

## LXR- $\alpha$ antibody (pAb)

---

**Catalog Nos:** 61175, 61176

**RRID:** AB\_2614981

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

**Reactivity:** Human, Mouse

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

**Purification:** Affinity Purified

**Host:** Rabbit

**Isotype:** IgG

**Molecular Weight:** 52 kDa

---

**Background:** LXR- $\alpha$  (Liver X receptor alpha, NR1H3) is an orphan nuclear receptor and a member of the NR1 receptor family. The LXR receptors (including LXR- $\alpha$  and LXR- $\beta$ ) bind oxysterols and play key roles in maintaining cholesterol homeostasis in macrophages, primarily by regulating multiple components of the reverse cholesterol transport. They also are potent inhibitors of inflammation and are capable of repressing cytokine and chemokine production by Toll-like receptor (TLR)-activated macrophages. LXR- $\alpha$  is expressed mainly in liver and adipose tissue and interacts with RXR, shifting RXR to an active ligand-binding subunit mediating responses to retinoids.

**Immunogen:** This LXR- $\alpha$  antibody was raised against a peptide in the N-terminal region of human LXR- $\alpha$ .

**Buffer:** Purified rabbit IgG in 70 mM Tris, 105 mM NaCl, 0.07 mM EDTA, 31 mM Glycine, 0.035% Sodium Azide, and 30% glycerol.

**Application Notes:**

Applications Validated by Active Motif:

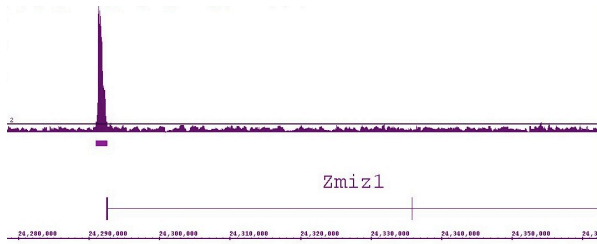
ChIP: 10  $\mu$ l per ChIP

ChIP-Seq: 10  $\mu$ l each

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

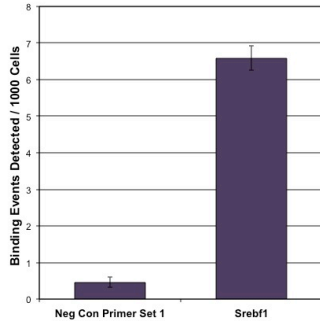
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



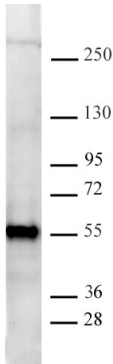
**LXR-α antibody (pAb) tested by ChIP-chip.**

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with chromatin from 2 million mouse embryonic fibroblasts. ChIP DNA was amplified by WGA, labeled and hybridized to a mouse tiling array. The image is zoomed in to show LXR-α binding at the promoter of the retinoic acid inducible gene Zmiz1. LXR binding at this gene is expected since LXR is known to form heterodimers with RXR (retinoic acid receptor).



**LXR-α antibody (pAb) tested by ChIP.**

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 20 µg of mouse embryonic fibroblast chromatin and 10 µl of LXR-α 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.



**LXR-α antibody (pAb) tested by Western blot.**

Whole cell extract of Hep G2 cells (30 µg) probed with LXR-α antibody (pAb) at a dilution of 1:500.