

## Androgen Receptor antibody (pAb)

**Catalog Nos:** 39781, 39081, 39782

**RRID:** AB\_2793341

**Isotype:** IgG

**Application(s):** ChIP, ChIP-Seq, WB

**Reactivity:** Human, Mouse

**Volumes:** 100 µl, 50 µl, 10 µl

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Concentration:** 1 µg/µl

**Molecular Weight:** 110 kDa

**Background:** Androgen Receptor (AR, NR3C4) is a nuclear receptor that facilitates signaling of testosterone and other androgenic hormones in the cytoplasm and translocates into the nucleus upon hormone binding. AR serves as a DNA binding activator of transcription of hormone responsive genes. There are two forms of the Androgen Receptor, the AR-A form, lacking the N-terminal 187 amino acids and the full length AR-B form. Androgen Receptor is involved in the development of both primary and secondary male sexual characteristics and abnormal expression of AR is linked to prostate cancer.

**Immunogen:** This Androgen Receptor antibody was raised against a peptide in the N-terminal region of human Androgen receptor.

**Buffer:** PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

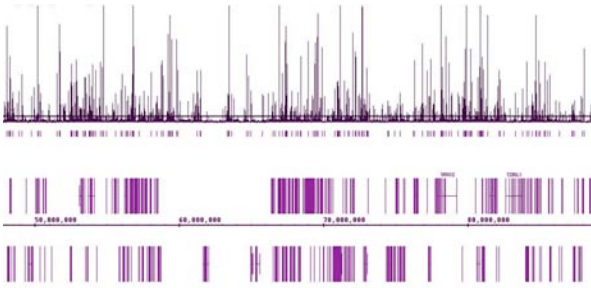
ChIP: 5 µg per ChIP

ChIP-Seq: 5 µg each

WB: 1 - 2 µg/ml dilution

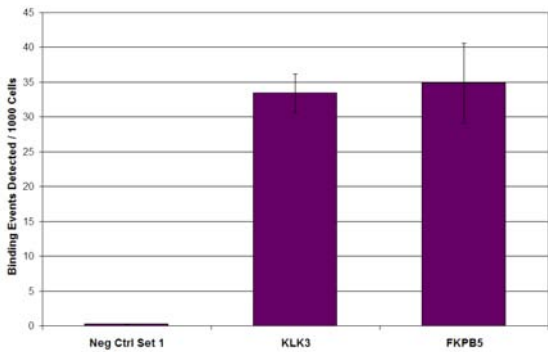
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



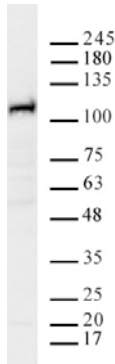
**Androgen Receptor antibody (pAb) tested by ChIP-Seq.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with chromatin from a human prostate cancer cell line (3 million cells). ChIP DNA was sequenced on the Illumina GAII and 22 million sequence tags were mapped to identify Androgen Receptor binding sites. The image shows hundreds of strong binding sites dispersed across the right arm of human chromosome 16.



**Androgen Receptor antibody (pAb) tested by ChIP.**

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 µg of VCAP60 cell chromatin and 10 µg of Androgen receptor antibody. ChIP DNA was used in qPCR with the negative control primer pairs or gene-specific primer pairs as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.



**Androgen Receptor antibody (pAb) tested by Western blot.**

Nuclear extract (20 µg) of LnCaP cells probed with Androgen Receptor antibody (pAb) (2 µg/ml).