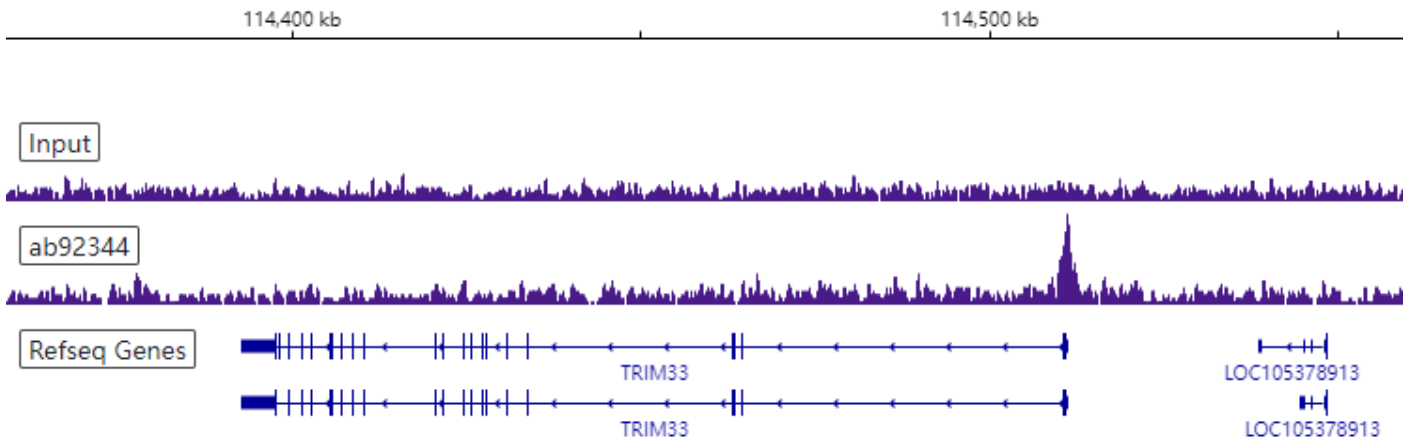


Anti-RBBP4 antibody [EPR3412] (ab92344)

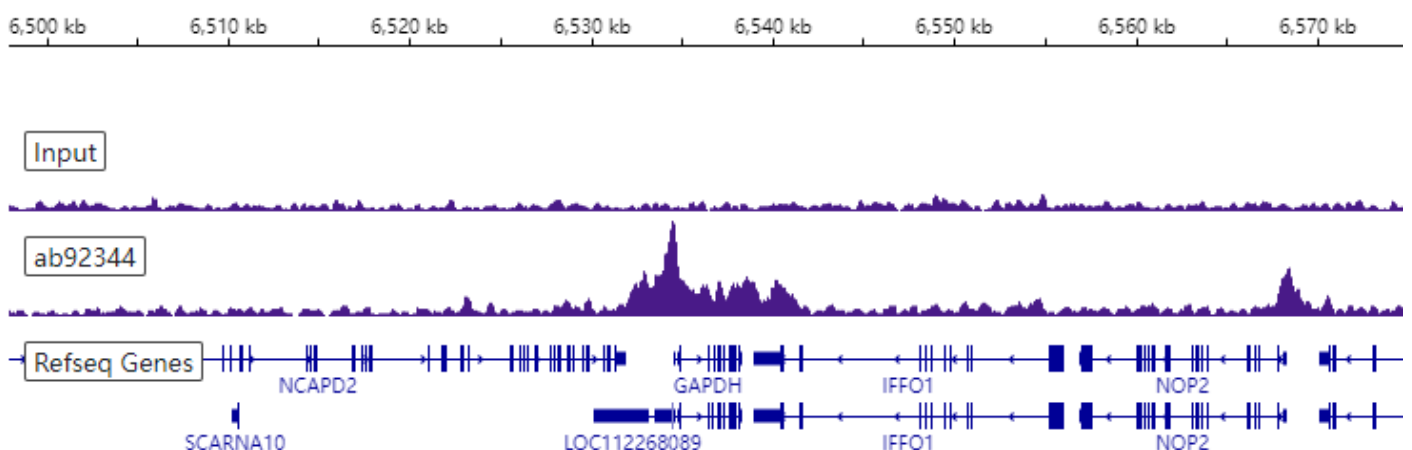
Target	RBBP4
Sample	HeLa
Treatment	None
Antibody amount	8 µg
Cell number	1×10 ⁷
Sequencing	NovaSeq 6000 (PE 150)
Peak calling	MACS 2.1.0
Reference genome	hg38

Chromosome 1 (active region)



