Proteins
ELISA Kits
G-LISA® Kits
Pull-down Assays
Motor Proteins
Small G-Proteins
Tubulin & FtsZ Proteins

Contact Us

P: 1 (303) 322.2254

F: 1 (303) 322.2257

E: cserve@cytoskeleton.com

W: cytoskeleton.com

Mitochondrial Acetylation: Emerging Concepts and Therapeutic Potential

General Overview of Lysine Acetylation in Health and Disease

Acetylation of the epsilon amino group of lysine residues (N^ε-acetylation) is an ancient, highly conserved post-translational modification (PTM)^{1,2} that links acetyl coenzyme A (acetyl-CoA) metabolism and cellular signaling. This occurs largely through the opposing activities of lysine acetyl transferases (KATs) and lysine deacetylases (KDACs)³. In humans, there are 3 major KAT families (GCN5, CBP/p300, and MYST) that all use acetyl-CoA as an essential cofactor to donate an acetyl group to target lysine residues. There are two KDAC families, the zinc-dependent histone deacetylases (HDAC1-11) and the NAD⁺-dependent sirtuins (SIRT1-7)¹⁻³.

It is well documented that acetylation of nuclear histones plays a major role in regulating chromatin compaction and transcriptional activity wherein acetylation favors a more open, transcriptionally active chromatin¹. Recent proteomic studies have identified over 4,500 non-histone proteins as targets of acetylation, thereby establishing lysine acetylation as a major global PTM. This PTM is present in many, if not all, cellular compartments, including the nucleus, cytoplasm, cell membrane, and mitochondria⁴. Functionally, reversible lysine acetylation has been shown to regulate enzyme activity⁴⁻⁶, protein-protein interactions⁷, and protein localization and stability^{8,9}. In addition, it plays critical regulatory roles in many cellular processes, including gene expression, cellular metabolism, apoptosis, cytoskeleton regulation, and membrane trafficking¹⁰.

In 2006, Vorinostat became the first FDA approved drug targeting histone epigenetics through HDAC inhibition for the treatment of cutaneous T-cell lymphomas (CTLC)¹¹. There are now four FDA approved HDAC inhibitors for the treatment of CTLC, peripheral T-cell lymphoma, and multiple myeloma¹², which clearly demonstrates the potential of targeting lysine acetylation for therapeutic intervention. Clinical success, coupled to the vast expansion in identification of non-histone acetyl lysine targeted proteins, has enhanced interest in exploring the regulatory functions and therapeutic potential of this PTM beyond epigenetics. The remainder of this Newsletter focuses on lysine acetylation in the mitochondria and its possible value as a therapeutic target.

Mitochondrial Lysine Acetylation

Lysine acetylation is abundant in mitochondria, with an estimated 700 proteins (63%) undergoing this modification, three fold higher than phospho-modifications in this organelle¹³. Elucidating the mechanisms that regulate mitochondrial acetylation and identifying potential therapeutic targets is a new and rapidly growing area of investigation¹⁰. In recent years, some general concepts have begun to emerge which are briefly outlined below:

- 1) Lysine acetylation is highly prevalent in enzymes of the major mitochondrial metabolic pathways and many of these proteins are hyper-acetylated⁴.
- 2) There is some sequence selectivity for mitochondrial protein acetylation as shown in a preference for negatively charged residues in the immediate vicinity of the acetylation site and a strong preference for hydrophobic residues at position +2⁴.
- 3) Mitochondrial protein acetylation changes rapidly in response to cellular nutrient availability or energy status and is likely a major mechanism regulating energy homeostasis¹⁴. It also appears that the levels of mitochondrial KATs and KDACs are tightly controlled by nutrient availability¹⁵. The link between pathological disruption of energy homeostasis and mitochondrial acetylation holds great therapeutic potential¹⁴.
- 4) Lysine acetylation has broadly been correlated with inhibition of mitochondrial enzyme activity, particularly in oxidative metabolism¹³. It is, however, noteworthy that of the 700 or so acetylated mitochondrial proteins, less than 40 have been characterized at the mechanistic level¹³, and there are cases where acetylation may either activate or inhibit the same protein depending on the tissue type^{6,15}.
- 5) Several groups have proposed a non-enzymatic mechanism for mitochondrial lysine acetylation, and the relative contribution of enzymatic vs non-enzymatic regulation remains to be determined¹⁶⁻¹⁸.
- 6) Low stoichiometry (< 1-5%) that translates to critical and regulated biological responses is a norm for PTMs and is one of the reasons that it is technically challenging to study these

News

Publications

Research Tools



Helping advance science,
one protein at a time.

Signal Seeker™ PRODUCTS

modifications^{19,20}. Low stoichiometry often reflects the highly dynamic nature and subcellular localization of the modified proteins and their role in cell signaling^{19,20}. Stoichiometry of acetylation in mitochondrial proteins appears to follow this trend^{23,24}. Understanding how sub-stoichiometric levels of PTMs can have significant effects on mitochondrial function remains an open question^{13,23,24}.

