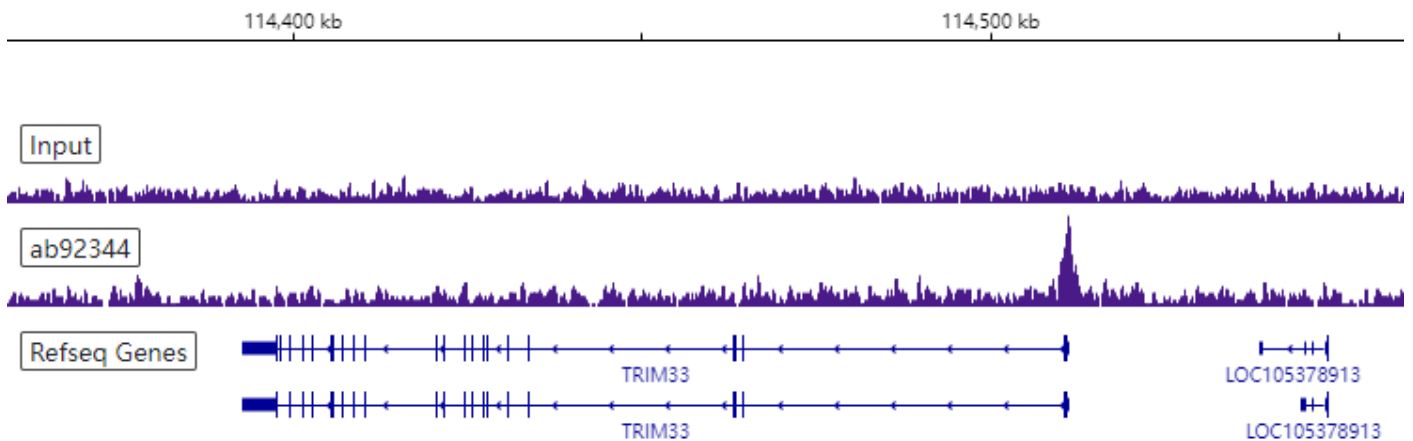


**Anti-RBBP4 antibody [EPR3412] (ab92344)**

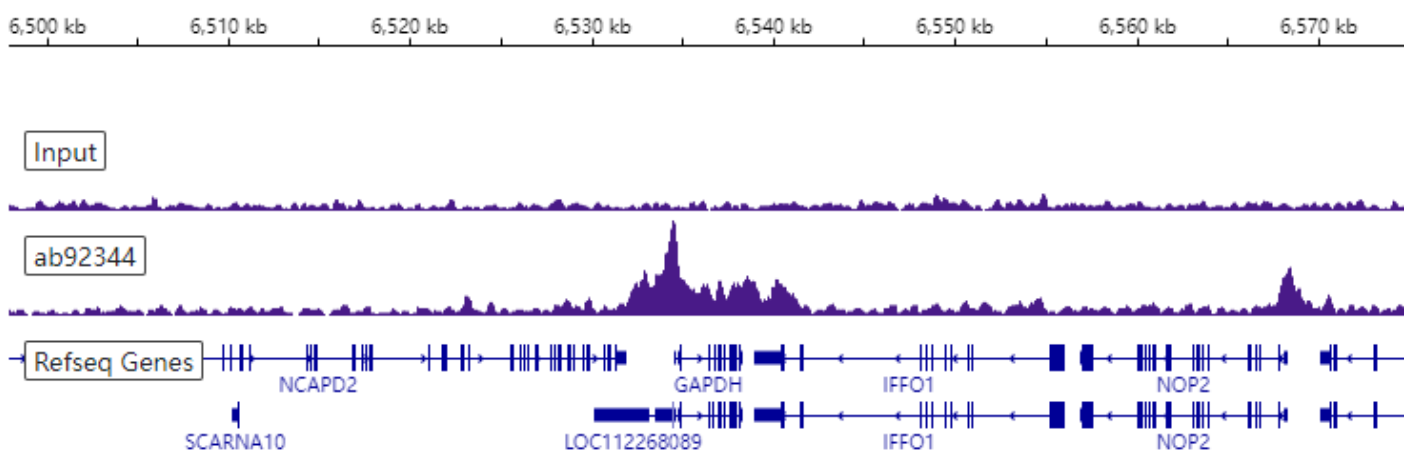
|                  |                       |
|------------------|-----------------------|
| Target           | RBBP4                 |
| Sample           | HeLa                  |
| Treatment        | None                  |
| Antibody amount  | 8 µg                  |
| Cell number      | 1×10 <sup>7</sup>     |
| Sequencing       | NovaSeq 6000 (PE 150) |
| Peak calling     | MACS 2.1.0            |
| Reference genome | hg38                  |

