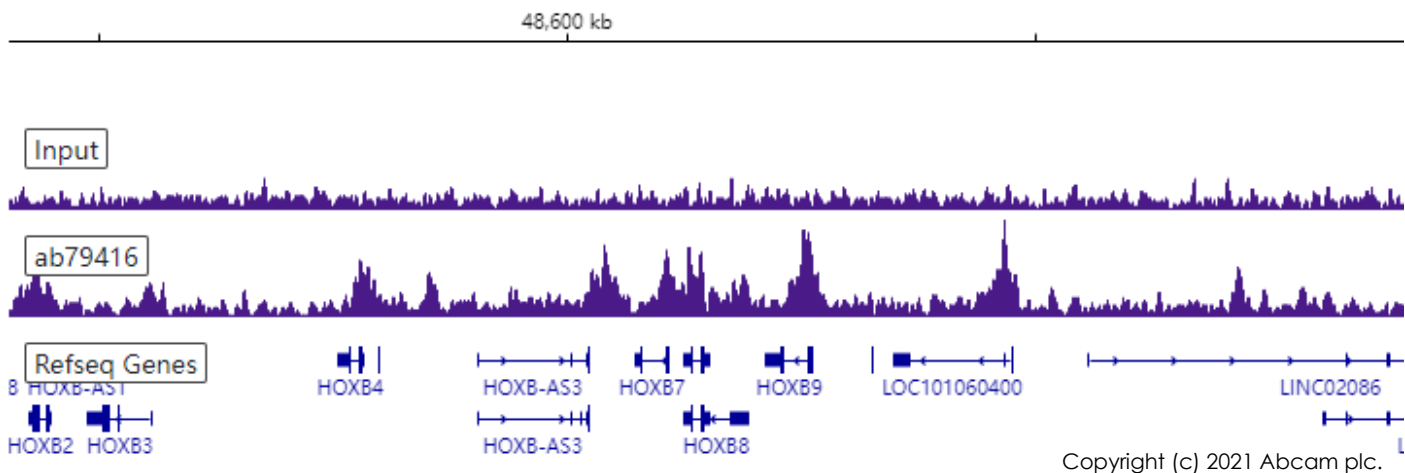


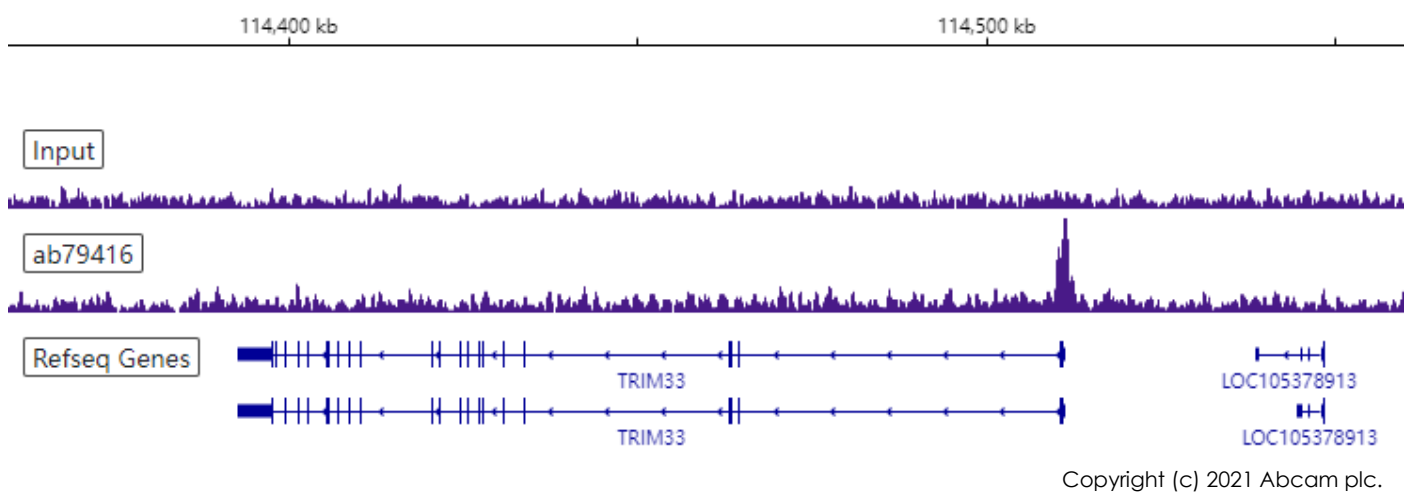
Anti-RBBP4 antibody [EPR3411] - ChIP Grade (ab79416)

