

# A monoclonal antibody for transcriptome-wide N<sup>6</sup>-methyladenosine analysis

Ryan T. Fuchs<sup>1</sup>, Laurence Ettwiller<sup>1</sup>, Curtis Desilets<sup>2</sup>, Christopher J. Fry<sup>2</sup>, and G. B. Robb<sup>1</sup>

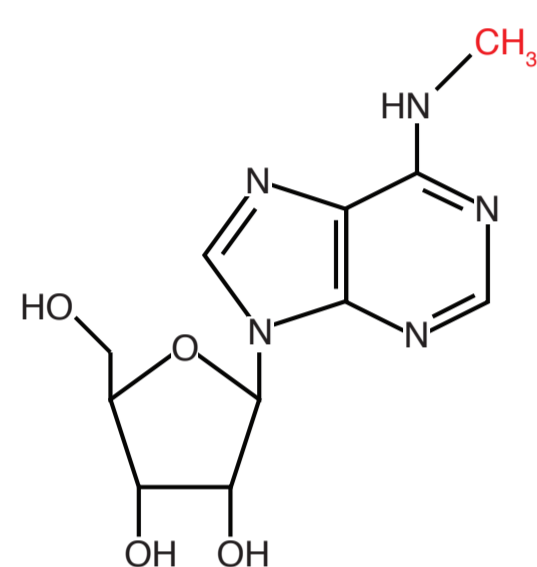
<sup>1</sup>New England Biolabs, Ipswich, MA 01938, USA

<sup>2</sup>Cell Signaling Technology, Beverly, MA 01915, USA

## Abstract

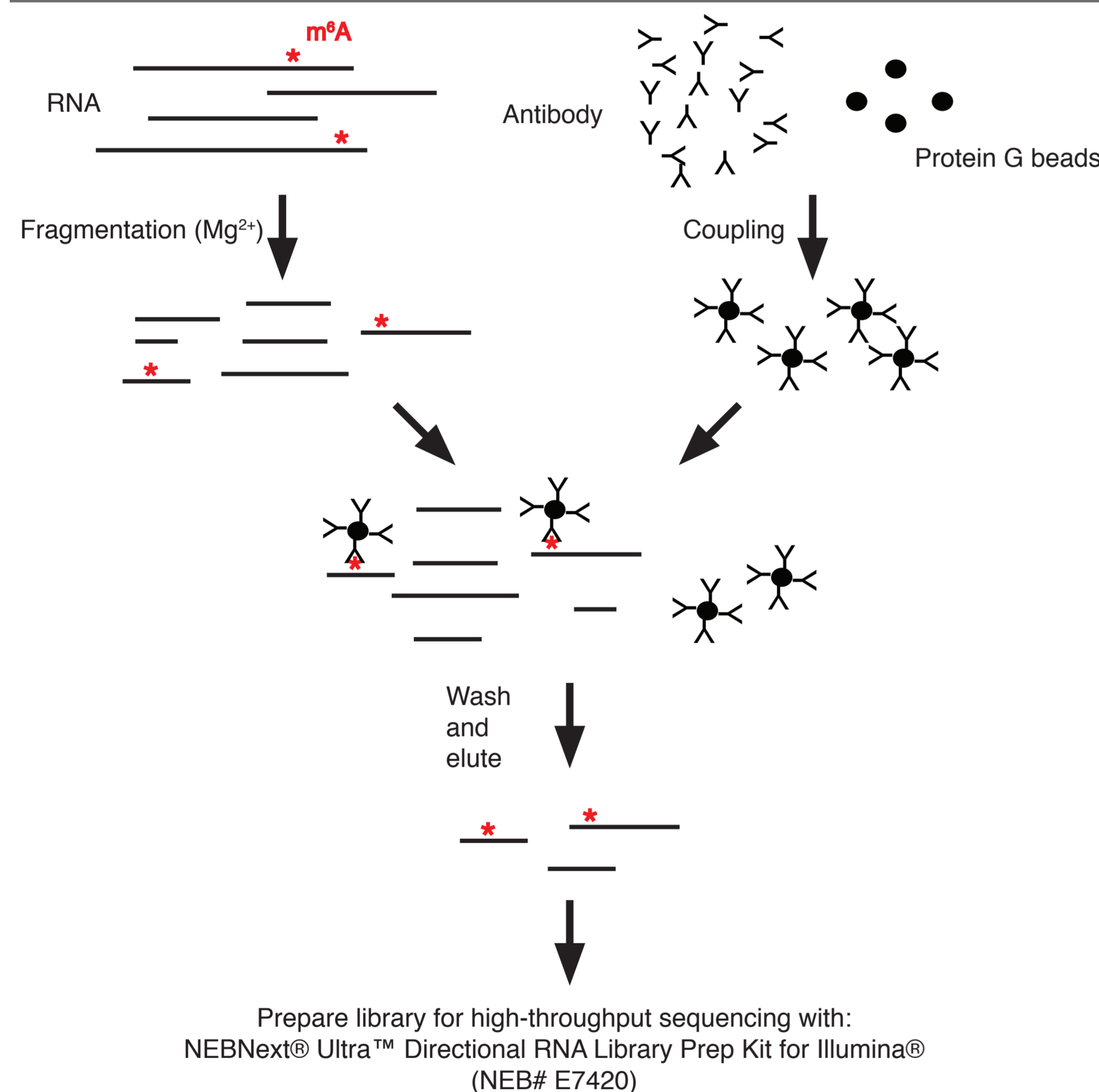
N<sup>6</sup>-methyladenosine (m<sup>6</sup>A) has been shown to be the most common base modification in eukaryotic messenger RNA (mRNA) other than the 7-methylguanosine cap. Recent m<sup>6</sup>A-RNA immunoprecipitation (m<sup>6</sup>A-RIP) with polyclonal antibodies combined with RNA high-throughput sequencing (RNA-seq) studies have identified the location of m<sup>6</sup>A sites in a transcriptome-wide manner in a variety of tissues and have started to analyze the function of m<sup>6</sup>A in mRNA (1-8). In humans m<sup>6</sup>A is most commonly associated with a sequence motif in the 3' UTR of mRNAs near stop codons and m<sup>6</sup>A modification is dependent upon a complex consisting of the methyltransferases METTL3+METTL14 and accessory proteins such as WTAP and KIAA1429. To further advance our understanding of m<sup>6</sup>A in RNA, it is important to continue improving the tools needed for m<sup>6</sup>A research. Here we present the generation of a new m<sup>6</sup>A-specific rabbit monoclonal antibody and its use in m<sup>6</sup>A-RIP-seq experiments.

