



PDF poster download

Biochemical and genomic approaches for high throughput drug discovery in chromatin remodeling research



Lu Sun¹, Tessa Firestone¹, Hannah Willis¹, Matthew R. Marunde¹, Vishnu U. Sunita Kumary¹, Matthew J. Meiners¹, Saarang Gopinath¹, Jonathan M. Burg¹, Bryan J. Venters¹, Allison Hickman¹, Zu-Wen Sun¹, Martis W. Cowles¹, Pierre Esteve², Hang Gyeong Chin², Chaithanya Ponnaluri², Sriharsa Pradhan² & Michael-Christopher Keogh¹

¹ EpiCypher Inc, Durham, NC 27709
² New England Biolabs, Ipswich, MA 01938

EpiDyne[®] remodeling assay and genomic approaches in remodeler research

Chromatin remodeling is mediated by ATP-dependent enzymes that play key roles regulating gene expression and genome replication / repair. Aberrant nucleosome organization from dysregulated chromatin remodeling can severely alter chromatin accessibility and disrupt these important processes, thereby driving various cancers. Remarkably, nearly 20% of all human cancers contain mutations in subunits from the SWI/SNF family of chromatin remodeling complexes, making them of great interest to basic research and therapeutic intervention^{1,2}.

In vitro studies on the remodeling enzymes such as SMARCA2/4/5 (and their multi-subunit complexes) are challenging, partially due to the strong preference for nucleosome-based substrates (the physiological target of these enzymes). We have created the EpiDyne[®] nucleosome portfolio to examine chromatin remodeler activity in biochemical assays, and here present the development of novel readouts (-PicoGreen[™] and -TR-FRET). These nonradioactive plate-based assays are automation adaptable, ready for high-throughput inhibitor screening, and can be customized for various remodeling enzymes that exhibit preferences in nucleosome composition (e.g. histone type or DNA linker length).

For parallel *in vivo* studies we note that genome-wide remodeler localization and open chromatin mapping are fundamental for understanding the function / activity of these enzymes in cancer development and inhibitor responses. However, traditional genomic approaches have significant issues: e.g. ChIP-seq demands high cell numbers and sequencing depths, and is unable to effectively map ATPases without heavily modified protocols; while ATAC-seq to map open regions deals poorly with cross-linking that could stabilize transient states of interest. To these ends, we optimized the CUTANA[™] CUT&RUN approach to efficiently capture the localization of all major classes of chromatin remodelers with high signal to background in multiple tissue cultures, enabling evaluation of specificity and/or efficiency from remodeler targeting inhibitors or degraders. We have also adopted NicE-seq for chromatin accessibility profiling in cross-linked material that captures landscape changes in response to remodeling disruptions. As complementary tools to the EpiDyne platform, CUT&RUN and NicE-seq facilitate epigenomic research on chromatin remodelers in cancer therapeutic intervention.