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

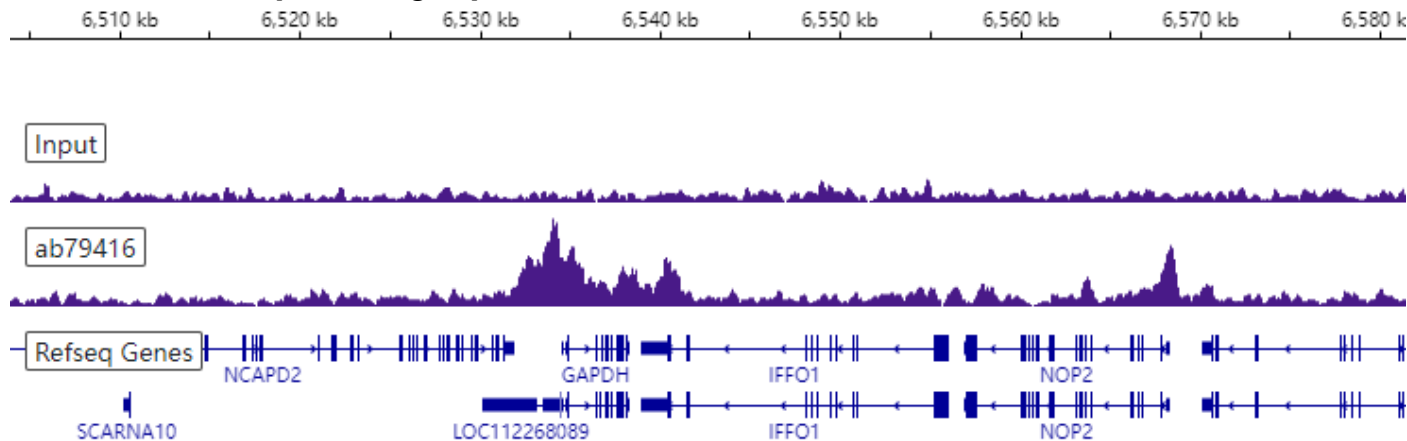
Chromosome 17 (active region)



Chromosome 1 (active region)

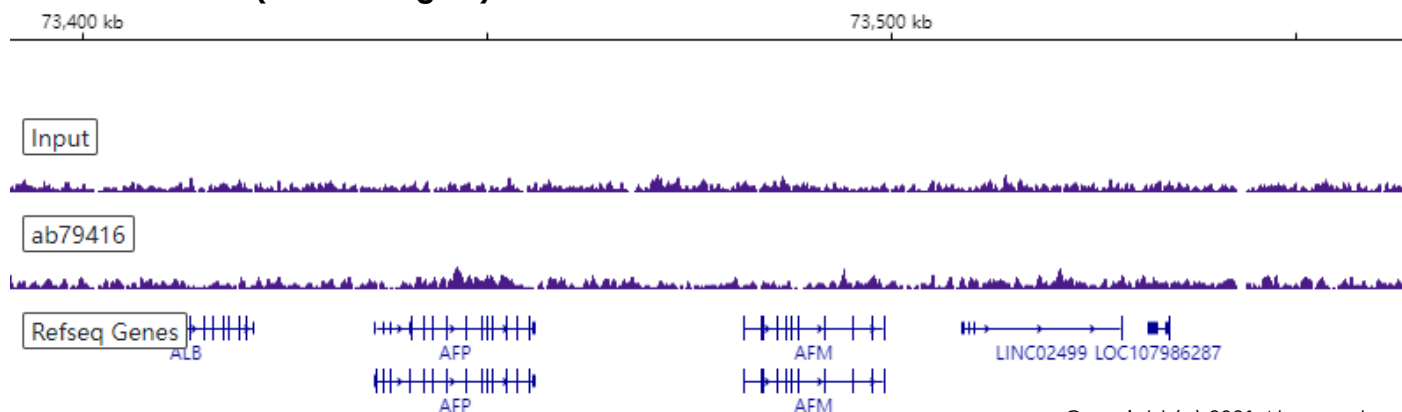


Chromosome 12 (active region)



