

## PXR antibody (pAb)

**Catalog Nos:** 39589, 39590

**RRID:** AB\_2793269

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 200 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 54 kDa

**Background:** PXR (NR1I2, Pregnane X receptor, nuclear receptor subfamily 1 group I member 2) belongs to a family of transcription factors characterized by a ligand-binding domain and a DNA-binding domain. PXR is an orphan receptor that regulates transcription of the cytochrome P450 gene CYP3A4, binding to the response element of the CYP3A4 promoter as a heterodimer with the 9-cis retinoic acid receptor RXR (RXR- $\alpha$  and RXR- $\beta$ ). PXR is activated by a range of compounds that induce CYP3A4, including dexamethasone and rifampicin.

**Immunogen:** This PXR antibody was raised against a peptide containing amino acid residues 57-71 of human PXR.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

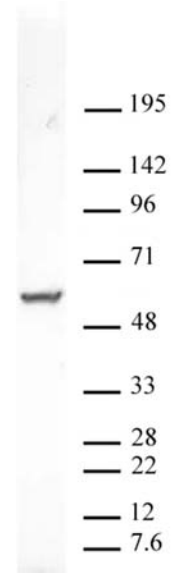
### Application Notes:

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.



### PXR pAb tested by Western blot.

Hep G2 nuclear extract (20 µg per lane) was probed with PXR pAb at a dilution of 1:1,000.