

## NCoA4 / ELE1 antibody (pAb)

**Catalog Nos:** 39895, 39896

**RRID:** AB\_2615019

**Isotype:** Serum

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** None

**Host:** Rabbit

**Molecular Weight:** 85 kDa

**Background:** NCoA4 (Nuclear receptor co-activator 4; also known as ELE1 and ARA70) is a transcriptional co-activator that enhances the transcriptional activity of several nuclear receptors, including androgen receptor and peroxisome proliferator-activated receptor (PPAR) gamma. It also interacts with the androgen receptor and the retinoid X receptor (RXR) in a ligand-dependent manner.

**Immunogen:** This NCoA4 / ELE1 antibody was raised against a peptide derived from the C-terminal region of human NCoA4.

**Buffer:** Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

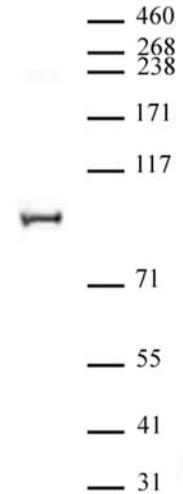
**Application Notes:**

Applications Validated by Active Motif:

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



**NCoA4 / ELE1 antibody (pAb) tested by Western blot.**

Nuclear extract (20 µg) of LNCaP cells probed with NCoA4 / ELE1 antibody at a dilution of 1:2,000.