## Antibody Development

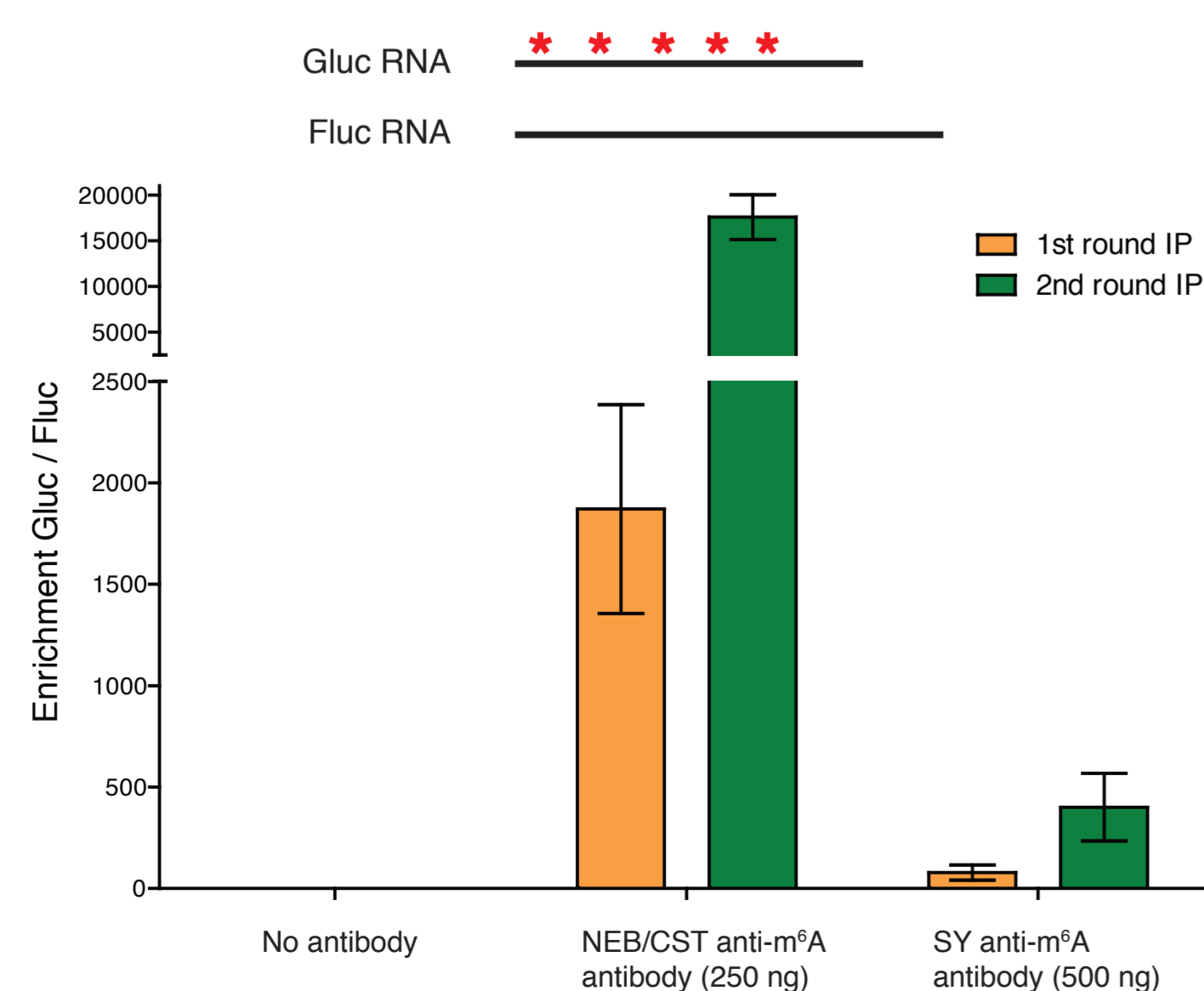


- N<sup>6</sup>-methyladenosine (m<sup>6</sup>A) was coupled to hemocyanin and introduced into rabbits to produce antibodies
- Dot blots and ELISA assays were used to show reactivity towards m<sup>6</sup>A, but no reactivity with unmodified adenosine or other modified nucleosides (m<sup>1</sup>A, 2'-OMe A)