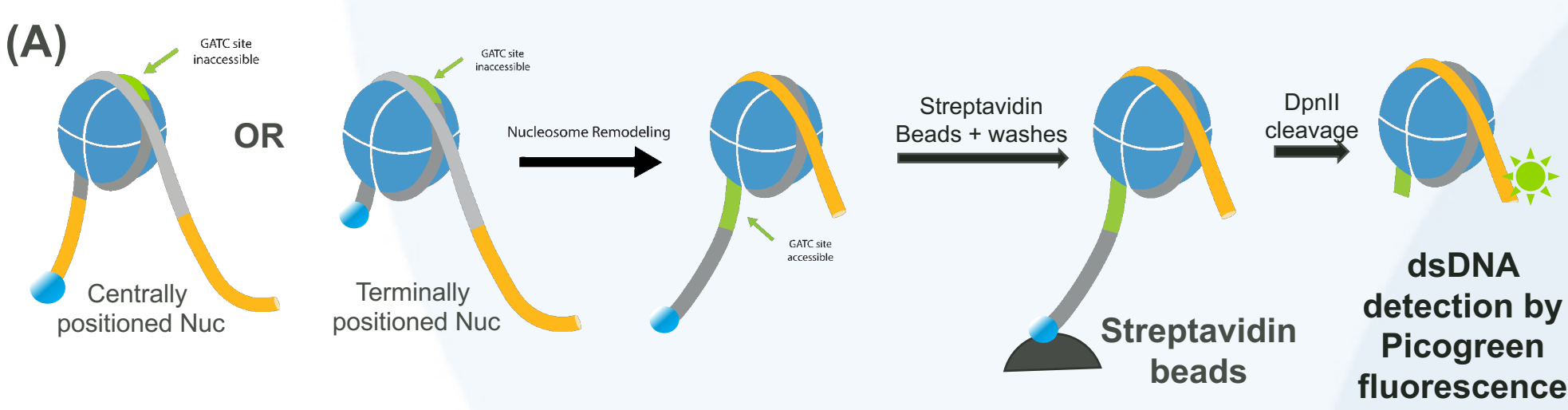


Figure 1. EpiDyne Nucleosome remodeling Assay Designs. (A) PicoGreen³ (B) TR-FRET

