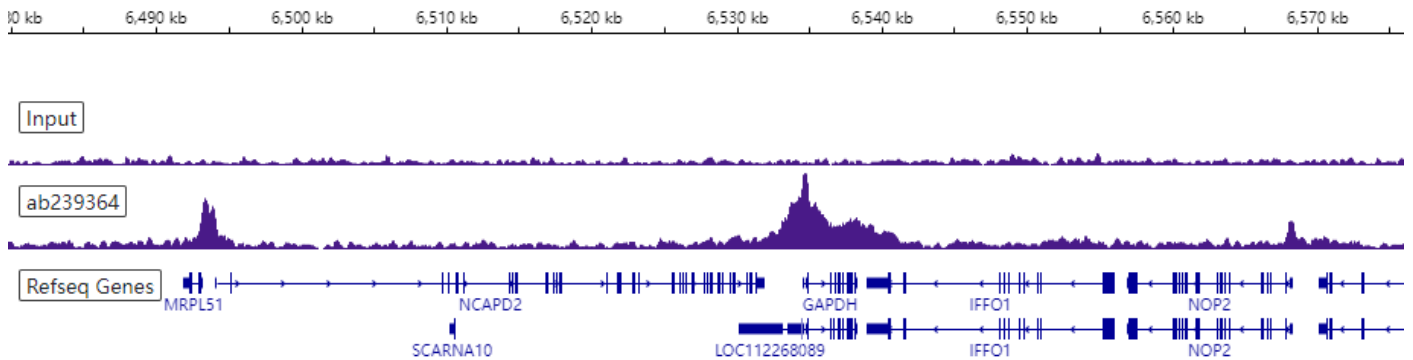


Anti-Cdk9 antibody [EPR22956-37] - ChIP Grade (ab239364)

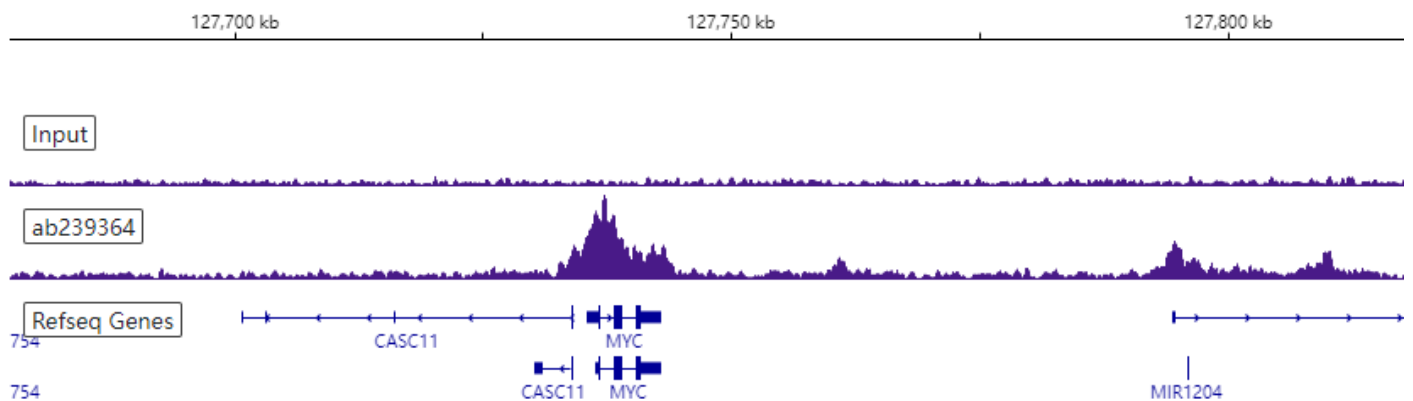
Target	Cdk9
Sample	HeLa
Treatment	None
Antibody amount	8 µg
Cell number	1×10 ⁷
Sequencing	NovaSeq 6000 (PE 150)
Peak calling	MACS 2.1.0
Reference genome	hg38

Chromosome 12 (active region)



Copyright (c) 2021 Abcam plc.