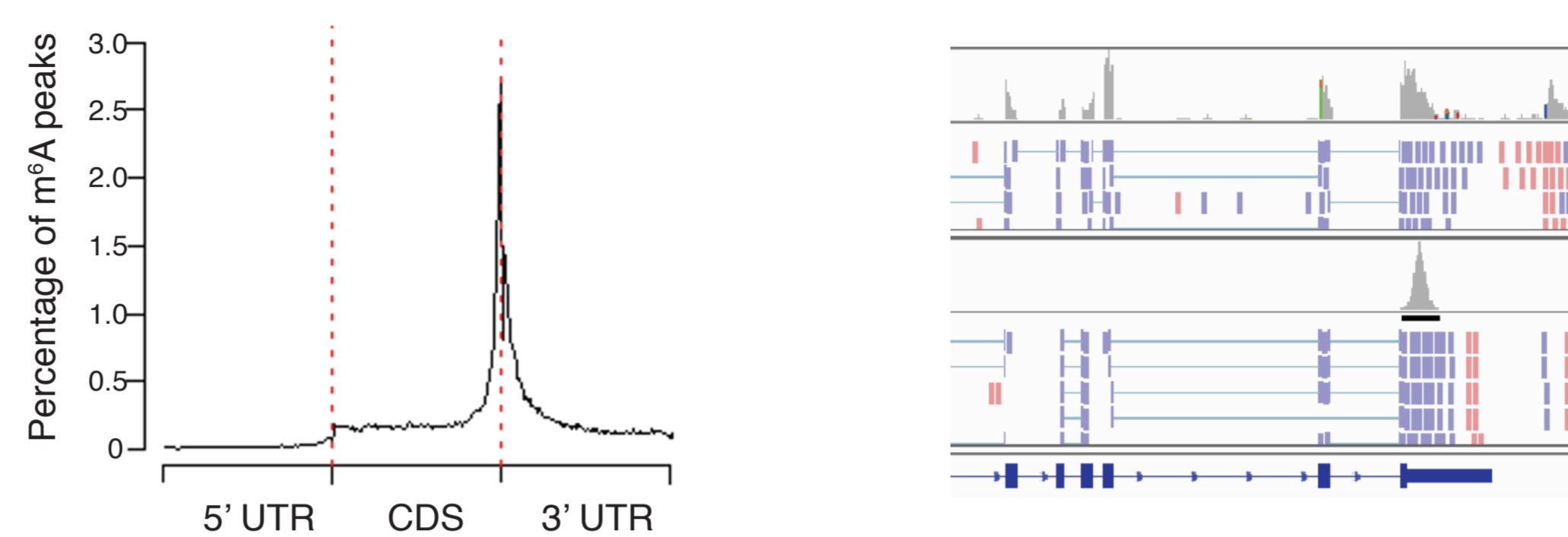
## RNA IP procedure



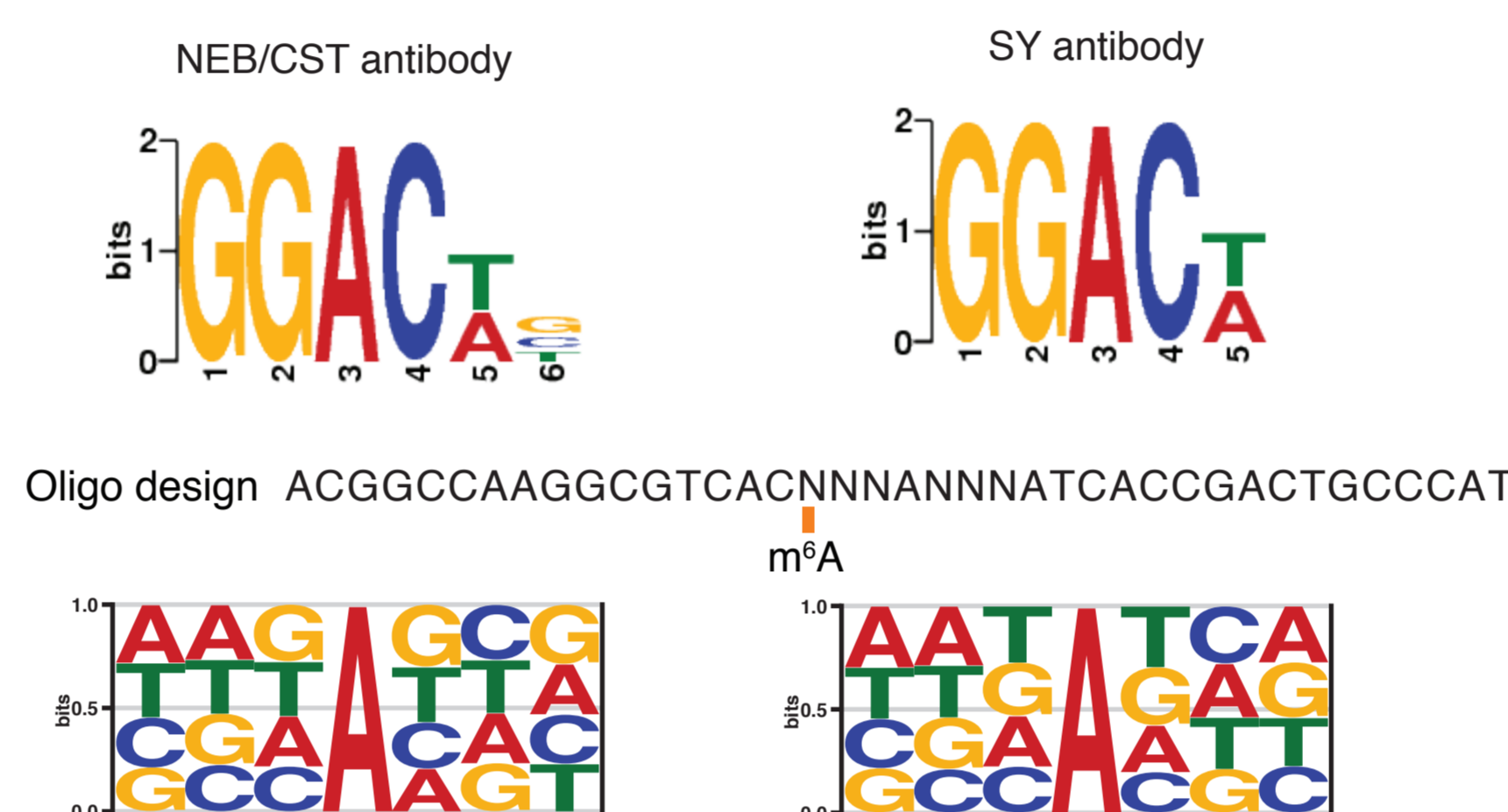