EpiDyne-picogreen (HTS compatible) reveals remodeler substrate preference

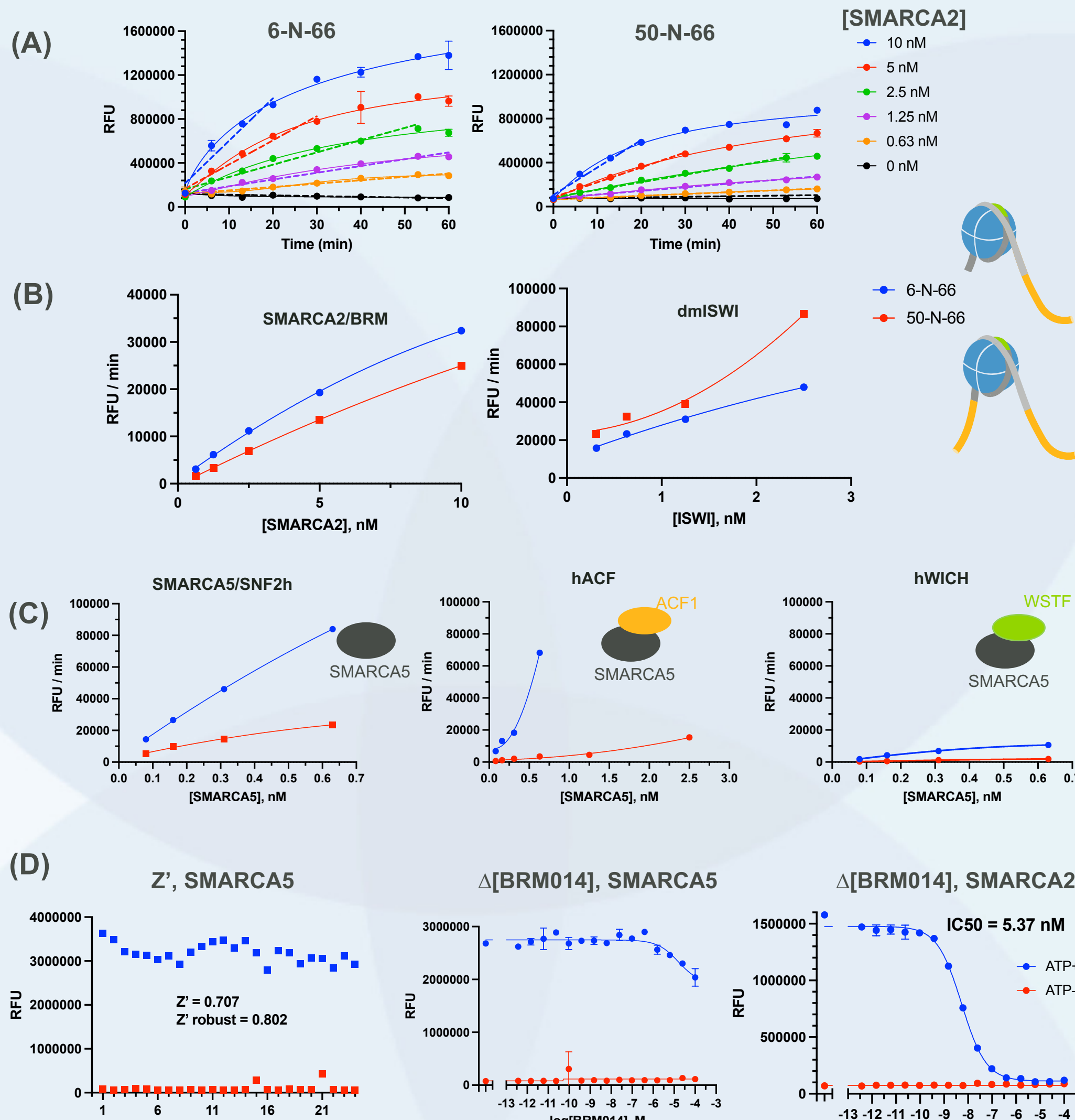


Figure 2. EpiDyne-PicoGreen remodeling assay. (A) Enzyme-, ATP- and time-dependent remodeling reactions by SMARCA2 (BRM). (B) Initial rates for terminally [6-N-66] or centrally [50-N-66] positioned nucleosomes by various ATPases⁴, plotted for reactions within linear ranges. (C) Substrate preference for various SMARCA5 remodeling complexes in EpiDyne-picogreen. (D) Z' analysis and tool compound^{5,6} dose response for SMARCA5/SNF2H compared to SMARCA2/BRM remodeling with EpiDyne-PicoGreen.

