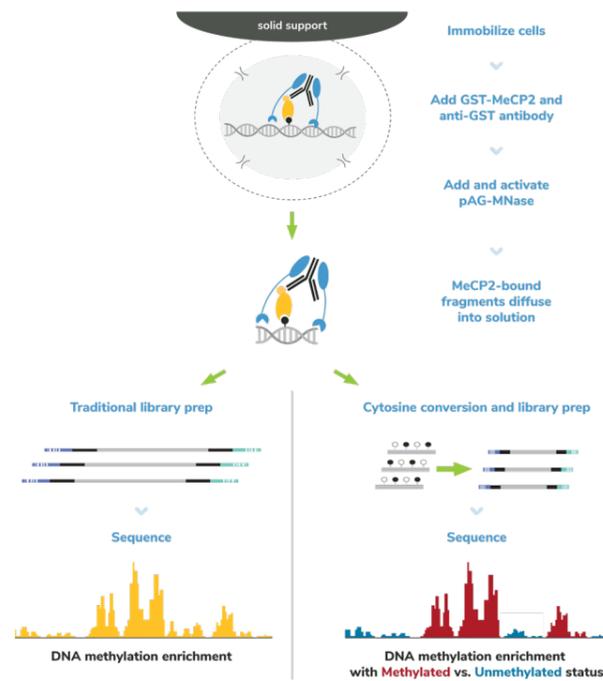


CUTANA™ meCUT&RUN

Genome-Wide
DNA Methylation Sequencing



Part 1: Enrich for methylated DNA

- A GST-tagged methylation-binding domain derived from human MeCP2 is used to capture methylated DNA fragments.

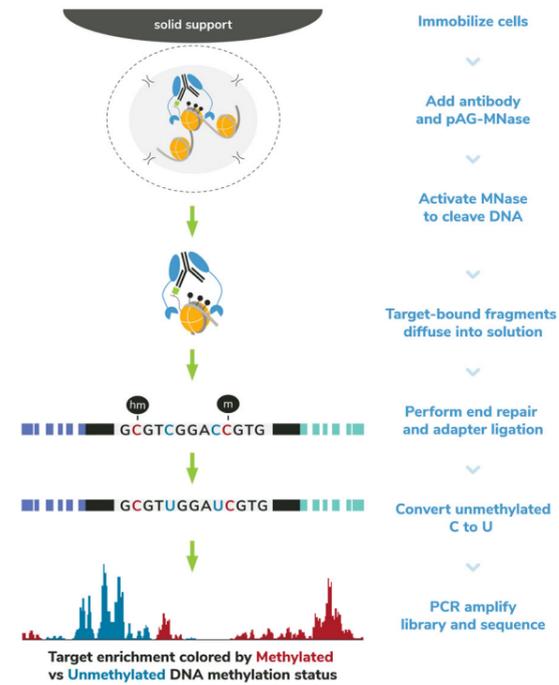
Part 2: Library prep and sequencing

Users have two options, providing maximal flexibility:

- Option 1 (meCUT&RUN):** Traditional library prep provides an overview of 5mC enrichment. Requires **15-20 M** total sequencing reads.
- Option 2 (meCUT&RUN-EM):** Cytosine conversion and library prep provides base-pair resolution 5mC analysis. Requires **30-50 M** total reads.

CUTANA™ Multiomic CUT&RUN

Simultaneous analysis of DNA
methylation and chromatin proteins



Part 1: CUT&RUN to enrich target chromatin

- CUT&RUN technology is used to first enrich for antibody-bound chromatin.
- Robust for histone PTMs, transcription factors, and other chromatin protein targets.

Part 2: Cytosine conversion & library prep

- Methylated CpGs are detected at target proteins using either bisulfite conversion or Enzymatic Methyl-seq (NEB® EM-seq™).
- EM-seq is the preferred method for Multiomic CUT&RUN because it is less harsh than bisulfite, preserves DNA integrity, and is optimized for low-input samples — making it ideal for CUT&RUN applications.



CUTANA™ DNA Methylation Sequencing Assays

meCUT&RUN



Low-cost, Genome-Wide
DNA Methylation Mapping

Multiomic CUT&RUN



Simultaneously map
DNA methylation and
chromatin proteins



epicypher.com | info@epicypher.com
855.374.2461



CUTANA™ meCUT&RUN

An Affordable Solution for Genome-Wide DNA Methylation Sequencing

CUTANA™ meCUT&RUN is a novel enrichment-based method for DNA methylation (DNAm) sequencing that delivers **high-quality, genome-wide profiles** at a **fraction of the cost and input required** for whole-genome bisulfite sequencing (WGBS).

