

EpiCypher 2018

Biological & Clinical Frontiers in Epigenetics

November 6-10, 2018

Nassau, Bahamas

Keynote Speakers: **C. David Allis**
Rockefeller University

Kristian Helin
University of Copenhagen

This exciting four-day conference at the oceanfront luxury all-inclusive Meliá Nassau Beach Resort will highlight recent breakthroughs in epigenetics research. Many talks will be selected from submitted poster abstracts.



Confirmed Speakers:

Scott Armstrong
Dana-Farber Cancer Institute

Emily Bernstein
Icahn School of Medicine at Mount Sinai

Taiping Chen
U Texas MD Anderson Cancer Center

Xiaodong Cheng
U Texas MD Anderson Cancer Center

Katrin Chua
Stanford University School of Medicine

Wolfgang Fischle
King Abdullah University of Sci. & Tech.

Ben Garcia
University of Pennsylvania

Nada Jabado
McGill Univ. Montreal Children's Hospital

Steve Jacobsen
Univ. of California Los Angeles

Bill Janzen
Epizyme, Inc

Cigall Kadoch
Dana-Farber Cancer Institute

Bob Kingston
Massachusetts General Hospital

Ryan Kruger
GlaxoSmithKline

Ross Levine
Memorial Sloan Kettering Cancer Center

En Li
Novartis Inst. for Biomedical Research

Haitao Li
Tsinghua University

Karen Maegley
Pfizer, Inc

Tom Misteli
National Cancer Institute - NIH

Bing Ren
University of California San Diego

Stephane Richard
Lady Davis Inst. - McGill University

Alex Ruthenburg
Univ. of Chicago

Yang Shi
Harvard Medical School

Xiaobing Shi
U Texas MD Anderson Cancer Center

Rob Sims
Constellation Pharmaceuticals

Marc Timmers
University of Freiburg

Jerry Workman
Stowers Inst. for Medical Research

Bing Zhu
Chinese Academy of Sciences

Organizers:

Mark Bedford
University of Texas
MD Anderson Cancer Center

Or Gozani
Stanford University

Dan Levy
Ben Gurion University

Brian Strahl
University of North Carolina
Lineberger Cancer Center

Email us at epicypher2018@epicypher.com for more information.

Register at epicypher.com/epi2018

855-374-2461



EpiCypher™

Bringing Epigenetics to Life