

LXR- β antibody (pAb)

Catalog Nos: 61177, 61178

RRID: AB_2614980

Isotype: IgG

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

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.5 μ g/ μ l

Molecular Weight: 55 kDa

Background: LXR- β (Liver X receptor beta, NR1H2, NER, UNR) is an orphan nuclear receptor and a member of the NR1 receptor family. The LXR receptors (including LXR- α and LXR- β) 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- β is more widely expressed than LXR- α , particularly in the brain.

Immunogen: This LXR- β antibody was raised against a peptide in the N-terminal region of human LXR- β .

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

Application Notes:

Applications Validated by Active Motif:

ChIP: 10 μ l per ChIP

ChIP-Seq: 10 μ l each

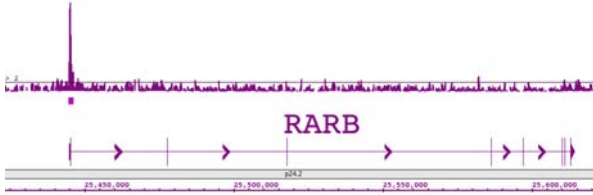
WB: 1:500 - 1:1,000 dilution

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.

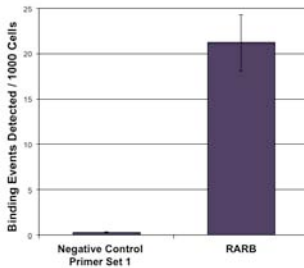
LXR-β pAb tested by ChIP-Chip.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with chromatin from 2.5 million primary human liver cells. ChIP DNA was amplified by WGA, labeled and hybridized to a human tiling array. The image is zoomed in to show LXR-β binding at a known LXR binding site in the RARB promoter.



LXR-β pAb tested by ChIP-qPCR.

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 25 μg of human liver chromatin and 10 μl of LXR-β antibody. ChIP DNA was used in qPCR with the negative control primer pairs and primers against the known LXR binding site in the RARB promoter. 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-β pAb tested by Western blot.

Detection of LXR-β by Western blot analysis. HeLa whole-cell extract (20 μg) was probed with LXR-β pAb at a 1:1000 dilution.