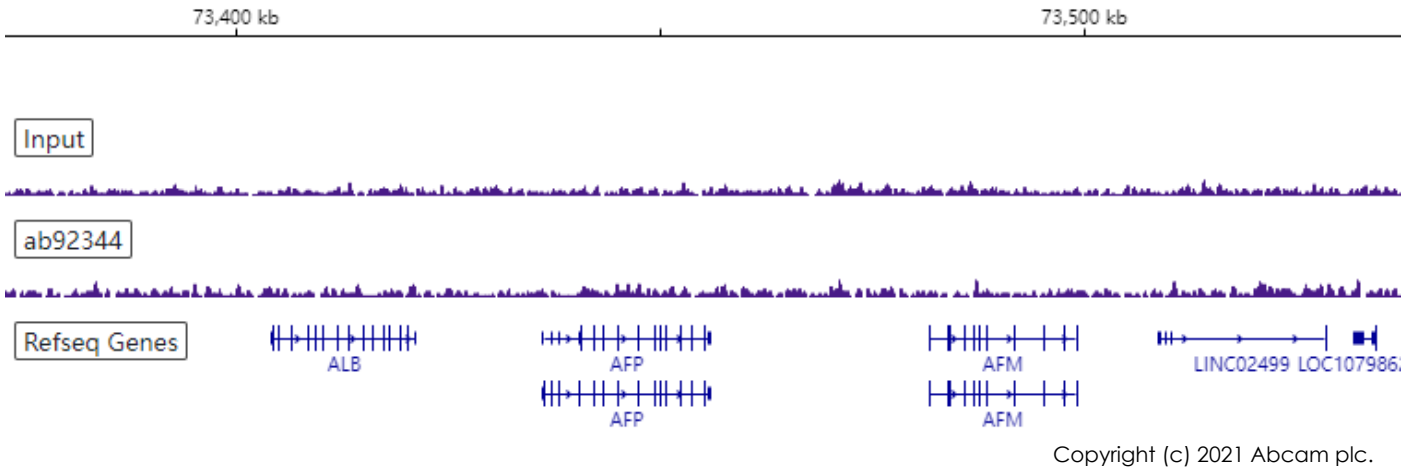
Copyright (c) 2021 Abcam plc.

Chromosome 12 (active region)

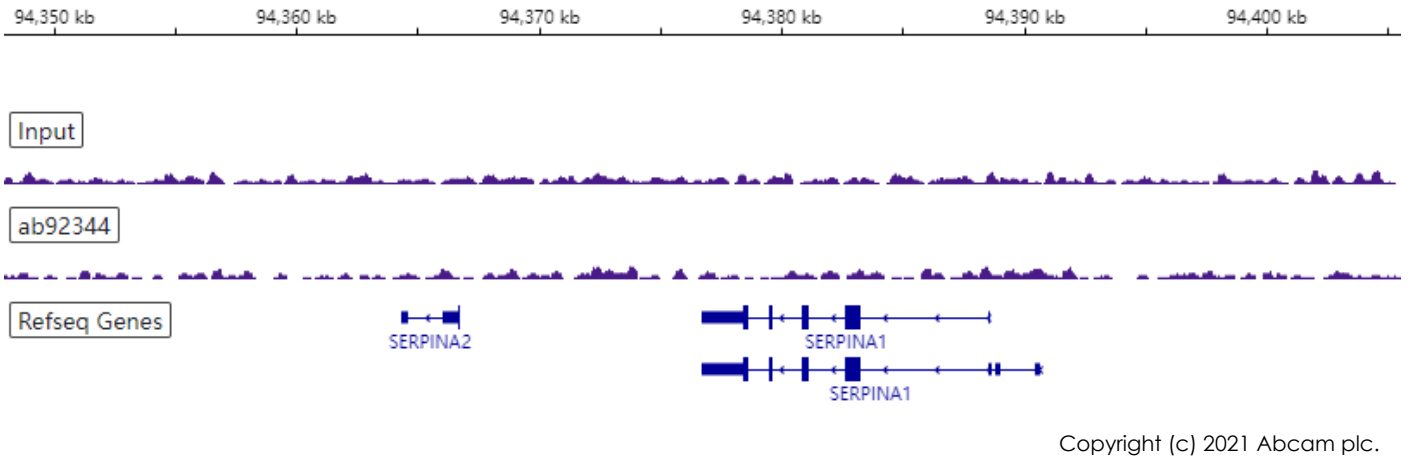


Copyright (c) 2021 Abcam plc.

Chromosome 4 (inactive region)



Chromosome 14 (inactive region)



Chromosome 11 (inactive region)

