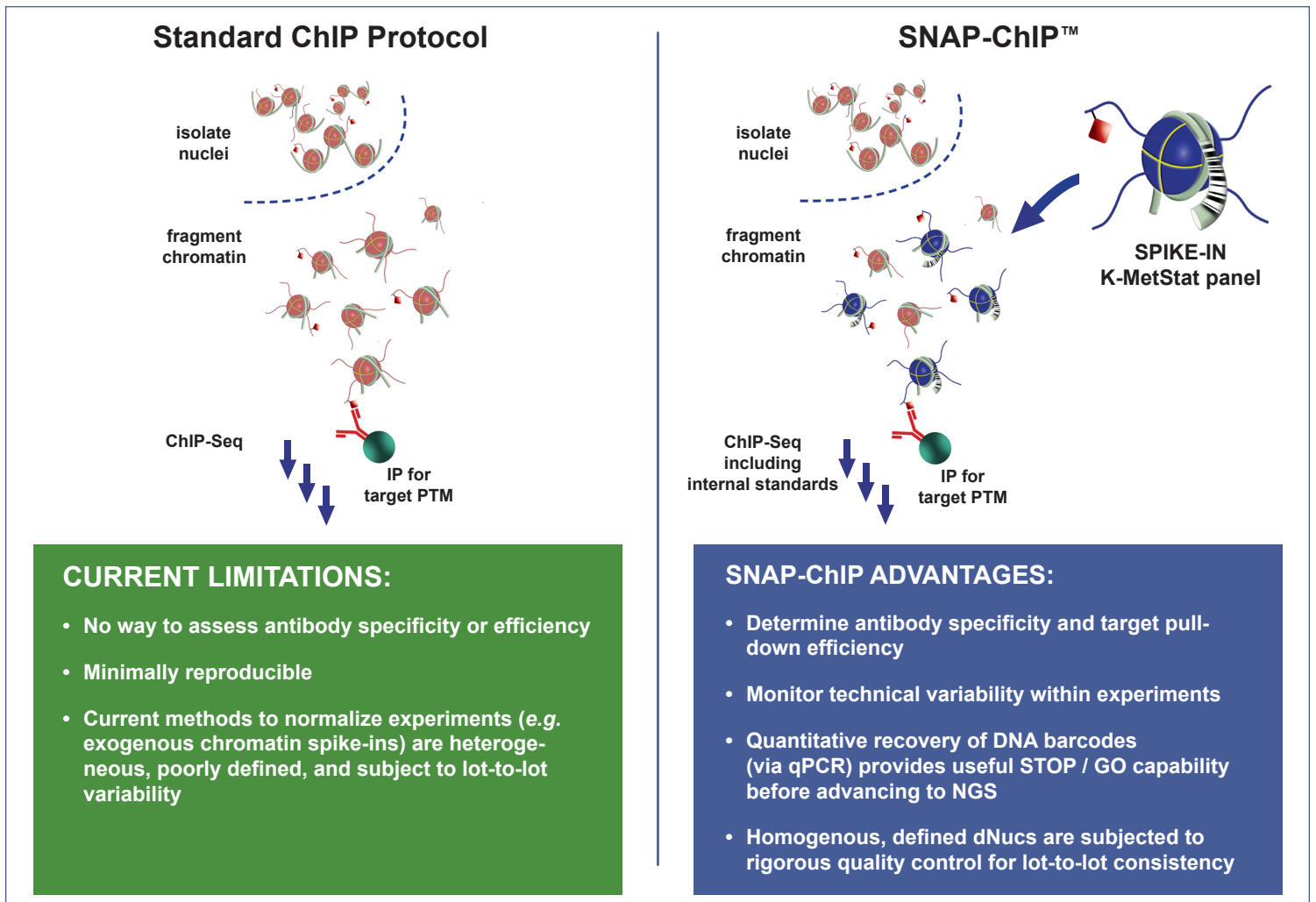


# SNAP-ChIP™ for histone lysine methylation

Sample Normalization & Antibody Profilng for ChIP using EpiCypher's **K-MetStat™** panel

**SNAP-ChIP** uses DNA barcoded recombinant designer nucleosomes (dNucs) as next-generation spike-ins for chromatin immunoprecipitation (ChIP). The first product in this family consists of a panel of dNucs carrying well-studied histone lysine methyl marks (**K-MetStat panel**). EpiCypher's K-MetStat panel can easily be added to any ChIP workflow to standardize and normalize samples (across experiments) and rigorously validate antibody specificity. For the first time, users can monitor antibody specificity and evaluate technical variability within a ChIP experiment, setting SNAP-ChIP apart from any other spike-in controls currently available on the market.

**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 the panel (H3K4, H3K9, H3K27, H3K36, H4K20) is represented by all lysine methylation states (*i.e.* me0, me1, me2, and me3). This allows for maximum user flexibility and provides the ability to gather detailed antibody cross-reactivity data. Additional modification-specific dNuc panels are currently in development (*e.g.* arginine methylation, lysine acetylation, etc.) and slated for release in 2018.