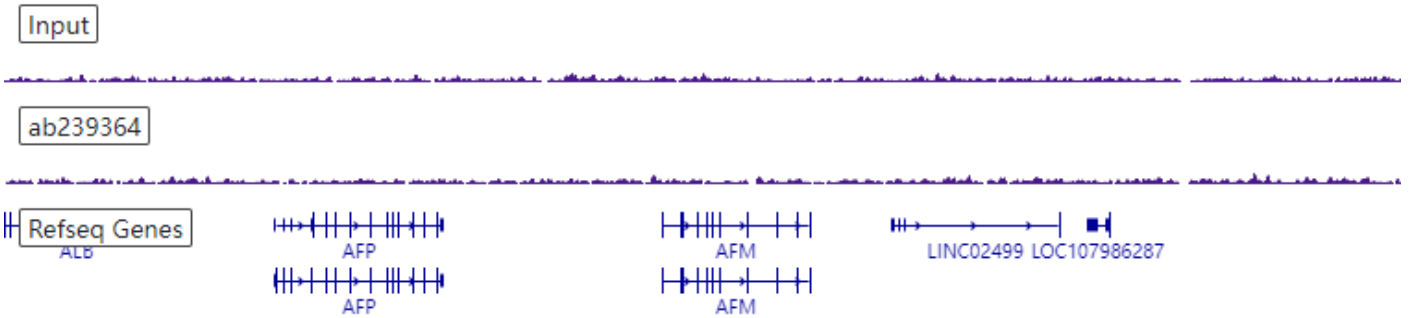
Chromosome 8 (active region)



Copyright (c) 2021 Abcam plc.

Chromosome 4 (inactive region)

73,500 kb

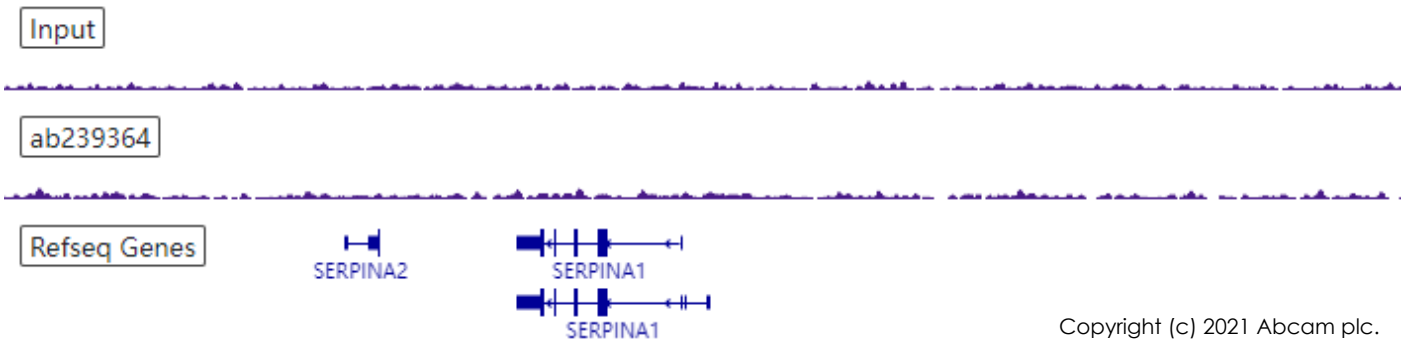


Copyright (c) 2021 Abcam plc.

Chromosome 14 (inactive region)

94,350 kb

94,400 kb



Copyright (c) 2021 Abcam plc.

Chromosome 11 (inactive region)

17,710 kb

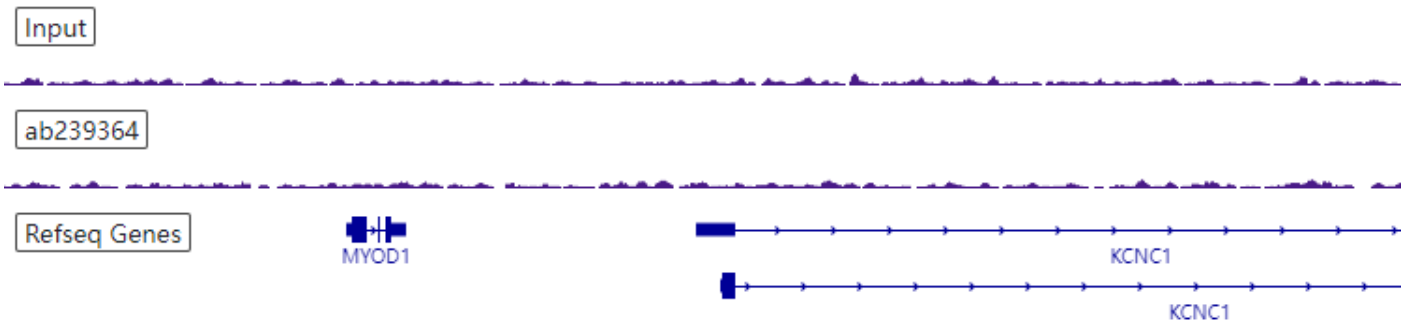
17,720 kb

17,730 kb

17,740 kb

17,750 kb

17,760 kb



Copyright (c) 2021 Abcam plc.