



EpiCypher 2025 CONFERENCE

Biological and Clinical Frontiers in Epigenetics

Nov 2-7 | Grand Fiesta Americana Coral Beach | Cancún, MX



Organizers:

Mark Bedford | MD Anderson

Or Gozani | Stanford University Stanford

Brian Strahl | University of North Carolina Chapel Hill

Michael-Christopher Keogh | EpiCypher

Sunday, Nov 2 CANCUN TIME	Speakers & Event	Location
12:00pm – 7:00pm	Poster Boards Available for Set Up <i>Poster Numbers TBD</i>	Grand Coral Ballroom
2:00pm – 5:00pm	Conference Check-in <i>If you arrive later, please see the marketing table outside the Grand Coral Ballroom.</i>	Main Lobby
2:00pm – 4:00pm	Session # 1 – 4 Speakers: Presentation Practice <i>Please give your presentation materials to AV desk on a USB/USB-C. AV will have access to Mac and PC computers with Microsoft PowerPoint.</i>	Grand Coral Ballroom
4:00pm – 4:50pm	Coffee Bar Available	Grand Coral Foyer
4:50pm – 5:00pm	Welcome & Opening Remarks from Organizers Or Gozani, Brian Strahl, Mark Bedford & Michael-Christopher Keogh	Grand Coral Ballroom
5:00pm – 7:00pm	Session #1 Grand Coral Ballroom Chair: Liling Wan , University of Pennsylvania	
5:00pm – 5:20pm	Speaker 1: Johnathan Whetstine , Fox Chase Cancer Center <i>"The Face-off: Epigenetics and Genetics Talk to One another"</i>	
5:20pm – 5:40pm	Speaker 2: Greg Wang , Duke University <i>"Chromatin dysregulation driving oncogenesis"</i>	
5:40pm – 6:00pm	Speaker 3: Scott Rothbart , Van Andel Research Institute <i>"Revealing novel cancer therapeutic strategies through molecular studies of chromatin modification crosstalk"</i>	
6:00pm – 6:20pm	Speaker 4: Alexander Ruthenburg , University of Chicago <i>"Promiscuous RNA binding by WDR5 remodels the KMT2A (MLL1) histone methyltransferase complex to an inactive state"</i>	
6:20pm – 6:40pm	Speaker 5: Karim-Jean Armache , New York University <i>"Visualizing the Mechanisms of Epigenetic Inheritance and Regulation"</i>	
6:40pm – 7:00pm	Speaker 6: Lucas Farnung , Harvard Medical School <i>"Mechanisms of Transcription and Replication Through Chromatin"</i>	
7:30pm – 10:00pm	Welcome Reception (Dinner Included)	Sunset Terrace

Monday, Nov 3	Speakers and Events	Location
9:00am – 11:00am	Session # 2 Grand Coral Ballroom Chair: Erin Green , University of Maryland, Baltimore County	
9:00am – 9:20am	Speaker 7: Haitao Li , Tsinghua University <i>“A CMiDAC Deacylation Complex Pillars an Epigenetic Landscape to Prevent Heart Failure”</i>	
9:20am – 9:40am	Speaker 8: Pawel Mazur , MD Anderson Cancer Center <i>“Mouse models to investigate the targetable protein methylome in cancer pathogenesis”</i>	
9:40am – 10:00am	Speaker 9: Marc Timmers , Albert-Ludwigs-Universität Freiburg <i>“TGF-β activated transcription is controlled by a MLL4-JUNB feed-forward loop”</i>	
10:00am – 10:20am	Speaker 10: Kunal Rai , MD Anderson Cancer Center <i>“Precision epigenetic engineering to improve T cell therapy”</i>	
10:20am – 10:40am	Speaker 11: Andrew Stergachis , University of Washington <i>“A single-cell and haplotype-resolved view of normal and pathogenic gene regulation”</i>	
10:40am – 11:00am	Speaker 12: Ari Melnick , Josep Carreras Leukemia Research Institute <i>“Epigenetic and architectural mechanisms driving immune neoplasms”</i>	
11:00am – 11:20am	Coffee Break	Grand Coral Foyer
11:20am – 1:00pm	Session # 3 Grand Coral Ballroom Chair: Isaac Hilton , Rice University	
11:20am – 11:40am	Speaker 13: Danesh Moazed , Harvard Medical School <i>“Regulation of Suv39h read-write activity by H3K14 ubiquitination”</i>	
11:40am – 12:00pm	Speaker 14: Serena Sanulli , Stanford School of Medicine <i>“Allosteric regulation of the nucleosome core biases cell state”</i>	
12:00pm – 12:20pm	Speaker 15: Gregory Young , PacBio <i>“Multiomics with HiFi sequencing: new methylation callers, new insights”</i>	
12:20pm – 12:40pm	Speaker 16: Jessica Downs , The Institute of Cancer Research <i>“Functional profiling of SWI/SNF chromatin remodeling complexes through base editing screens”</i>	
12:40pm – 1:00pm	Speaker 17: Capucine Van Rechem , Stanford School of Medicine <i>“Replication stress and cholesterol biosynthesis as therapeutic vulnerabilities in Diffuse Midline Glioma”</i>	