EpiDyne-TR-FRET (HTS compatible) for real time remodeler studies

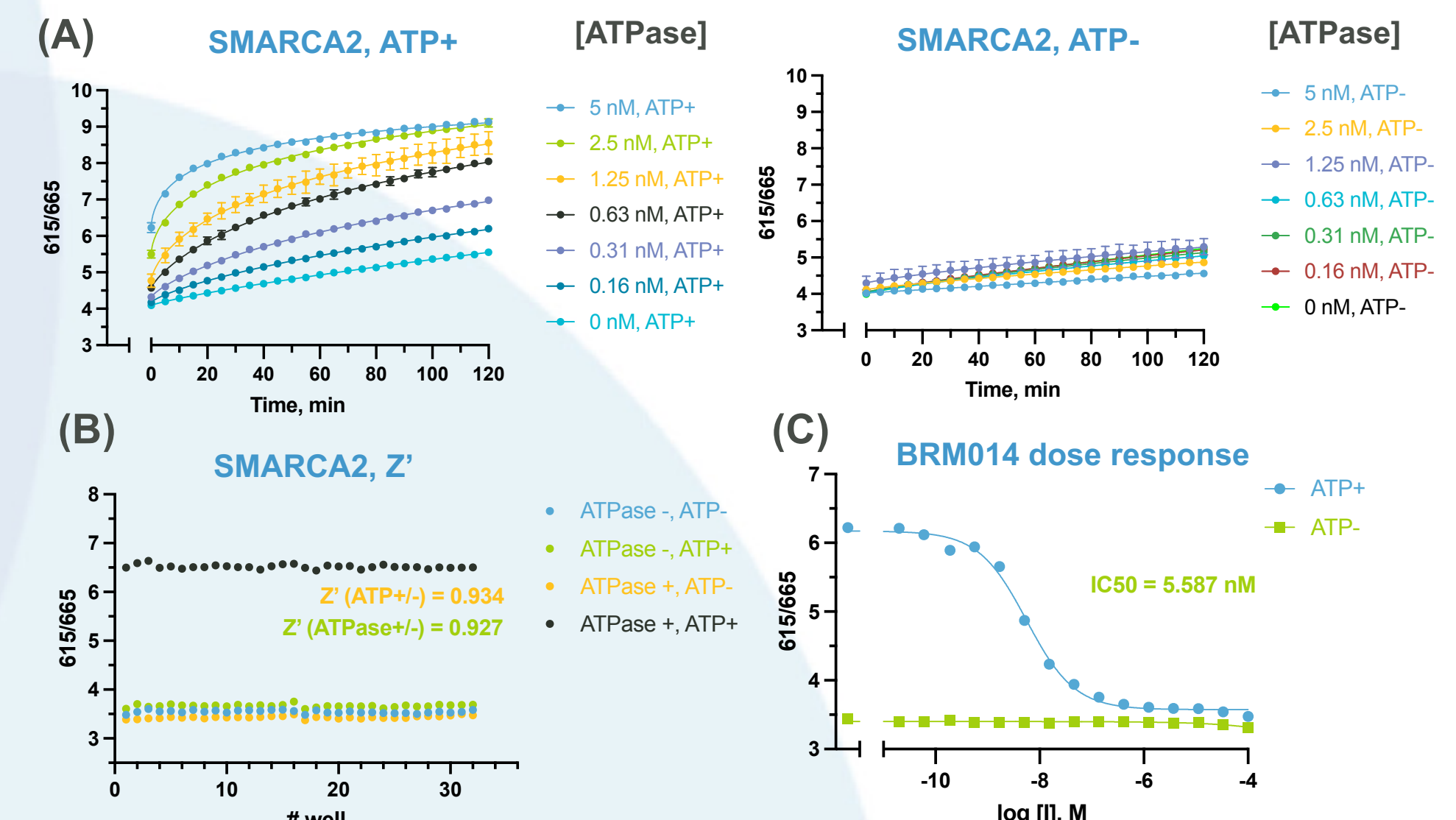


Figure 3. EpiDyne-TR-FRET remodeling assay. (A) Enzyme-, ATP- and time-dependent remodeling reactions by SMARCA2. (B) Z' analysis and (C) tool compound^{5,6} inhibition of SMARCA2/BRM remodeling in EpiDyne-TR-FRET.

| | PicoGreen | TR-FRET |
|---------------------------|----------------------|-----------------------|
| Suitable Nucleosome | terminal/central | terminal |
| Assay readout | end point | real time / end point |
| Operation time | 4-5 hr | ~3 hr |
| | [2 hr hands on] | [0.5 hr hands on] |
| S/B window | 5 ~ 20 | 1.5 ~ 2.5 |
| Z' with SMARCA2/4/5 | > 0.7 | > 0.8 |
| Fluorescence interference | Very low | low |
| Assay format | 384-well plate based | 384-well plate based |

