

EpiCypher 2018

Biological & Clinical Frontiers in Epigenetics

November 6-10, 2018

Nassau, Bahamas

Keynote Speakers: C. David Allis

Kristian Helin

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
Emily Bernstein
Taiping Chen
Xiaodong Cheng
Katrin Chua
Nada Jabado
Steve Jacobsen

Bill Janzen
Cigall Kadoch
Bob Kingston
Ryan Kruger
En Li
Haitao Li
Hiten Madhani

Karen Maegley
Tom Misteli
Bing Ren
Stephane Richard
Alex Ruthenburg
Yang Shi
Xiaobing Shi

Rob Sims
Marc Timmers
Jonathan Whetstine
Jerry Workman
Bing Zhu

Organizers:

Mark Bedford
Or Gozani
Dan Levy
Brian Strahl

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

Register at epicypher.com/epi2018

855-374-2461



EpiCypher™

Bringing Epigenetics to Life