1:00pm – 4:00pm	Free Time	
2:00pm – 4:00pm	Session # 5 – 7 Speakers: Presentation Practice <i>Please give your presentation materials to AV desk on a USB/USB-C. AV will have access to Mac and PC computers with Microsoft PowerPoint.</i>	Grand Coral Ballroom
4:00pm – 6:00pm	Session #4 Grand Coral Ballroom Chair: Mitchell Vollger , University of Utah	
4:00pm – 4:20pm	Speaker 18: Nick Young , Baylor College of Medicine <i>“A Needed Nomenclature for Nucleosomes and Histone Proteoform Mechanisms”</i>	
4:20pm – 4:40pm	Speaker 19: Neil Kelleher , Northwestern University <i>“Targeted enrichment of endogenous nucleosomes for characterization by tri-level”</i>	
4:40pm – 5:00pm	Speaker 20: Sriharsa Pradhan , New England Biolabs, Inc. <i>“RNA pol I inhibition remodels chromatin and alters nuclear architecture in mammalian cells”</i>	
5:00pm – 5:20pm	Speaker 21: Hong Wen , Van Andel Research Institute <i>“The histone acetylation reader ENL in development and cancer”</i>	
5:20pm – 5:40pm	Speaker 22: Yadira Soto-Feliciano , MIT <i>“Mechanisms of Transcriptional Regulation by Chromatin Adaptors in Cancer”</i>	
5:40pm – 6:00pm	Speaker 23: Liling Wan , University of Pennsylvania <i>“Chromatin-associated oncogenic condensates”</i>	
6:00pm – 8:00pm	Free Time	
8:00pm – 10:00pm	Poster Session 1 – Even Numbers <i>Drinks & Snacks will be provided. Please enjoy dinner beforehand.</i>	Grand Coral Ballroom/ Foyer

Tuesday, Nov 4	Speakers & Event	Location
9:00am – 11:00am	Session #5 Grand Coral Ballroom Chair: Kanishk Jain , University of Minnesota	
9:00am – 9:20am	Speaker 24: Jian Jin , Icahn School of Medicine at Mount Sinai <i>“Novel Degradators and Stabilizers of Epigenetic Targets”</i>	
9:20am – 9:40am	Speaker 25: Ryan Kruger , SK Life Science Labs <i>“Selective Degradation of p300 Results in a Potent Antitumor Effect in Prostate and CBP Loss of Function Cell Lines”</i>	
9:40am – 10:00am	Speaker 26: Yael David , Memorial Sloan Kettering Cancer Center <i>“Targeting a new metabolism-epigenetic axis towards improved CAR T therapy in Multiple Myeloma”</i>	
10:00am – 10:20am	Speaker 27: Irfan Asangani , University of Pennsylvania <i>“NSD2-Driven Enhancer Reprogramming as a Therapeutic Target in Prostate Cancer”</i>	
10:20am – 10:40am	Speaker 28: Ho Man Chan , AstraZeneca <i>“Epigenetic Drug Development: PRMT5 Inhibitor AZD3470 for MTAP-deficient cancers”</i>	
10:40am – 11:00am	Speaker 29: Karen Maegley , Pfizer <i>“Targeting the SETDB1 Methyltransferase with Covalent Inhibitors”</i>	
11:00am – 11:20am	Coffee Break	Grand Coral Foyer
11:20am – 1:00pm	Session #6 Grand Coral Ballroom Chair: Steve Josefowicz , Weill Cornell Medicine	
11:20am – 11:40am	Speaker 30: Stephane Richard , McGill University <i>“Understanding arginine methylation and RNA processes to design new treatments for cancer.”</i>	
11:40am – 12:00pm	Speaker 31: David Shechter , Albert Einstein College of Medicine <i>“PRMT5 as a Central Regulator of both RNA Processing and Histone Homeostasis”</i>	
12:00pm – 12:20pm	Speaker 32: Yanzhong (Frankie) Yang , City of Hope Cancer Center <i>“Arginine Methylation Facilitates the Recruitment of FACT complex to Chromatin to Maintain Active Transcription”</i>	
12:20pm – 12:40pm	Speaker 33: Y. George Zheng , University of Georgia <i>“Discovery of a non-catalytic function of PRMT1”</i>	

