

# CUT&RUN for advanced MeRIP applications

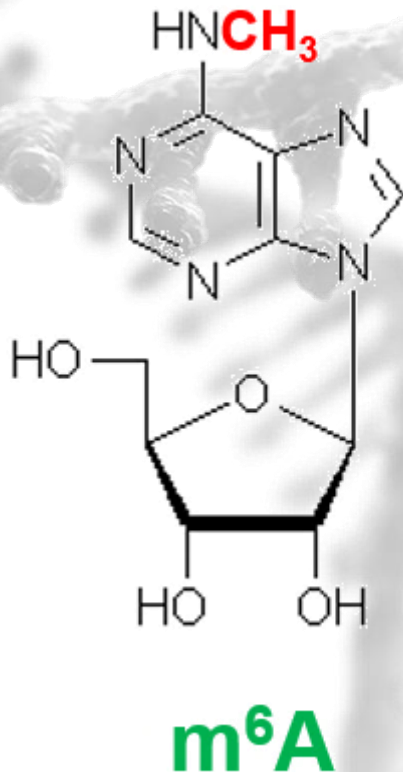


EpigenTek Group Inc.  
July 2023

# Product Summary

	<u>CUT&amp;RUN MeRIP Kit</u>	<u>CUT&amp;RUN MeRIP-Seq Kit</u>
<b>Catalog No.</b>	P-9018	P-9016
<b>Purpose</b>	Enrich short RNA fragments containing m <sup>6</sup> A methylation sites for downstream analysis (PCR, NGS)	m <sup>6</sup> A RNA enrichment + library preparation for NGS
<b>Kit Size</b>	24 rxns	12 and 24 rxns
<b>Starting Material</b>	Isolated RNA	
<b>Input Amount</b>	1 - 20 µg per rxn 10 µg optimal 500 ng minimum	
<b>Protocol Time</b>	< 3 h	< 6 h
<b>Antibody Notes</b>	Highly-specific and MeRIP-grade anti-m6A antibody is included with kit and not cross-reactive to unmethylated A	

# m<sup>6</sup>A RNA Modification



## N<sup>6</sup>-Methyladenosine

- “The fifth RNA base”
- Most common and abundant eukaryotic RNA modification
- Accounts for >80% of all RNA methylation
- Found mainly in mRNA; also observed in non-coding species (tRNA, rRNA, miRNA, etc.)
- Functions
  - RNA metabolism
  - Cell differentiation
  - Immunity
  - Inflammation
  - Circadian clock
  - Viral life cycle

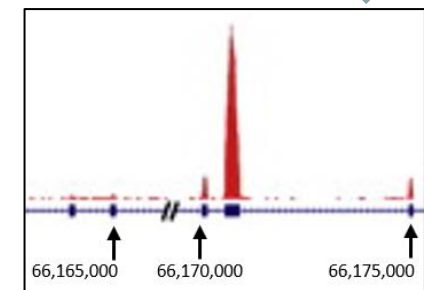
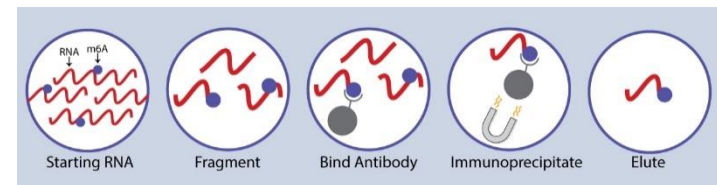
# MeRIP-seq

Methylated RNA ImmunoPrecipitation Sequencing

- First method to detect m<sup>6</sup>A on transcriptome-wide level
- Couples m<sup>6</sup>A RNA immunoprecipitation with NGS
- Allows for high-throughput localization of modified sites from enriched m<sup>6</sup>A-containing RNA fragments

