

## FOR IMMEDIATE RELEASE

### EpiCypher SNAP-ChIP® technology poised to transform market for histone PTM antibodies

**Research Triangle Park, NC – Oct. 4<sup>th</sup>, 2018** – A recently published report in *Molecular Cell* validates the use of EpiCypher’s proprietary SNAP-ChIP® technology for profiling antibody specificity and sensitivity. SNAP-ChIP® (**S**ample **N**ormalization and **A**ntibody **P**rofilng for Chromatin Immunoprecipitation) utilizes pools of DNA-barcoded designer modified nucleosomes (dNucs) as quantitative spike-in controls for ChIP-Seq normalization and antibody specificity testing.

Using this new approach, many commonly used ‘ChIP-grade’ antibodies failed due to significant cross-reactivity with other histone PTMs, including some of those recommended by the ENCODE consortium. These results highlight the need for internal controls to monitor antibody performance in each ChIP experiment, and position SNAP-ChIP as the new gold standard for antibody validation in ChIP.

*‘ The results are striking and will have an immediate impact on the field ’*, said Dr. Michael-Christopher Keogh, Chief Scientific Officer at EpiCypher and a corresponding author on the study. *‘ Our ongoing investigations demonstrate that ChIP-incapability, despite what their datasheets promise, is a widespread issue for histone PTM antibodies, and highlights a critical need for re-evaluating these reagents ’*.

EpiCypher is using SNAP-ChIP® technology to screen hundreds of commercially available antibodies to identify those with best-in-class performance. These “SNAP-ChIP® certified antibodies” are available now from EpiCypher and their use will be essential to provide a renewed confidence in data generated using ChIP.

**About EpiCypher** - A pioneer in the field of epigenetics and chromatin biology, EpiCypher is a biotechnology company developing transformative technologies and delivering superior products to researchers worldwide. EpiCypher manufactures and sells recombinant “designer” modified nucleosomes (dNucs), including the EpiDyne™ family of substrates for nucleosome remodeling assays, as well as recombinant histone binding proteins, peptides and antibodies, and offers a broad range of custom substrate manufacturing and assay development services.

For more information about SNAP-ChIP®, visit [EpiCypher.com/SNAP-ChIP](http://EpiCypher.com/SNAP-ChIP). For more information about SNAP-ChIP® certified antibodies visit [epicypher.com/snap-chip-certified-antibodies](http://epicypher.com/snap-chip-certified-antibodies)

#### Media Contact:

Ana Aza, PhD  
EpiCypher, Inc.  
[ana@epicypher.com](mailto:ana@epicypher.com)