12:40pm – 1:00pm	Speaker 34: Sung Hee Baek , Seoul National University <i>“Epigenetic Code in Health and Disease”</i>	
1:00pm – 8:00pm	Free time <i>Please enjoy dinner before Trainee Session.</i>	
2:00pm – 3:30pm	Session # 8 – 9 Speakers: Presentation Practice <i>Please give your presentation materials to AV desk on a USB/USB-C. AV will have access to Mac and PC computers with Microsoft PowerPoint.</i>	Grand Coral Ballroom
8:00pm – 10:00pm	Session #7: C. David Allis Trainee Session Grand Coral Ballroom Session Introduction: Brian Strahl , UNC Chapel Hill Chair: Emily Bernstein , Icahn School of Medicine at Mount Sinai	
8:00pm – 8:15pm	Speaker 35: Abid Khan , UNC Chapel Hill <i>“Epigenetic Moonlighting by SETD2: From Chromatin to Cell Cycle”</i>	
8:15pm – 8:30pm	Speaker 36: Jaemin Kim , Seoul National University <i>“Regulation of HIF transcriptional programs promoting tumor progression in clear cell renal cell carcinoma”</i>	
8:30pm – 8:45pm	Speaker 37: Ben Mallory , University of Washington <i>“Principles and functional consequences of plasmid chromatinization in mammalian cells”</i>	
8:45pm – 9:00pm	Speaker 38: Alexia Martinez de Paz , Weill Cornell Medicine <i>“Signaling induced biophysical disruption of repressed chromatin domains drives immune cell fate”</i>	
9:00pm – 9:15pm	Speaker 39: Claudia Negrón-Lomas , McGill University <i>“The oncohistone H3.3 G34W impairs differentiation programs in development and cancer through epigenetic remodeling”</i>	
9:15pm – 9:30pm	Speaker 40: Deepak Saini , McGill University <i>“Divergent roles of DNA methylation, TRIM28, and p53 Surveillance in human embryonic and trophoblast stem cells”</i>	
9:30pm – 9:45pm	Speaker 41: Rebecca Smith , Fox Chase Cancer Center <i>“Histone Acetylation Differentially Modulates CTCF-CTCF Loops and Intra-TAD Interactions”</i>	
9:45pm – 10:00pm	Speaker 42: Adrienne Stampley , The University of Alabama at Birmingham <i>“The epigenome fine tunes transcriptional pause release across developmental potential”</i>	

