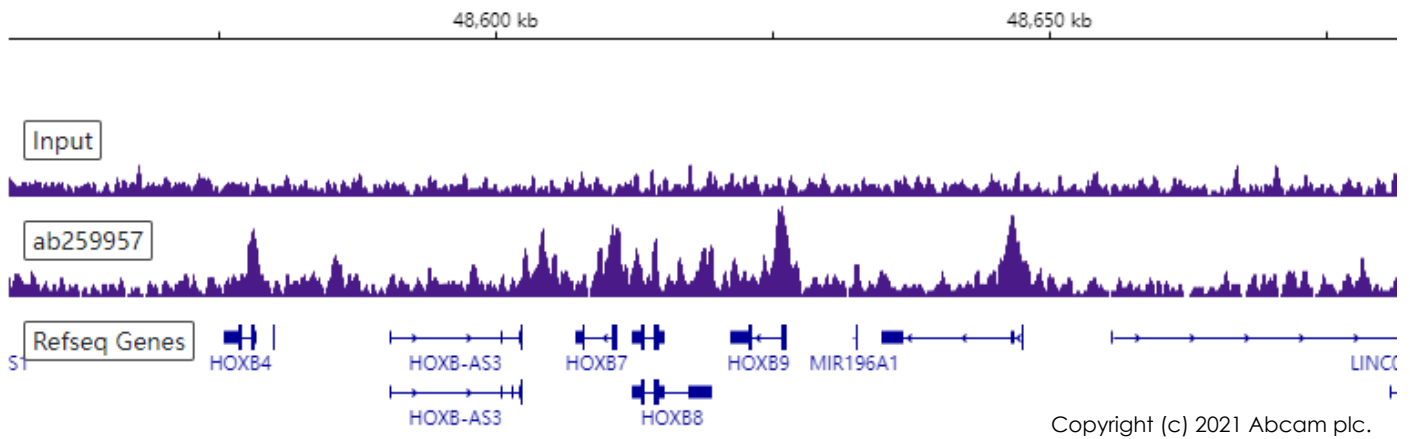


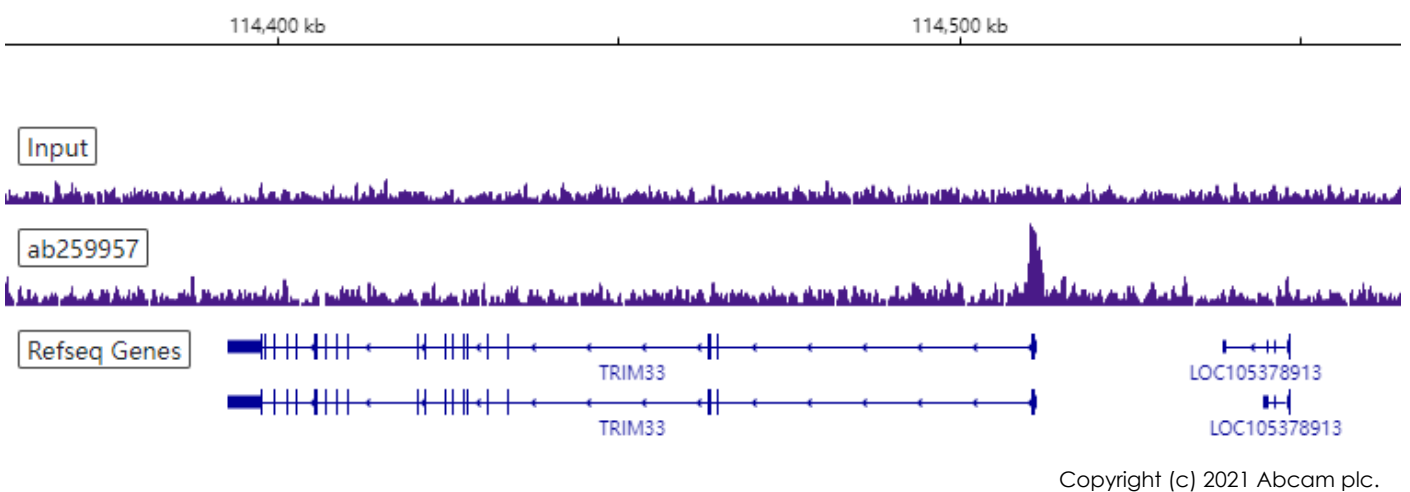
Anti-RBBP7 antibody [EPR23796-74] - ChIP Grade (ab259957)