WGBS require deep sequencing, while targeted methods sacrifice coverage and resolution. DNAm researchers thus face a tough choice between expensive whole-genome bisulfite sequencing (WGBS) or limited insights from reduced representation bisulfite sequencing (RRBS) and other assays.

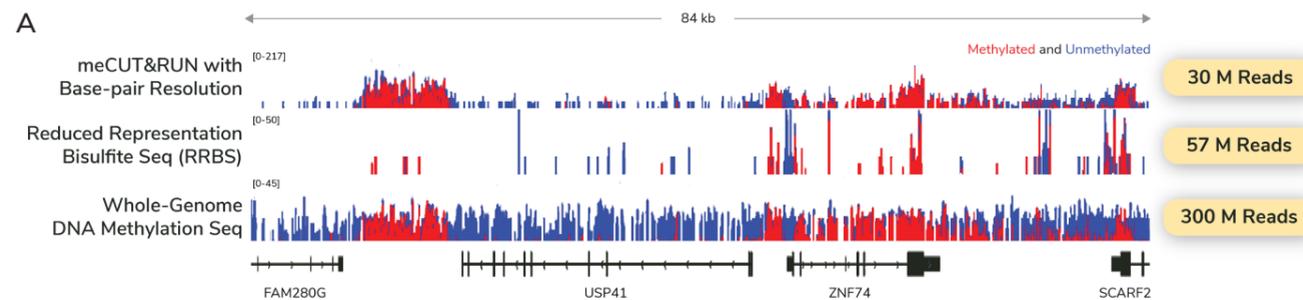
meCUT&RUN is a low-cost DNAm mapping solution that doesn't sacrifice coverage

| Technology | meCUT&RUN | RRBS | Microarrays | Hybridization Panels | WGBS |
|------------------------------|---------------------|----------------------|--------------------|----------------------|------------------|
| Input | 10k - 500k cells | 10 - 500 ng DNA | 250 ng DNA | 200 ng DNA | 50 ng - 5 µg DNA |
| Coverage | 80% Methylated CpGs | Limited (5-15% CpGs) | Limited (~3% CpGs) | Limited (~15% CpGs) | Whole genome |
| Recommended Sequencing Depth | >15 M reads | 50-100 M reads | N/A | >100 M reads | >800 M reads |
| Turnaround Time | Days | Weeks | Days | Weeks | Weeks |
| Base-Pair Conversion | Optional | Required | Required | Required | Required |
| Overall Cost | \$ | \$\$ | \$\$ | \$\$ | \$\$\$ |

Why Choose meCUT&RUN:

- High coverage**
Captures 80% of methylated CpGs.
- Low cost**
20-fold reduced sequencing costs vs. WGBS.
- Compatible with precious samples**
Robust down to 10,000 cells.
- Minimal bias**
Avoids destructive bisulfite conversion.

meCUT&RUN delivers genome-wide signal with low sequencing requirements



meCUT&RUN provides comprehensive coverage across functional elements, far exceeding RRBS

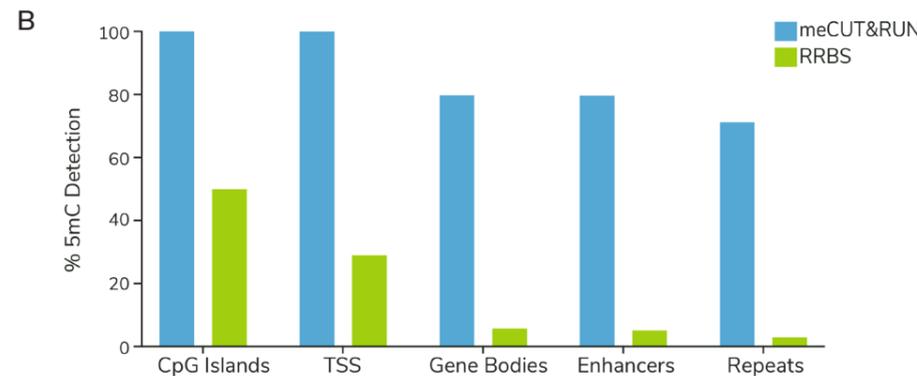


FIGURE 1
CUTANA™ meCUT&RUN was paired with enzymatic conversion to thoroughly evaluate its performance against targeted bisulfite-conversion assays (RRBS) and whole-genome, enzymatic-based approaches (EM-seq; 300 M reads) in K562 cells.
(A) Representative sequencing tracks.
(B) Percent 5mC coverage calculated relative to EM-seq.

