

## PELP1 antibody (pAb)

**Catalog Nos:** 61263, 61264

**RRID:** AB\_2615014

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 150 kDa

**Background:** The PELP1 gene (proline, glutamic acid and leucine rich protein 1, Modulator of non-genomic activity of estrogen receptor, HMX3, MNAR, P160) encodes a protein that is a coactivator of estrogen receptor-mediated transcription and a corepressor of other nuclear hormone receptors and sequence-specific transcription factors. Plays a role in estrogen receptor (ER) genomic activity when present in the nuclear compartment by activating the ER target genes in a hormonal stimulation dependent manner. PELP1 interacts with KDM1 (LSD1) and may act as reader of histone methyl modifications for ER mediated gene activation.

**Immunogen:** This PELP1 antibody was raised against a peptide within the internal region of human PELP1.

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

### Application Notes:

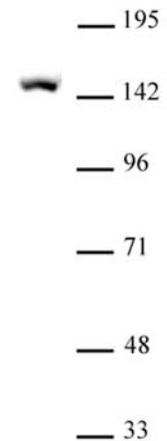
Applications Validated by Active Motif:

WB: 1:500 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. 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.



### PELP1 antibody (pAb) tested by Western blot

Whole cell extract of MCF-7 cells (30  $\mu$ g) probed with PELP1 antibody (1:500).