



EpiCypher 2016

Biological & Clinical Frontiers in Epigenetics

Keynote Speaker: Yang Shi

April 20-23, 2016

San Juan, Puerto Rico

This exciting four-day conference at the luxury oceanfront Caribe Hilton will highlight recent breakthroughs in epigenetics research.

Confirmed Speakers

<i>Scott Armstrong</i>	<i>Ben Garcia</i>	<i>Alex Ruthenburg</i>
<i>Cheryl Arrowsmith</i>	<i>Steve Jacobsen</i>	<i>Julien Sage</i>
<i>Emily Bernstein</i>	<i>Ryan Kruger</i>	<i>Xiaobing Shi</i>
<i>Xiaodong Cheng</i>	<i>Bing Li</i>	<i>Jesse Smith</i>
<i>Katrin Chua</i>	<i>Minkui Luo</i>	<i>Paul Wade</i>
<i>Luciano Di Croce</i>	<i>Hiten Madhani</i>	<i>Jerry Workman</i>
<i>Wolfgang Fischle</i>	<i>Dinshaw Patel</i>	

Many talks will be selected from submitted poster abstracts.

Early Reg. Deadline: **March 1st, 2016**

Abstract deadline: **March 1st, 2016**

Please email abstracts to:
epicypher2016@epicypher.com

Space is limited - Register Soon!

Organizers

<i>Mark Bedford</i>	<i>Or Gozani</i>	<i>Brian Strahl</i>
---------------------	------------------	---------------------

Sponsored in part by 
PerkinElmer
For the Better

www.epicypher.com/epicypher-2016

855-374-2461

 **EPICYPHER**TM
Pioneering the Science of EpigeneticsTM