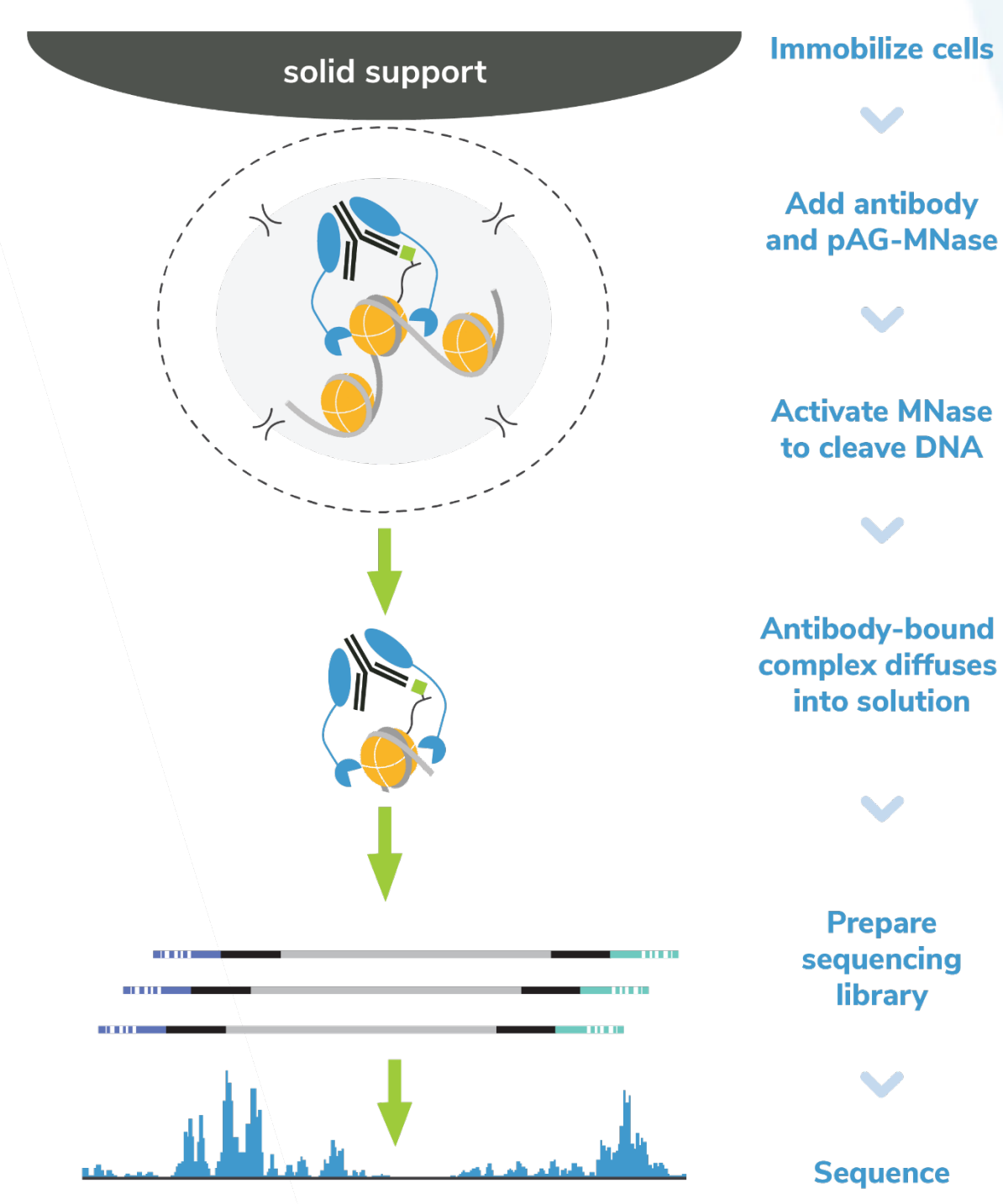
Figure 4. Performance comparisons of EpiDyne remodeling assays.

Acknowledgement

We are grateful to Dr. Dmitry Fyodorov (*Albert Einstein College of Medicine*) for the kind gift of dm ISWI⁴ and human SMARCA5 complexes. This work was supported by multiple awards from the National Institutes of Health (R44GM116584, R44GM123869, R44DE029633, R44HG010640, and R44HG011006).

