



FOR IMMEDIATE RELEASE

EpiCypher Announces EpiCypher2016 as Major Epigenetics Conference

PerkinElmer among sponsors of EpiCypher 2016.

Research Triangle Park, NC - 10 February 2016 – EpiCypher, Inc. has announced [EpiCypher2016](#), a conference focused on Biological and Chemical Frontiers in Epigenetics, to be held April 20-23 in San Juan, Puerto Rico. The conference is being organized by Drs. Brian Strahl, University of North Carolina at Chapel Hill; Or Gozani, Stanford University; and Mark Bedford, Univ. Texas MD Anderson Cancer Center.

The four-day conference at the luxury oceanfront Caribe Hilton in San Juan will highlight recent breakthroughs in epigenetics and chromatin research and will foster the collaboration and exchange of scientific expertise and knowledge in one of the most vibrant areas of life sciences. Through scientific presentations, a wide spectrum of topics in epigenetics and chromatin biology will be explored. The conference has over 25 confirmed speakers with a lineup of luminaries and thought leaders from industry and academic centers. PerkinElmer, Inc. is a sponsor for the event.

“EpiCypher2016 represents a significant commitment to the field and the thought leaders that are shaping the future of epigenetics. We’re delighted to have PerkinElmer as part of this inaugural event in light of our recent collaboration yielding a panel of bromodomain binding assays utilizing their highly sensitive AlphaLISA® detection technology.” said Sam Tetlow, CEO, EpiCypher. “We look forward to noteworthy announcements and scientific findings from EpiCypher2016.”

[EpiCypher2016](#) will feature keynote addresses by Dr. C. David Allis and Dr. Yang Shi. Dr. Allis is the Joy and Jack Fishman Professor and Head of the Laboratory of Chromatin Biology and Epigenetics at The Rockefeller University. Dr. Shi is the Merton Bernfield Professor of Neonatology at Boston Children’s Hospital / Harvard Medical School.

The plenary session talks will cover broad aspects of epigenetics and chromatin biology, ranging from basic molecular mechanisms to therapeutic efforts. There will be discussions around the elucidation of novel mechanisms, new tools, and other breaking topics in epigenetics and chromatin research.

About EpiCypher - A pioneer in the field of epigenetics and chromatin biology, EpiCypher is a bioscience company developing transformative technologies and delivering superior products to researchers worldwide. EpiCypher’s patented EpiTitan™ Histone Peptide Array is the most advanced histone array currently available. EpiCypher also sells nucleosomes, recombinant histone binding proteins, peptides and antibodies, as well as peptide synthesis and peptide array screening services.

For more information about EpiCypher, visit [EpiCypher.com](#).

Media Contact:

Sam C. Tetlow, CEO

EpiCypher, Inc.

919-627-7882 (direct)

sam@epicypher.com

###