## **Procedure**

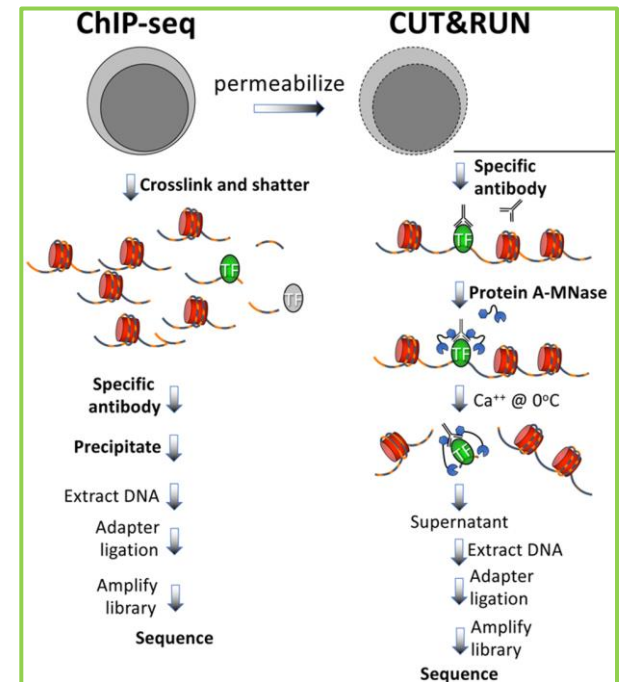
- RNA extraction
- RNA fragmentation
- Binding of magnetic bead-coupled antibody to sheared RNA
- Immunoprecipitation of m<sup>6</sup>A RNA fragments
- RNA separation from bound antibody and elution
- Reverse transcription to generate cDNA
- Library construction
- Sequencing



# CUT&RUN

Cleavage Under Targets and Release Using Nuclease

- Recent innovation in protein/DNA interaction studies
- Originally developed to resolve the shortcomings of ChIP
  - Large amount of starting material
  - Long protocol time (several days)
  - Limited resolution



Kaya-Okur et al., 2020

- **EpigenTek** has developed a **new MeRIP method** using CUT&RUN technology for improved m6A enrichment

# CUT&RUN m<sup>6</sup>A MeRIP

- **Advantages**

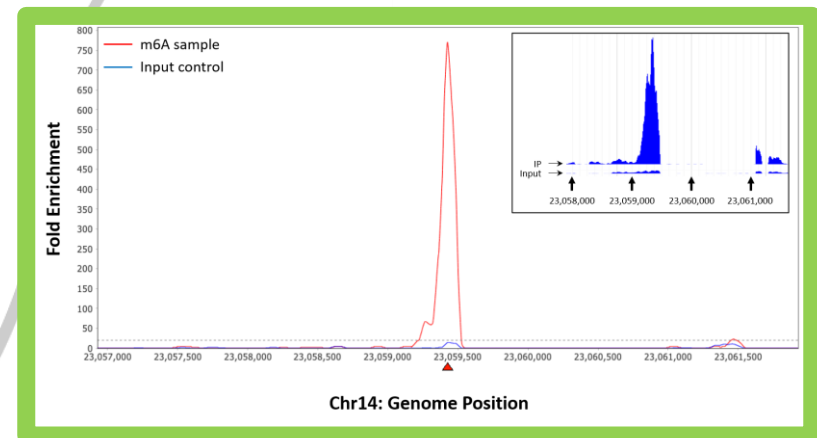
- Higher enrichment
- Lower input
- Reduced background
- Faster, more streamlined procedure

- **Cleavage under targets**

- Simultaneous fragmentation of immunocaptured RNA and cleavage/removal of any sequences in both ends of target m<sup>6</sup>A-containing sequences
- RNA regions occupied by antibody are unaffected

- **Generation of short RNA fragments only bound with anti-m<sup>6</sup>A antibody**