Functional epigenomic approaches in remodeler research

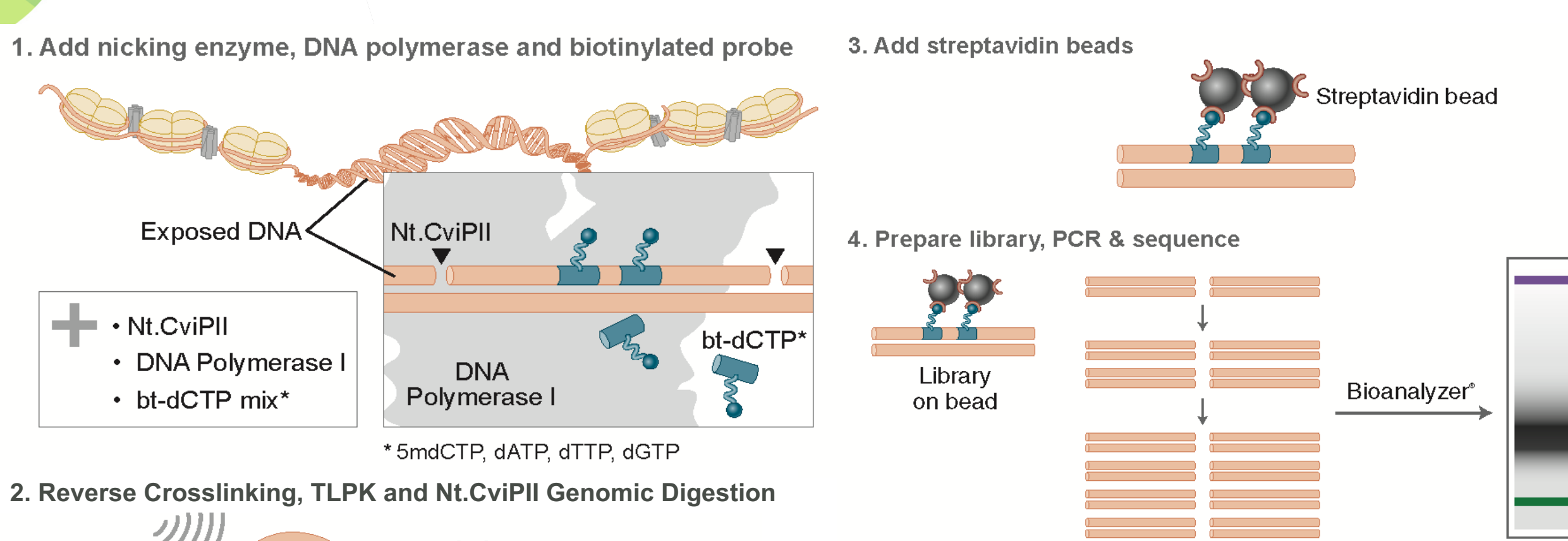
CUT&RUN Workflow



CUT&RUN vs. ChIP-seq

| Platform Comparison | ChIP-seq | CUTANA [™] CUT&RUN |
|--------------------------|-------------------|--|
| Required Cells | >1 million | 5,000-500,000 |
| Ideal for Profiling | Histone PTMs, TFs | Histone PTMs, TFs & chromatin remodelers |
| Sequencing Depth (Reads) | >30 million | 3-8 million |
| Experimental Throughput | Low | High |
| Signal-to-Noise | Low | High |
| Assay Automation | Difficult | Yes |

One-pot Uni-NicE-seq vs. ATAC-seq



| Platform Comparison | ATAC-seq | NicE-seq |
|--------------------------|---|--|
| Required Cells | Single-50k nuclei | 25-50k nuclei |
| Sequencing Depth (Reads) | >50 million | 30-35 million |
| Strengths & limitations | <ul style="list-style-type: none"> Field standard Sensitive | <ul style="list-style-type: none"> Robust and sensitive Low mitochondrial contamination (<5%) Works with fixed and FFPE tissue |

