

RXR- α antibody (mAb)

Catalog No: 61029

RRID: AB_2793479

Clone: 4RX-3A2

Isotype: IgG1

Application(s): EMSA, IP, WB

Reactivity: Human

Quantity: 100 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 0.49 μ g/ μ l

Molecular Weight: 52 kDa

Background: RXR- α (Retinoid X Receptor alpha, NR2B1) is a nuclear receptor that mediates the biological effects of retinoids by its involvement in retinoic acid-mediated gene activation. RXR- α exerts its action by binding, as a homodimer or a heterodimer, to specific sequences in the promoters of target genes and regulating their transcription. RXR- α is a member of the steroid and thyroid hormone receptor superfamily of transcriptional regulators.

Immunogen: This RXR-alpha antibody was raised against a recombinant protein corresponding to the C-terminal region of mouse RXR-alpha.

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:

Published Applications:

IP

EMSA

WB

See references for more information. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.