7) PTM crosstalk is recognized as a major cell regulatory mechanism²⁵. Mitochondrial acetylation has been shown to participate in PTM crosstalk as exemplified by the regulation of the pyruvate dehydrogenase complex via the hierarchical coordination of kinase, phosphatase, acetyltransferase, and deacetylase activities²¹.

8) Heart disease and heart failure are strongly correlated with increased acetylation of mitochondrial proteins in this organ²², and multiple reports have shown that the major mitochondrial deacetylase, Sirtuin 3 (Table 1), exerts a protective function against heart disease^{26,27}. Currently, there is intense interest in targeting SIRT3 and the mitochondrial acetylome for therapeutic intervention^{28,29}.

9) As shown in Table 1, the presence of mitochondrial KATs has only recently been described; however, they are considered to have strong therapeutic potential as targets for a variety of human diseases including cancer, heart disease, diabetes, and obesity (see references in Table 1).

Conclusion

Therapeutic exploitation of the mammalian mitochondrial acetylome offers great potential for the treatment of human disease; however, it is clear that much is still to be learned about the basic biology of this PTM beyond epigenetics. The role of lysine acetylation in the cytoplasm, plasma membrane, and other cellular compartments is also an area of high interest. In this regard, the development of an expanded toolset to aid in this research is urgently needed.

Table 1: Mammalian Mitochondrial Acetyltransferases and Deacetylases

Protein	Gene Name	Activity	Disease Linkage	Year Activity Described	Refs.
Acetyl-CoA acetyltransferase 1	ACAT1	acetyl transferase	C	2014	21
Males absent on the first	MOF	acetyl transferase	H	2016	30
General control of amino acid synthesis 5-like 1	GCN5L1	Lacks catalytic domain, essential component of acetyl transferase complex	H, O, D	2012	31,32
Sirtuin 3	SIRT3	deacetylase	H, O, D	2002	33
Sirtuin 4	SIRT4	Deacylase and very low deacetylase	D	2014	34
Sirtuin 5	SIRT5	deacylase	na	na	35

C-cancer, H-heart disease, O-obesity, D-diabetes

Signal-Seeker™ PTM Kits, Antibodies, Beads, Etc

Kit or Affinity Beads	Type	Size	Cat. #
Acetyl-Lysine Antibody Mouse Monoclonal (7B5A1)	MAB	2 x 100 ul 1 x 25 ul	AAC02 AAC02-S
Acetyl-Lysine Antibody Mouse Monoclonal (19C4B2.1)	MAB	2 x 100 ul 1 x 25ul	AAC03 AAC03-S
Acetyl-Lysine-HRP Antibody Mouse Monoclonal (19C4B2.1)	MAB	1 x 100 ul 1 x 25 ul	AAC03-HRP AAC03-HRP-S
Acetyl-Lysine Affinity Beads	Beads	4 x 500 ul	AAC04-beads
Signal-Seeker™ Acetyl-Lysine Enrichment Kit	Kit	30 assays 10 assays	BK163 BK163-S