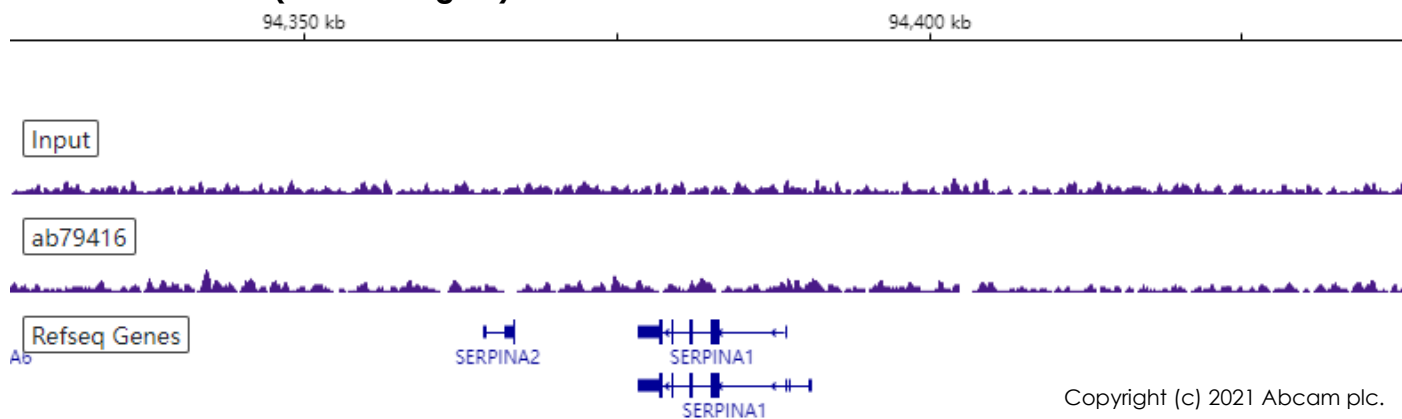
Copyright (c) 2021 Abcam plc.

Chromosome 4 (inactive region)



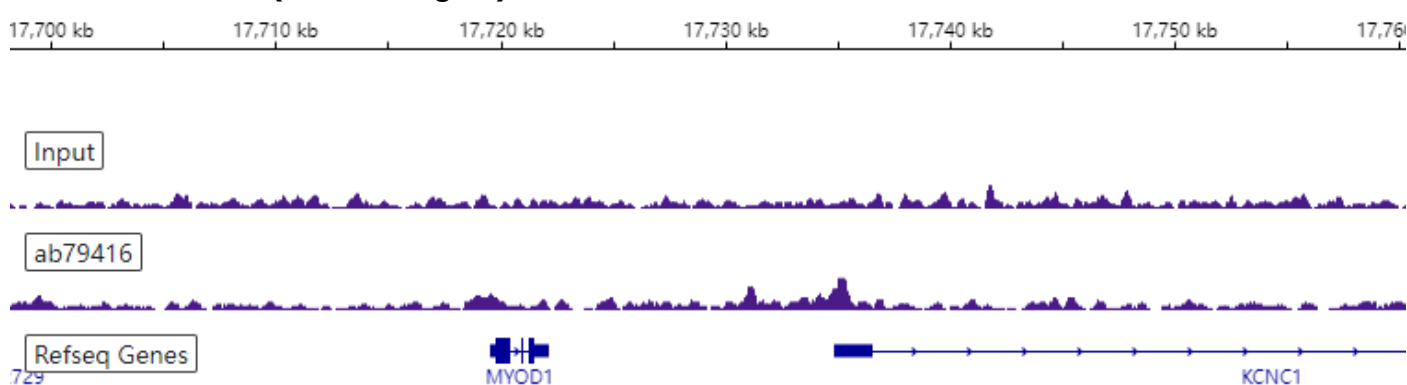
Copyright (c) 2021 Abcam plc.

Chromosome 14 (inactive region)



Copyright (c) 2021 Abcam plc.

Chromosome 11 (inactive region)



Copyright (c) 2021 Abcam plc.