Wednesday, Nov 5	Speakers and Events	Location
9:00am – 11:00am	Session #8 Grand Coral Ballroom Chair: Leah Gates , Case Western Reserve University	
9:00am – 9:20am	Speaker 43: Emily Bernstein , Icahn School of Medicine at Mount Sinai <i>"Mechanisms of chromatin remodeling mutations in cancer"</i>	
9:20am – 9:40am	Speaker 44: Chao Lu , Columbia University <i>"Genetic Dissection of Chromatin Regulatory Network"</i>	
9:40am – 10:00am	Speaker 45: Dan Levy , Ben-Gurion University <i>"Lysine Methylation Signaling at Chromatin"</i>	
10:00am – 10:20am	Speaker 46: Jacek Majewski , McGill University <i>"The Role of Histone 3 Lysine 36 Methyltransferases in Maintaining Epigenome Integrity"</i>	
10:20am – 10:40am	Speaker 47: Shehzad Z Sheikh , UNC Chapel Hill <i>"Chromatin Accessibility as a Predictors and Modifiers of Chronic Intestinal Inflammation"</i>	
10:40am – 11:00am	Speaker 48: Nicolas Reynoird , Université Grenoble Alpes <i>"RNF113A methylation dynamics controls Small Cell Lung Cancer sensitivity to alkylation damage"</i>	
11:00am – 11:20am	Coffee Break	Grand Coral Foyer
11:20am – 1:00pm	Session #9 Grand Coral Ballroom Chair: Capucine Van Rechem , Stanford School of Medicine	
11:20am – 11:40am	Speaker 49: Erin Green , University of Maryland, Baltimore County <i>"Pathways of chromatin regulation in hypoxia"</i>	
11:40am – 12:00pm	Speaker 50: Jessica Tyler , Weill Cornell Medicine <i>"Harnessed or hurtful heterologous engagement with the nucleosome acidic patch"</i>	
12:00pm – 12:20pm	Speaker 51: Xiaobing Shi , Van Andel Research Institute <i>"Epigenetic regulation of drug resistance in cancer"</i>	
12:20pm – 12:40pm	Speaker 52: Leah Gates , Case Western Reserve University <i>"Microbial metabolites regulate the chromatin landscape and gene expression"</i>	
12:40pm – 1:00pm	Speaker 53: Gerry Brien , University of Edinburgh <i>"Therapeutic targeting of oncogenic transcriptional biology in synovial sarcoma"</i>	

1:00pm – 4:00pm	Free Time	
2:00pm – 3:30pm	Session # 10 – 12 Speakers: Presentation Practice <i>Please give your presentation materials to AV desk on a USB/USB-C. AV will have access to Mac and PC computers with Microsoft PowerPoint.</i>	Grand Coral Ballroom
2:00pm – 3:00pm	Workshop 1: Advance your chromatin insights with CUT&RUN: Principles, Progress, and Possibilities Presenter: Tessa M. Firestone, <i>EpiCypher</i> Kingdom 3	
3:00pm – 4:00pm	Workshop 2: One assay, many insights: getting started with the amazingly capable long-read epigenomic assay, Fiber-seq Presenter: Bryan J. Venters, <i>EpiCypher</i> Kingdom 3	
4:00pm – 6:00pm	Session #10 Grand Coral Ballroom Chair: Nick Young , Baylor College of Medicine	
4:00pm – 4:20pm	Speaker 54: David Lahr , Foghorn Therapeutics, Inc. <i>“Proposal to use serial correlation to eliminate need for peak calling from analysis of epigenomic data”</i>	
4:20pm – 4:40pm	Speaker 55: Andrew Farmer , Takara Bio <i>“Dissecting gene regulatory programs through single-cell mapping of histone modifications”</i>	
4:40pm – 5:00pm	Speaker 56: Colin Ng , AtlasXomics <i>“Advancing Spatial Epigenomics: High-Resolution Mapping of Chromatin States in Intact Tissues”</i>	
5:00pm – 5:20pm	Speaker 57: Andres Blanco , University of Pennsylvania <i>“Targeting epigenetic regulators to induce therapeutic differentiation in acute myeloid leukemia.”</i>	
5:20pm – 5:40pm	Speaker 58: Xiaodong Cheng , MD Anderson Cancer Center <i>“Updated understanding of the protein–DNA recognition code used by C2H2 zinc finger proteins”</i>	
5:40pm – 6:00pm	Speaker 59: Makiko Iwafuchi , Cincinnati Children's Hospital Medical Center <i>“PRDM1 establishes super-poised enhancers in a pioneer factor dose-dependent manner to safeguard cell identity”</i>	
6:00pm – 8:00pm	Free Time	
8:00pm – 10:00pm	Poster Session 2- Odd Numbers <i>Drinks & Snacks will be provided. Please enjoy dinner beforehand.</i>	Grand Coral Ballroom/ Foyer