## RNA IP-qPCR with control RNAs



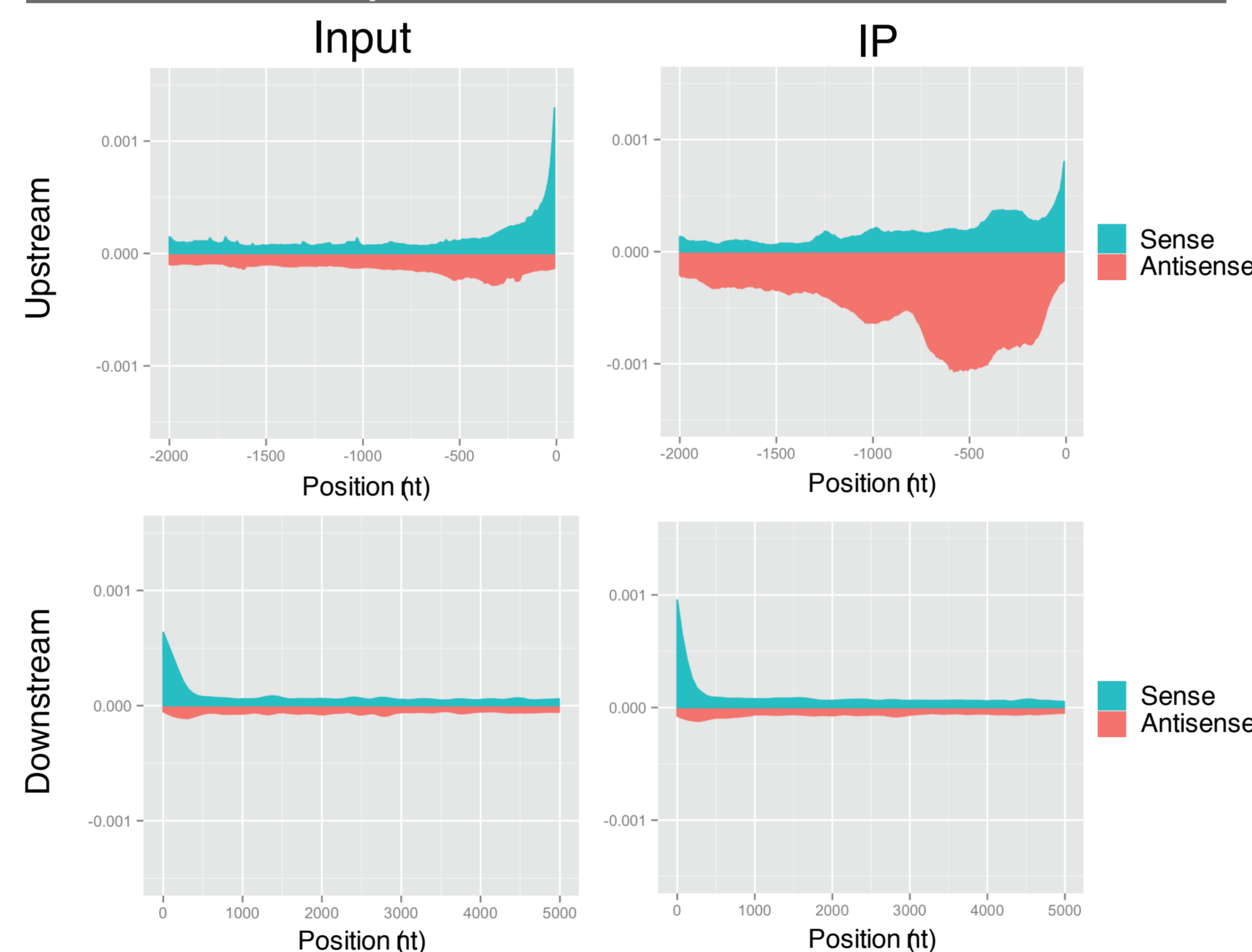
## RIP-seq using