Distinct remodeler mapping responses to SMARCA2/4 inhibitors in H1299 cells

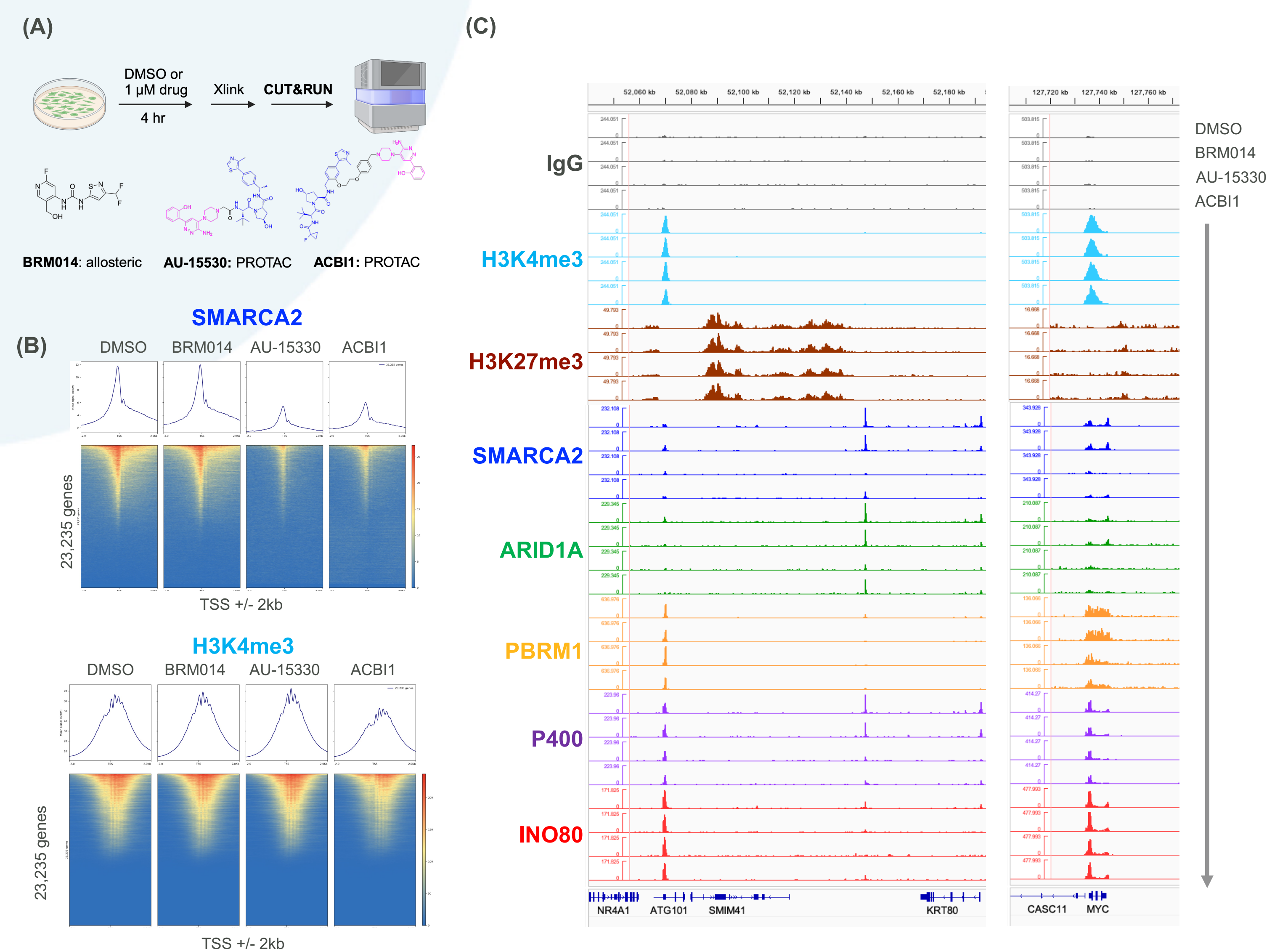


Figure 4. CUT&RUN validation of CUTANA compatible antibodies against remodelers in drug treatments. (A) Experimental workflow. Proliferating H1299 cells were treated with 1 μM of SMARCA2/4 targeting BRM014⁶, AU-15330⁹, ACB1¹⁰, or 0.1% DMSO vehicle control for 4 hours before light crosslinked and harvested for CUT&RUN mapping. (B) Heatmaps showing SMARCA2 and H3K4me3 enrichment at TSS and their responses to variant drug treatments. Rows are aligned, ranked and group scaled by intensity (top to bottom) in respective DMSO controls and colored such that red indicates high localized enrichment and blue denotes background signal. (C) Representative chromatin regions showing enrichment of core or accessory remodeler proteins in response to DMSO, various allosteric or PROTAC based drugs, group scaled by individual targets.