Target	RBBP7
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

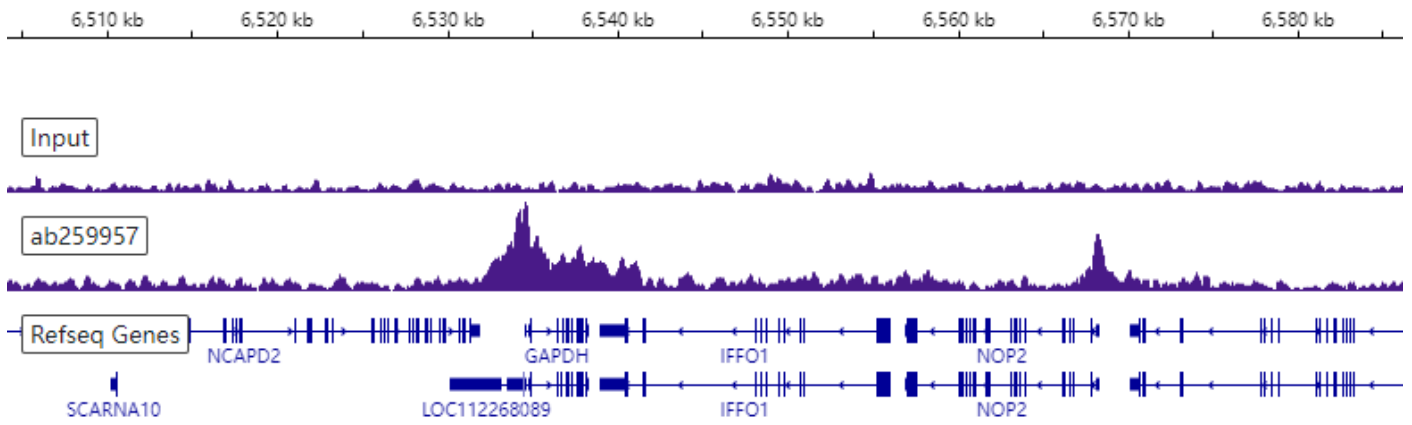
Chromosome 17 (active region)



Chromosome 1 (active region)