mouse brain polyA+ RNA



Motif analysis within m<sup>6</sup>A peaks near stop codons

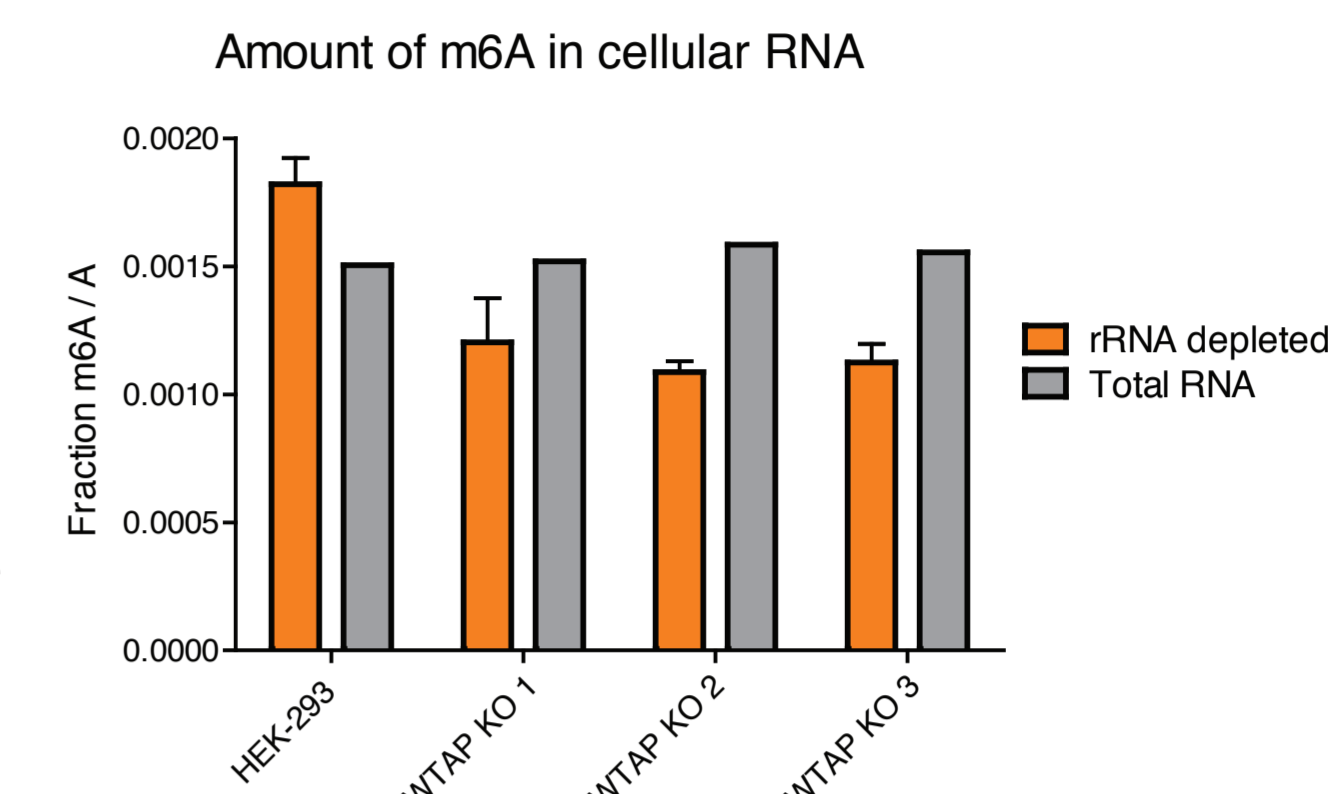


## RIP-seq reveals antisense enrichment



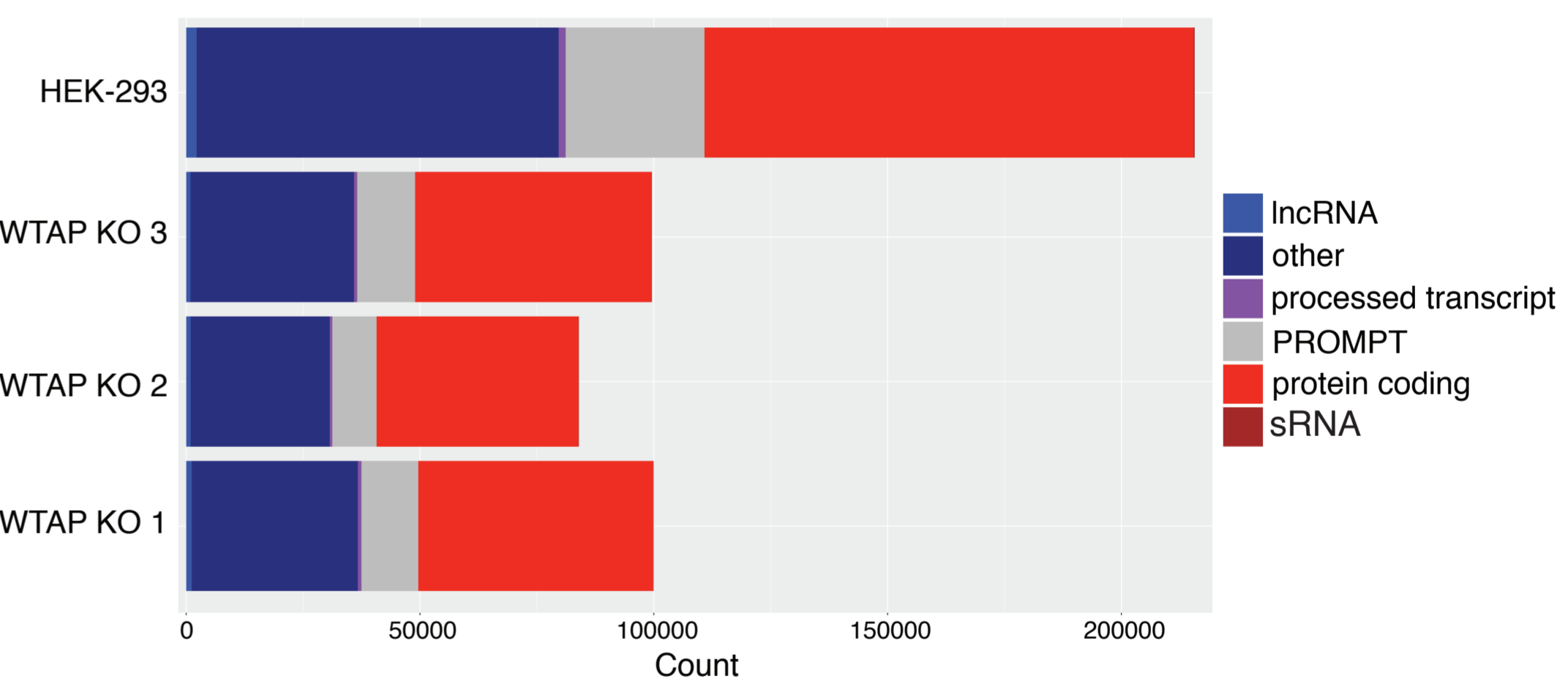
