

## NR2F2 / COUP2 antibody (pAb)

**Catalog Nos:** 61213, 61214

**RRID:** AB\_2614942

**Isotype:** IgG

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

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 50 kDa

**Background:** NR2F2 / COUP2 (ARP1, TFCOUP2, COT2, COUP-TF2) is a member of the nuclear receptor superfamily related to the COUP transcription factor. It interacts with several nuclear receptor co-activator proteins to activate transcription. NR2F2 / COUP2 plays important roles in signaling pathways in multiple endocrine systems and is crucial for differentiation of the trophoblast during development.

**Immunogen:** This NR2F2 / COUP2 antibody was raised against a peptide within the N-terminal region of human NR2F2.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Validated Applications:

ChIP: 5 µl per ChIP

ChIP-Seq: 5 µl each

WB: 1:500 - 1:2,000 dilution

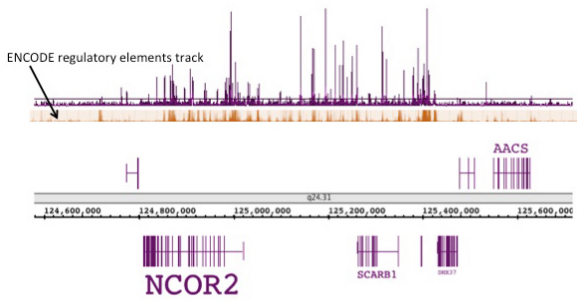
Published Applications:

ChIP

For optimal results, primary antibody incubations should be performed at 4°C. The addition of 0.05% Tween 20 to all Blotto solutions may also reduce background. Individual optimization may be required.

**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.



#### NR2F2 / COUP2 (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with chromatin from 4.5 million human endometrial stromal cells and 5  $\mu$ l of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 28 million sequence tags were mapped to identify NR2F2 binding sites. The purple track shows NR2F2 binding across a 1.2 million bp region on chromosome 12. Binding sites showed good correlation with the ENCODE active regulatory elements track (yellow).

#### NR2F2 / COUP2 (pAb) tested by Western blot.

Detection of NR2F2 / COUP2 by Western blot. The analysis was performed using 20  $\mu$ g MCF-7 nuclear extract and NR2F2 / COUP2 (pAb) at a 1:500 dilution.