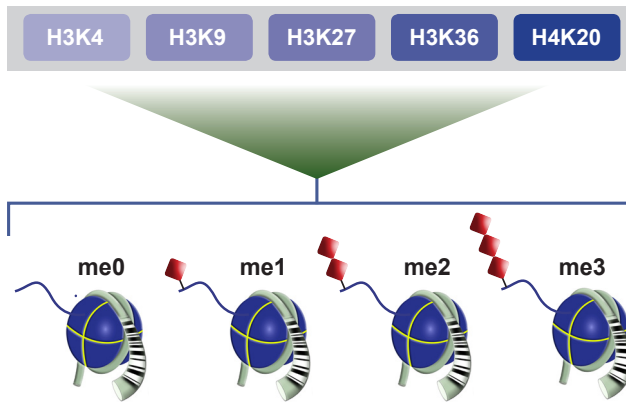


**Figure 1:** Overview of the SNAP-ChIP approach (adapted from ICe-ChIP technology: Grzybowski *et al. Mol Cell*, Vol. 58, Issue 5, 886 - 899, 2015). A pool of recombinant dNucs with defined post-translational modifications (PTMs) identified by unique DNA barcodes is added to sample chromatin prior to immunoprecipitation (IP). Capture of the barcoded nucleosomes (on / off target) allows the user to assess antibody specificity, monitor technical variability, and normalize experiments. Quantitative recovery of barcoded dNucs (via qPCR) provides a useful STOP / GO capability prior to advancing to next-generation sequencing.

**K-MetStat Panel details** ▶

## K-MetStat panel:

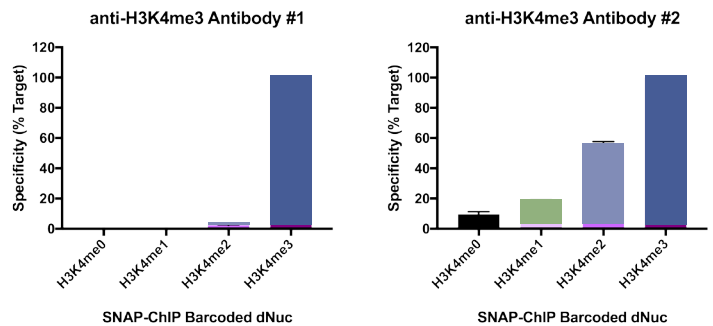
16 uniquely modified designer nucleosomes



**Figure 2:** Schematic depicting the 16 dNucs included in the K-MetStat Panel, each uniquely DNA barcoded.

## Why do I need to assess antibody specificity in ChIP?

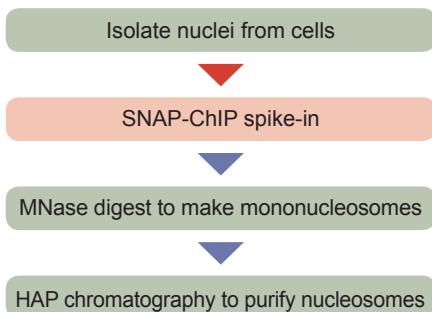
- Antibody cross-reactivity can lead to gross misinterpretation of biological findings
- Methylation states (me0/1/2/3) are challenging targets for antibodies due to high degree of structural similarities
- SNAP-ChIP addresses these limitations by enabling quantification of antibody specificity within every ChIP experiment:



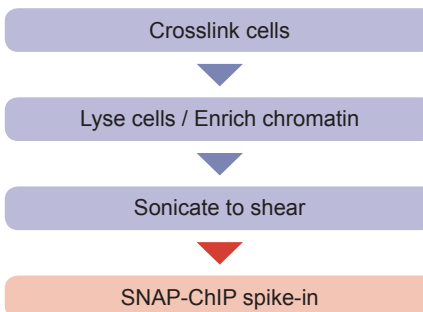
**Figure 3:** A SNAP-ChIP experiment (n = 3) using two H3K4me3 antibodies shows that Antibody #1 exhibits <3% cross-reactivity with alternate H3K4 methyl states; Antibody #2 shows ~60% cross-reactivity with H3K4me2, substantially compromising ChIP-studies.

## SNAP-ChIP seamlessly integrates into existing ChIP workflows

### Native ChIP Workflow:



### Crosslinked ChIP Workflow:



Immunoprecipitate nucleosomes using antibody against target histone PTM

Purify DNA

qPCR to determine **antibody specificity, & technical variability**

**STOP / GO decision**

Next Generation Sequencing (NGS) to identify epigenetic changes of interest

**Normalize data by equalizing SNAP-ChIP spike-ins across samples**

## ORDERING INFO:

### SNAP-ChIP K-MetStat

Catalog No. 19-1001  
Price: \$249 / 10 ChIP equivalents  
Website: [EpiCypher.com/SNAP-ChIP](http://EpiCypher.com/SNAP-ChIP)

### Related Products (Coming Soon):

Lysine acetylation status panel (K-AcStat)  
Arginine methylation status panel (R-MetStat)

### Related Nucleosome products:

EpiDyne™ Nucleosome Remodeling Substrates  
Recombinant Nucleosomes (rNucs)  
Designer Nucleosomes (dNucs)  
Variant Nucleosomes (vNucs)  
Oncogenic Nucleosomes (OncoNucs)  
Purified Nucleosomes (HeLa, Chicken)