SMARCA2/4 disruption alters chromatin accessibility

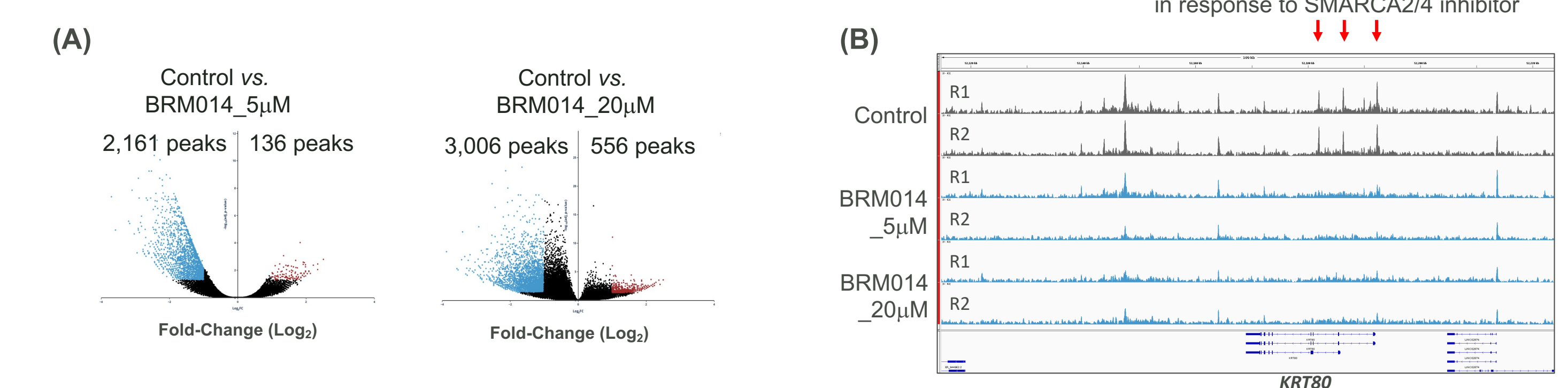


Figure 5. NicE-seq⁹ measures chromatin response to SMARCA2/4 remodeling inhibitor. Peaks were called from 100k formaldehyde-fixed NCI-H1299 cells (untreated / DMSO / + BRM014⁶). (A) Volcano plots of differential peak occupancy between untreated and H1299 drug treated cells. (B) Representative biomarker (*KRT80*); locus functionality is reliant on SMARCA2 (BRM) function⁹ showing open chromatin changes in response to BRM014 treatment.

References

- Hoffman *et al* (2014) Functional epigenetics approach identifies BRM/SMARCA2 as a critical synthetic lethal target in BRG1-deficient cancers. *PNAS* 111:3128.
- Hodges *et al* (2016) The Many Roles of BAF (mSWI/SNF) and PBAF Complexes in Cancer. *Cold Spring Harbor Perspect Med* 6:a026930.
- Andreyeva *et al* (2022) *Drosophila* SUMM4 complex couples insulator function and DNA replication control. *Elife* 11:e81828.
- Fyodorov DV & Kadonaga JT (2003) Chromatin assembly *in vitro* with purified recombinant ACF and NAP-1. *Methods Enzymol* 371:499.
- Papillon *et al* (2018) Discovery of Orally Active Inhibitors of Brahma Homolog (BRM)/SMARCA2 ATPase Activity for the Treatment of Brahma Related Gene 1 (BRG1)/SMARCA4-Mutant Cancers. *J Med Chem* 61:10155.
- Jagani *et al* (2019) In-Depth Characterization and Validation in BRG1-Mutant Lung Cancers Define Novel Catalytic Inhibitors of SWI/SNF Chromatin Remodeling. *bioRxiv* (10.1101/812628v1)
- Chin *et al* (2020) Universal NicE-seq for high-resolution accessible chromatin profiling for formaldehyde-fixed and FFPE tissues. *Clin Epigenet* 12: 143.
- Vishnu *et al* (2021) One-pot universal NicE-seq: all enzymatic downstream processing of 4% formaldehyde crosslinked cells for chromatin accessibility genomics. *Epigenetics & Chromatin* 14:53.
- Xiao *et al*. (2021) Targeting SWI/SNF ATPases in enhanced-addicted prostate cancer. *Nature*. 601, 434-439.
- Farnaby *et al* (2019) BAF complex vulnerabilities in cancer demonstrated via structure-based PROTAC design. *Nat. Chem Biol*, 15, 672-680.