Thursday, Nov 6	Speakers and Events	Location
9:00am – 11:00am	Session #11 Grand Coral Ballroom Chair: Dan Levy , Ben-Gurion University	
9:00am – 9:20am	Speaker 60: Ernesto Guccione , Icahn School of Medicine at Mount Sinai <i>“Targeting solid tumors addiction to the WNT pathway”</i>	
9:20am – 9:40am	Speaker 61: Nada Jabado , McGill University <i>“The Fire Exit: Redirecting H3K27M-Stalled OPC States in Gliomas via ASCL1 Hijacking”</i>	
9:40am – 10:00am	Speaker 62: Christopher Hartl , Epigenome Technologies <i>“Practical Single-Cell CUT&Tag: Antibodies, Spike-Ins, and the 12x Rule”</i>	
10:00am – 10:20am	Speaker 63: Robert Schneider , Helmholtz Munich <i>“New modulators of chromatin and gene expression”</i>	
10:20am – 10:40am	Speaker 64: François Fuks , Université Libre de Bruxelles <i>“RNA modifications in Health and Disease”</i>	
10:40am – 11:00am	Speaker 65: Winston Timp , Johns Hopkins University <i>“Increasing Versatility of Single Molecules Sequencing for Epigenetics”</i>	
11:00am – 11:20am	Coffee Break	Grand Coral Foyer
11:20am – 1:00pm	Session #12 Grand Coral Ballroom Chair: Nicolas Reynoird , Université Grenoble Alpes	
11:20am – 11:40am	Speaker 66: Chen Davidovich , Monash University <i>“H3K27me3 mimicry restricts the spread of Polycomb domains”</i>	
11:40am – 12:00pm	Speaker 67: Isaac Hilton , Rice University <i>“Advanced epigenome editing technologies for probing and programming chromatin states and gene expression”</i>	
12:00pm – 12:20pm	Speaker 68: Matthew Easterday , Dovetail Genomics <i>“Structural variants drive enhancer hijacking and oncogene expression in an ovarian adenocarcinoma”</i>	
12:20pm – 12:40pm	Speaker 69: Kami Ahmad , Fred Hutchinson Cancer Center <i>“Histone gene control in cancer”</i>	
12:40pm – 1:00pm	Speaker 70: Nicolas Altemose , Stanford University <i>“Long-read, single-molecule sequencing methods for studying protein-DNA interactions”</i>	
1:00pm – 5:00pm	Free Time	

4:00pm – 5:00pm	Workshop 3: From Bench to Browser: Unlocking CUT&RUN and CUT&Tag Analysis with CUTANA™ Cloud Presenter: Keith Maier, PhD, <i>EpiCypher</i> Board of Directors Room	
5:00pm – 7:00pm	Session #13 Grand Coral Ballroom Chair: Yanzhong (Frankie) Yang , City of Hope Cancer Center	
5:00pm – 5:20pm	Speaker 71: Maria Elena Torres-Padilla , Helmholtz Munich <i>“Epigenetic mechanisms in early mammalian development”</i>	
5:20pm – 5:40pm	Speaker 72: Mo Chen , Tsinghua University <i>“Spatiotemporal control of SMARCA5 by a MAPK-RUNX1 axis distinguishes mutant KRAS-driven pancreatic malignancy from tissue regeneration”</i>	
5:40pm – 6:00pm	Speaker 73: Jacques Côté , Laval University Cancer Research Center <i>“Acetylation and incorporation of H2A.Z by NuA4/TIP60 in genome expression and stability”</i>	
6:00pm – 6:20pm	Speaker 74: Steve Josefowicz , Weill Cornell Medicine <i>“A histone code for stimulation induced chromatin activation and transcription”</i>	
6:20pm – 6:40pm	Speaker 75: Jake Kirkland , Oklahoma Medical Research Foundation <i>“A Repressive Role for PBRM1 in Balancing Pluripotency and Differentiation via Chromatin State Control”</i>	
6:40pm – 7:00pm	Speaker 76: Hiten Madhani , University of California, San Francisco <i>“Evolution of the 53BP1-RIF1-shieldin pathway”</i>	
7:30pm – 10:00pm	Farewell Reception (Dinner Included)	Sunset Terrace