CUTANA™ Multiomic CUT&RUN

Next-Gen Multiomic Profiling of DNA Methylation and Chromatin Proteins

CUTANA™ Multiomic CUT&RUN is a next-generation epigenomic mapping strategy that **profiles chromatin proteins and DNA methylation together** in a **single reaction**, capturing novel insights to drive biological discovery.

Traditional multiomic DNA methylation approaches integrate ChIP-seq with whole genome bisulfite sequencing (WGBS) datasets. However, these assays are cost-prohibitive for many labs, require large numbers of cells, take weeks (or longer!) to complete, and generate data that are indirect, restricting biological interpretations.

Multiomic CUT&RUN provides insights of chromatin regulation in a single workflow

| Technology | Multiomic CUT&RUN | ChIP-seq + Whole Genome Bisulfite Sequencing | |
|------------------------------|-------------------|---|------------------|
| | 1 Simple Workflow | Workflow 1: ChIP-seq | Workflow 2: WGBS |
| Input | 5k-500k cells | >1 million cells | 50 ng - 5 µg DNA |
| Turnaround Time | 4-5 Days | Weeks | Weeks |
| Multiomic Data? | Yes | No | No |
| Recommended Sequencing Depth | >30 M reads | >30 M reads | >800 M reads |
| Computational Complexity | Low | Requires complex integration of two distinct datasets | |
| Overall Cost | \$ | \$\$\$ | |

Why Choose Multiomic CUT&RUN:

- Reveal the full regulatory landscape**
Map DNA methylation at protein binding sites with base-pair resolution.
- Streamline discovery with a single workflow**
Multiple -omic insights from one sample.
- Low cost, maximal sensitivity**
Low cell inputs and sequencing requirements provide flexibility.

Multiomic CUT&RUN maps DNA methylation at diverse chromatin targets

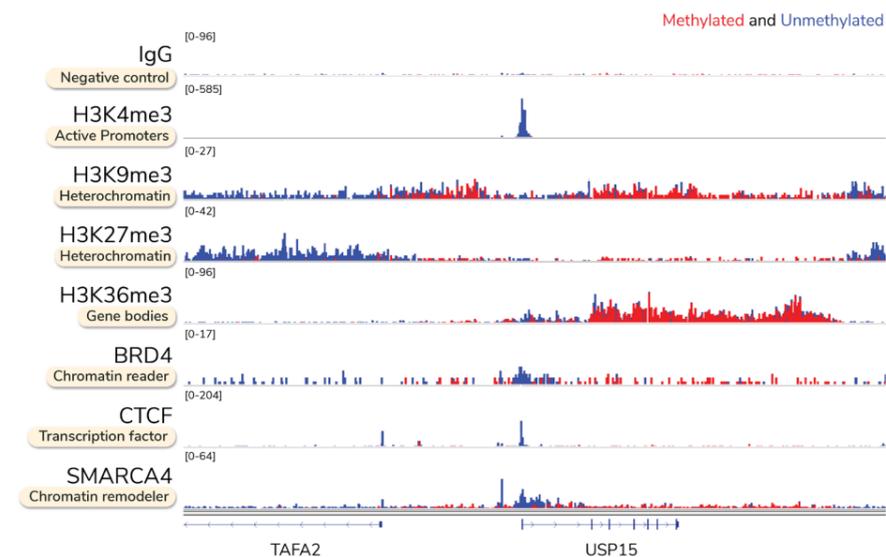


FIGURE 2
CUTANA™ Multiomic CUT&RUN provides base-pair resolution of CpG methylation at transcription factor binding sites, histone PTMs, and chromatin-modifying enzymes.
Each reaction used 500,000 K562 cells. Target chromatin was enriched with the CUTANA Multiomic CUT&RUN Workflow, followed by Enzymatic Methyl-seq (NEB® EM-seq™) conversion and library prep.
Key Applications:

- Gauge the impact of DNA methylation on transcription factor binding
- Resolve crosstalk between histone modifications and DNA methylation
- Understand the epigenetic mechanisms of gene regulation