**Chromosome 1 (active region)**



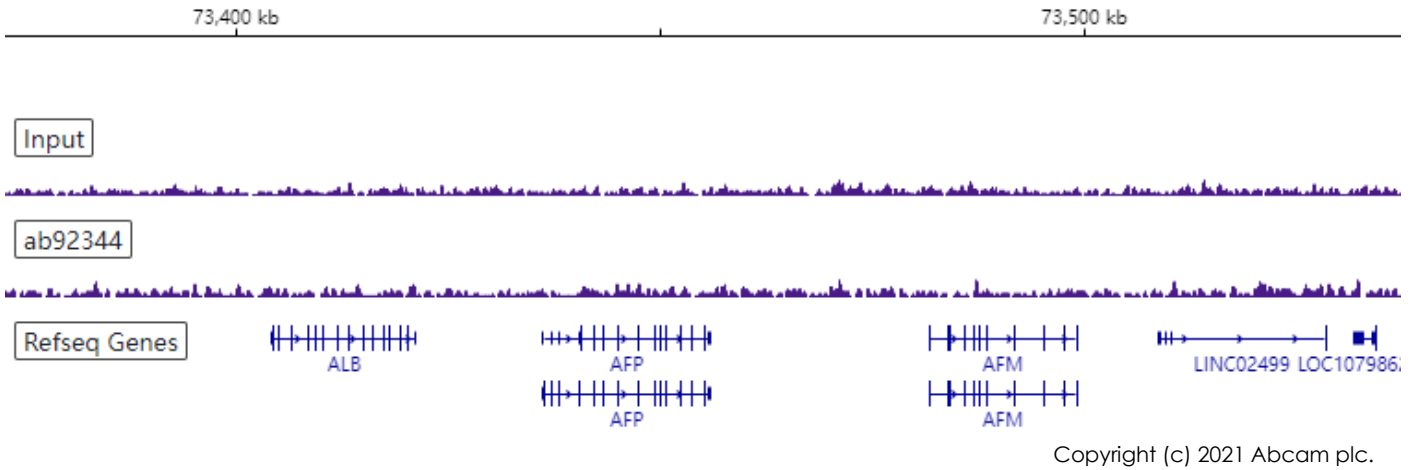
Copyright (c) 2021 Abcam plc.

**Chromosome 12 (active region)**

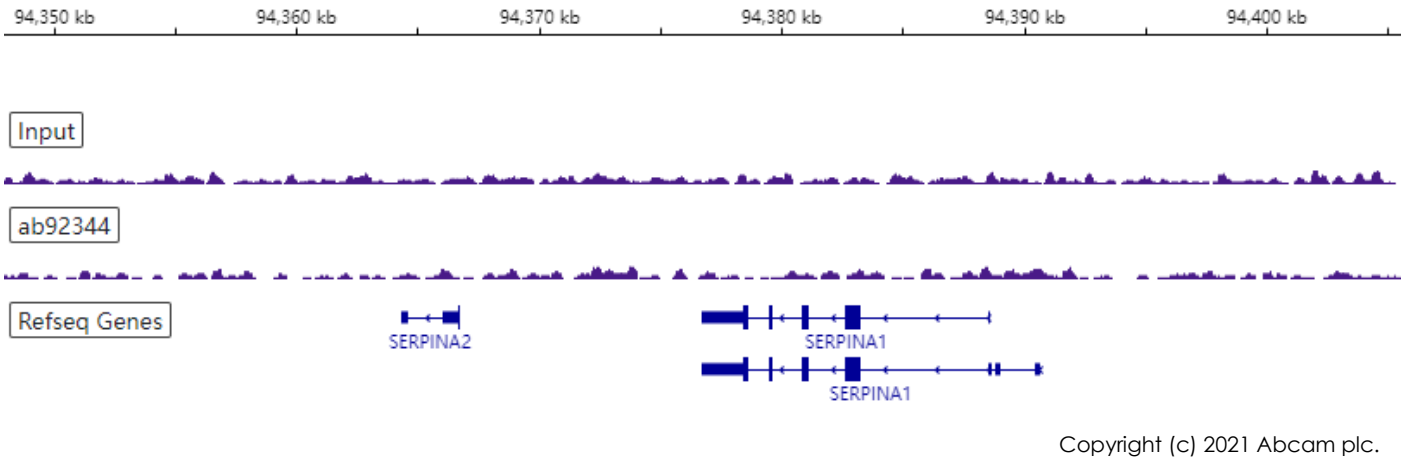


Copyright (c) 2021 Abcam plc.

## Chromosome 4 (inactive region)



## Chromosome 14 (inactive region)



## Chromosome 11 (inactive region)

