



## FOR IMMEDIATE RELEASE

### EpiCypher Launches SNAP-ChIP™ Product Family: Now Enabling Normalized Chromatin IP Experiments.

**Research Triangle Park, NC - 22 September 2017** – EpiCypher, Inc. has released **SNAP-ChIP** (Sample Normalization and Antibody Profiling for Chromatin Immunoprecipitation), a breakthrough product family that will dramatically improve researchers' ability to normalize ChIP experiments. SNAP-ChIP utilizes EpiCypher's proprietary DNA barcoded recombinant designer nucleosomes (dNucs) as next-generation spike-ins for ChIP. With SNAP-ChIP, researchers now have the ability to both normalize samples across experiments (enabling broad collaboration throughout the scientific community, and generate detailed data pertaining to antibody specificity and cross-reactivity).

The first offering in the SNAP-ChIP family, the SNAP-ChIP K-MetStat panel (lysine-methylation status panel), is comprised of 16 uniquely modified dNucs carrying disease-relevant lysine methylation modifications on histones H3 and H4. Each modification site in this pool (H3K4, H3K9, H3K27, H3K36, H4K20) is represented by all lysine methylation states (*i.e.* me0, me1, me2, and me3). "For the first time, scientists can both monitor technical variability and measure antibody specificity in the same experiment. This has not been possible with any other spike-in product or approach currently available," states EpiCypher's Chief Scientific Officer, Michael Keogh, PhD. Keogh further adds "The beauty of the SNAP-ChIP product line is that it fits seamlessly into existing ChIP workflows, allowing the epigenetics community immediate access to this breakthrough. In fact, we have already contracted with a number of antibody development companies to use our SNAP-ChIP K-MetStat Panel to directly analyze the specificity of antibodies in their portfolios in a ChIP setting."

Numerous additional modification-specific SNAP-ChIP dNuc panels are currently in development (*e.g.* arginine methylation, lysine acylation, etc.) and are slated for release in early 2018.

**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 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 EpiCypher, visit [EpiCypher.com](http://EpiCypher.com).

#### Media Contact:

Jim Bone, PhD

President, EpiCypher, Inc.

434-996-8107

[jim@epicypher.com](mailto:jim@epicypher.com)