- High resolution mapping



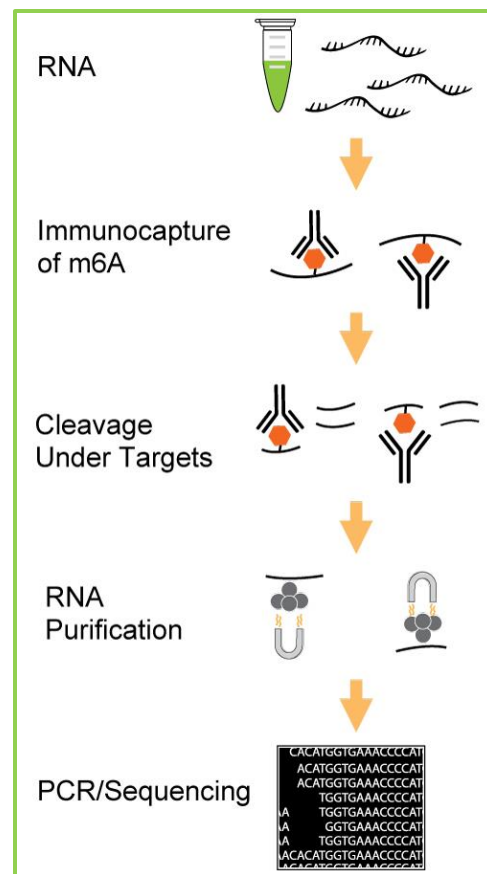
The m<sup>6</sup>A peak distribution within the human *ACIN1* transcript from samples processed with **CUT&RUN m<sup>6</sup>A MeRIP** (EpigenTek cat. #P-9018) correlates well with public database records (inset).



# P-9018 CUT&RUN MeRIP Procedure

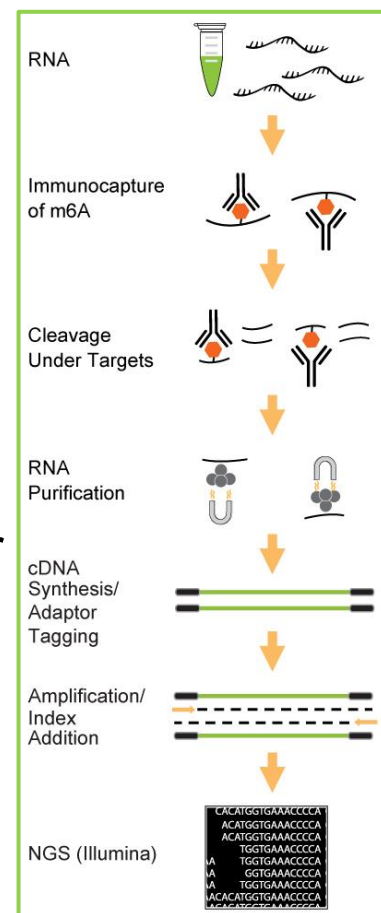
- Isolated and purified total RNA from any species
- m<sup>6</sup>A RNA immunocapture with highly-specific anti-m<sup>6</sup>A antibody (included in kit)
- Cleavage under targets (in situ)
- Wash away unbound RNA (via antibody affinity magnetic beads)
- Enriched m<sup>6</sup>A RNA release, purification (via RNA affinity magnetic beads), and elution

The enriched m<sup>6</sup>A RNA is now ready for downstream analysis (PCR, NGS)



# P-9016 CUT&RUN MeRIP-Seq Kit

- Same features as P-9018 **+ library prep**
- Reagents included for:
  - **Reverse transcription** of released enriched RNA
  - **Adaptor ligation**
  - **Amplification** of purified, adaptor-ligated cDNA for NGS library construction





# Key Features

- Fast, streamlined procedures
- Sonication-free fragmentation
- Cleavage and immunocapture processed **in same single-tube**
  - Minimized sample loss
  - Low input amount
- Unbound RNA cleavage/removal in situ using unique nucleic acid cleavage enzyme mix with low sequence bias
- RNA cleaved at both ends, directly adjacent to antibody binding site
  - Enrichment of shorter RNA fragments for more reliable identification of m<sup>6</sup>A RNA methylation sites
  - Minimized immunocapture/sequencing background
  - High-resolution mapping

For more information, please visit:

[www.epigentek.com](http://www.epigentek.com)

[www.epigentek.com/distributors](http://www.epigentek.com/distributors) (non-US customers)

Or contact your account associate today at:

**[sales@epigentek.com](mailto:sales@epigentek.com)**