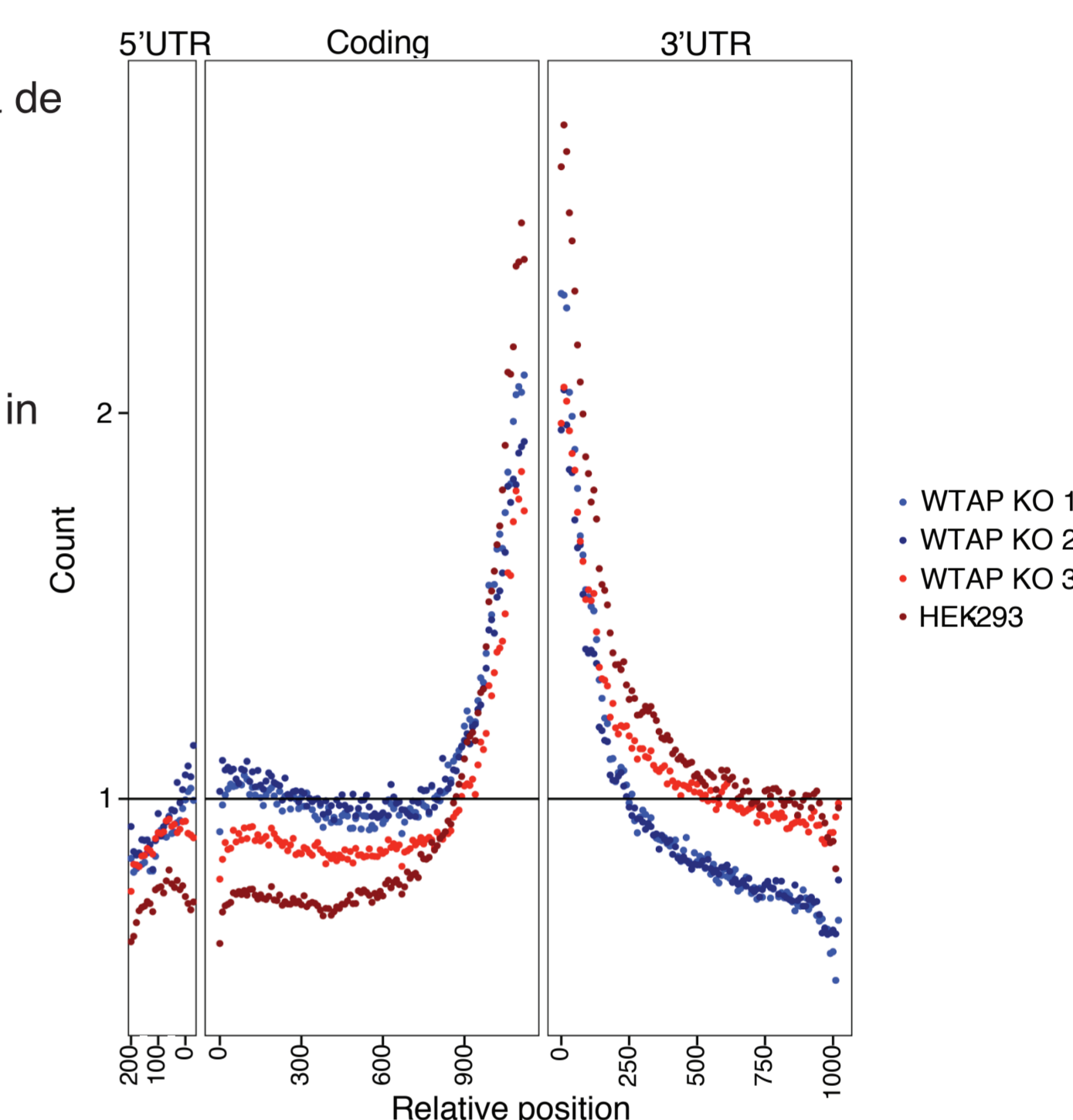
## RIP-seq using RNA from WTAP knockout cells