References

- Kim G.W. and Yang X.J. 2011. Comprehensive lysine acetylomes emerging from bacteria to humans. *Trends Biochem. Sci.* **36**, 211-220.
- Weinert B.T. et al. 2011. Proteome wide mapping of the Drosophila acetylome demonstrates a high degree of conservation of lysine acetylation. *Sci. Signal.* **4**, ra48
- Choudhary C. et al. 2014. The growing landscape of lysine acetylation links metabolism and cell signaling. *Nat. Rev. Mol. Cell. Bio.* **15**, 536-550.
- Lundby A. et al. 2012. Proteomic analysis of lysine acetylation sites in rat tissues reveals organ specificity and subcellular patterns. *Cell Rep.* **2**, 419-431.
- Samant S.A. et al. 2014. SIRT3 deacetylates and activates OPA1 to regulate mitochondrial dynamics during stress. *Mol. Cell. Biol.* **34**, 807-819.
- Bharathi S.S. et al. 2013. Sirtuin 3 (SIRT3) protein regulates long chain acyl-CoA dehydrogenase by deacetylating conserved lysines near the active site. *J. Biol. Chem.* **288**, 33837-33847.
- Wei L. et al. 2013. Oroxylin A induces dissociation of hexokinase II from the mitochondria and inhibits glycolysis by SIRT3-mediated deacetylation of cyclophilin D in breast cancer. *Cell Death Dis.* **4**, e601.
- Inuzuka H. et al. 2012. Acetylation-dependent regulation of Skp2 function. *Cell.* **105**, 179-193
- De Boor S. et al. 2015. Small GTP-binding protein Ran is regulated by post translational lysine acetylation. *Proc. Natl. Acad. Sci. USA.* **112**, E3679-E3688.
- Drazic A. et al. 2016. The world of protein acetylation. *Biochim. Biophys. Acta.* **1864**, 1372-1401.
- Mann B.S. et al. 2007. FDA approval summary: vorinostat for treatment of advanced primary cutaneous T-cell lymphoma. *Oncologist.* **12**, 1247-1252.
- Yoon S. and Eom G.H. 2016. HDAC and HDAC inhibitor: from cancer to cardiovascular diseases. *Chonnam Med. J.* **52**, 1-11.
- Baeza J. et al. 2016. Mechanisms and dynamics of protein acetylation in mitochondria. *Trends Bio. Sci.* **41**, 231-244.
- Anderson K.A. and Hirschev M.D. 2012. Mitochondrial protein acetylation regulates metabolism. *Essays Biochem.* **52**, 23-35.
- Thapa D. et al. 2017. Acetylation of mitochondrial proteins by GCN5L1 promotes enhanced fatty acid oxidation in the heart. *Am. J. Physiol. Heart Circ. Physiol.* **313**, H265-H274.
- Wagner G.R. and Payne R.M. 2013. Widespread and enzyme-independent Nε-acetylation and Nε-succinylation of proteins in the chemical conditions of the mitochondrial matrix. *J. Biol. Chem.* **288**, 29036-29045.
- Davies M.N. et al. 2016. The acetyl group buffering action of carnitine acetyltransferase offsets macronutrient-induced lysine acetylation of mitochondrial proteins. *Cell Rep.* **14**, 243-254.
- Hosp F. et al. 2017. Lysine acetylation in mitochondria: from inventory to function. *Mitochondrion.* **33**, 58-71.
- Johnson H. and Eyers C.E. 2010. Analysis of post-translational modifications by LC-MS/MS. *Methods Mol. Biol.* **658**, 93-108.
- Thompson J. et al. 2012. https://genome.duke.edu/sites/default/files/ABRF_Poster_JWT_2012_PTMs.pdf.
- Fan J. et al. 2014. Tyr phosphorylation of PDP1 toggles recruitment between ACAT1 and SIRT3 to regulate the pyruvate dehydrogenase complex. *Mol. Cell.* **53**, 534-548.
- Stram A.R. et al. 2017. Progressive mitochondrial protein lysine acetylation and heart failure in a model of Friedreich's ataxia cardiomyopathy. *PLoS One.* **12**, e0178345.
- Weinert B.T. et al. 2014. Acetylation dynamics and stoichiometry in Saccharomyces cerevisiae. *Mol. Systems Biol.* **10**, 716.
- Weinert B.T. et al. 2015. Analysis of acetylation stoichiometry suggests that SIRT3 repairs nonenzymatic acetylation lesions. *EMBO J.* **34**, 2620-2632.
- Venne S.A. et al. 2014. The next level of complexity: crosstalk of posttranslational modifications. *Proteomics.* **14**, 513-524.
- Chen T.S. et al. 2015. Mouse SIRT3 attenuates hypertrophy-related lipid accumulation in the heart through the deacetylation of LCAD. *PLoS One.* **10**, e0118909.
- Koentges C. et al. 2015. SIRT3 deficiency impairs mitochondrial and contractile function in the heart. *Basic Res. Cardiol.* **110**, 36.
- Lee C.F. et al. 2015. Mitochondrion as a target for heart failure therapy. *Circ. J.* **79**, 1863-1870.
- Pillai V.B. et al. 2010. Exogenous NAD blocks cardiac hypertrophic response via activation of SIRT3-LKB1-AMP-activated kinase pathway. *J. Biol. Chem.* **285**, 3133-3144.
- Chatterjee A. et al. 2016. MOF acetyl transferase regulates transcription and respiration in mitochondria. *Cell.* **167**, 722-738.
- Scott I. et al. 2012. Identification of a molecular component of the mitochondrial acetyltransferase programme: a novel role for GCN5L1. *Biochem. J.* **443**, 655-661.
- Wang L. et al. 2017. GCN5L1 modulates cross-talk between mitochondria and cell signaling to regulate FoxO1 stability and gluconeogenesis. *Nat. Comm.* **8**, 523-534.
- Schwer B. et al. 2002. The human silent information regulator (Sir)2 homologue hSIRT3 is a mitochondrial nicotinamide adenine dinucleotide-dependent deacetylase. *J. Cell Biol.* **158**, 647-657.
- Mathias R.A. et al. 2014. Sirtuin 4 is a lipamidase regulating pyruvate dehydrogenase complex activity. *Cell.* **159**, 1615-1625.
- Tan M. et al. 2014. Lysine glutarylation is a protein posttranslational modification regulated by SIRT5. *Cell Metab.* **19**, 605-617.