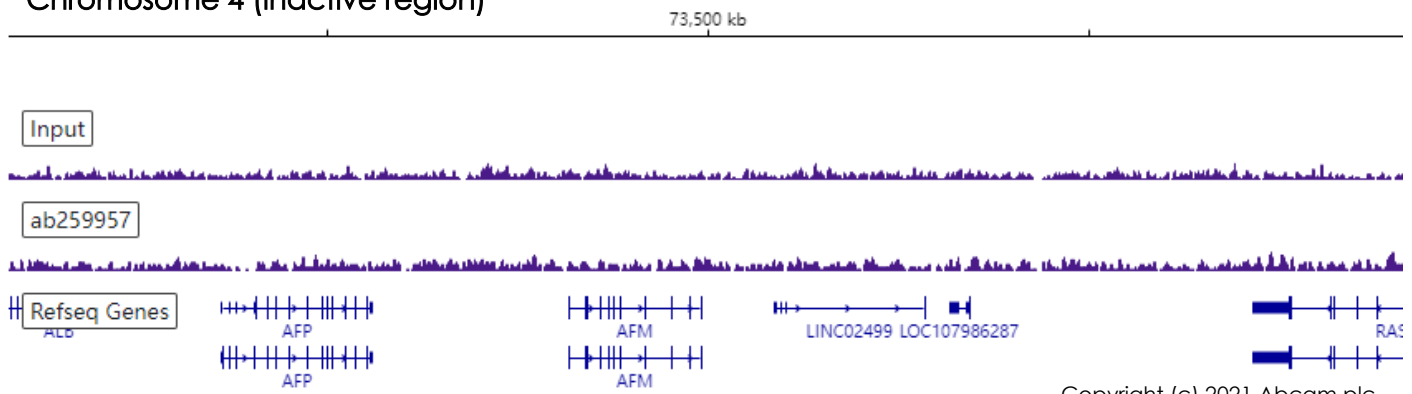


### Chromosome 12 (active region)



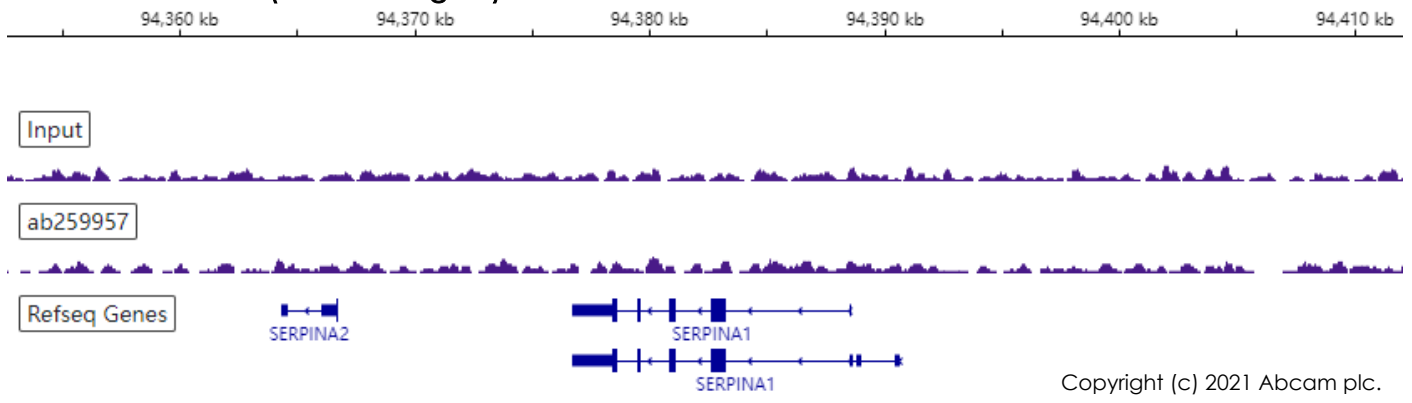
Copyright (c) 2021 Abcam plc.

### Chromosome 4 (inactive region)



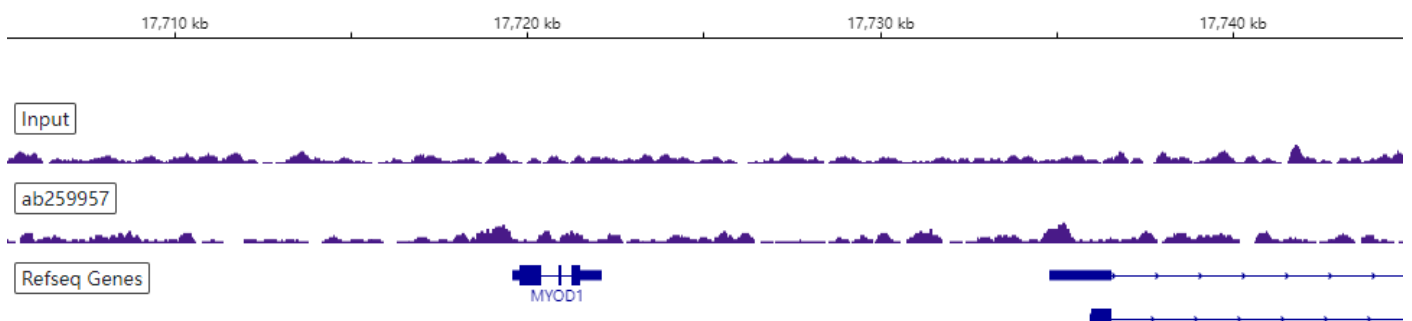
Copyright (c) 2021 Abcam plc.

### Chromosome 14 (inactive region)



Copyright (c) 2021 Abcam plc.

### Chromosome 11 (inactive region)



Copyright (c) 2021 Abcam plc.