- HPLC analysis of RNA showed ~40-50% decrease in m<sup>6</sup>A in mRNA in WTAP null cells



- m<sup>6</sup>A RIP-seq showed a decrease of m<sup>6</sup>A near stop codons

- The distribution of m<sup>6</sup>A in classes of RNAs is unchanged in WTAP null cells



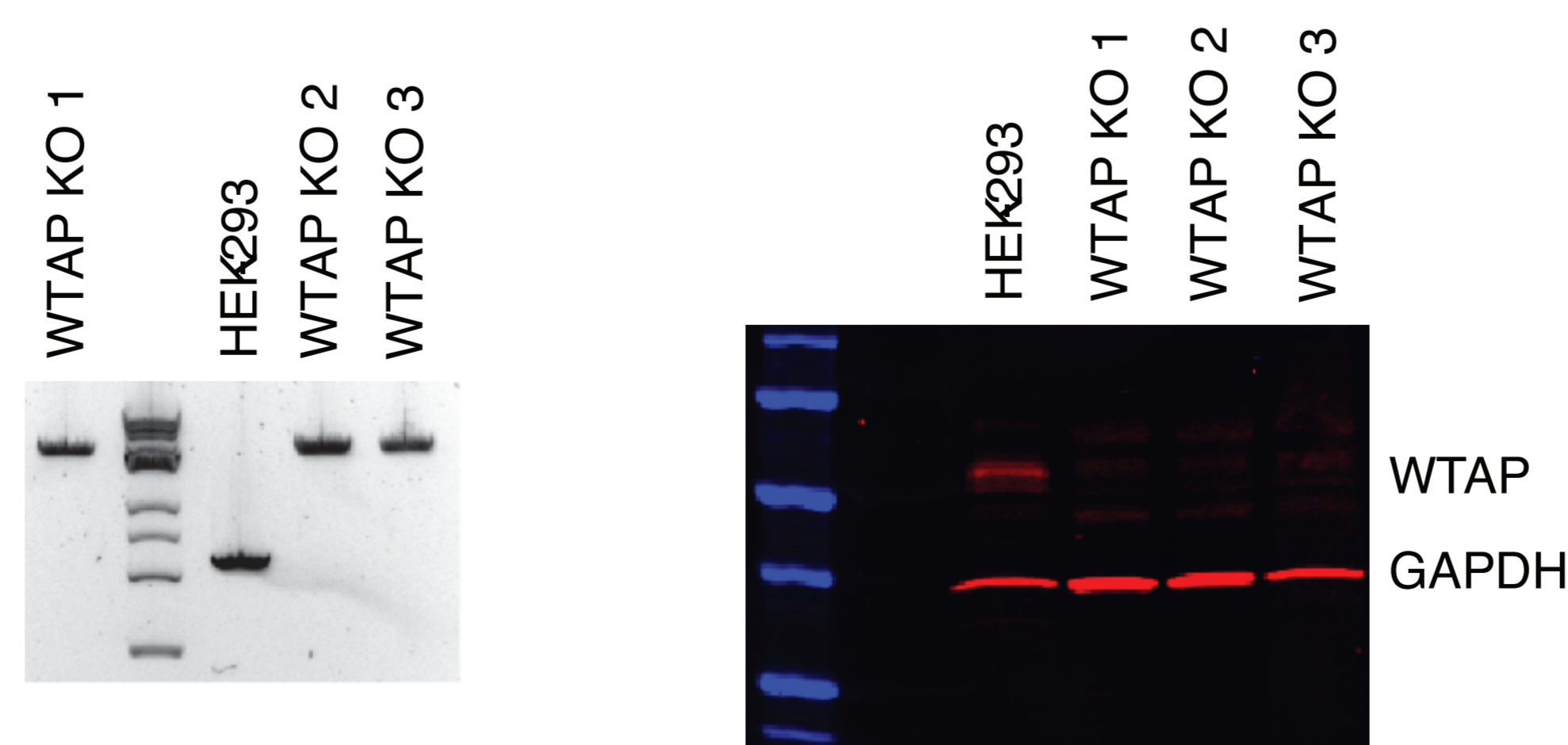
## Conclusions

- We have developed and validated a monoclonal antibody specific for m<sup>6</sup>A.
- The monoclonal antibody results in better enrichment of m<sup>6</sup>A-containing control RNA than a currently available polyclonal antibody.
- The monoclonal antibody performs well in m<sup>6</sup>A-RIP-seq and produces results consistent with published m<sup>6</sup>A-RIP-seq data. In addition, our analysis revealed enrichment of antisense RNAs.
- WTAP knockout strains show a reduction in RNA m<sup>6</sup>A levels and reduced enrichment of m<sup>6</sup>A modified sites after RIP-seq, but the overall pattern of m<sup>6</sup>A sites does not change.
- This suggests that WTAP is necessary for full methyltransferase complex activity, but not significantly involved in determining which sites are methylated.

## References

- Schwartz S, Mumbach MR, Jovanovic M, et al. Cell Rep. 2014. 8:284-296.
- Schwartz S, Agarwala SD, Mumbach MR, et al. Cell. 2013. 155:1409-1421.
- Dominissini D, Moshitch-Moshkovitz S, Schwartz S, et al. Nature. 2012. 485:201-206.
- Meyer KD, Saletore Y, Zumbo P, Elemento O, Mason CE, and Jaffrey SR. Cell. 2012. 1635-1646.
- Wang X, Lu Z, Gomez A, et al. Nature. 2014. 505:117-120.
- Dominissini D, Moshitch-Moshkovitz S, Salmon-Divon M, et al. Nat. Protoc. 2013. 8:176-189.
- Fustin J, Doi M, Yamaguchi Y, et al. Cell. 2013. 155:793-806.
- Wang Y, Li Y, Toth JI, Petroski MD, Zhang Z, and Zhao JC. Nat. Cell Biol. 2014. 16:191-198.

## Disruption of WTAP via Cas9 genome editing



- PCR from genomic DNA shows rearrangement at the WTAP locus
- Western blot analysis shows complete loss of WTAP protein expression