

RXR- β antibody (mAb)

Catalog No: 61031

RRID: AB_2793480

Clone: 16RX-2E1

Isotype: IgG1

Application(s): IHC, WB

Reactivity: Human

Quantity: 100 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 56 kDa

Background: RXR- β (retinoid X receptor beta, NR2B2) is a member of the nuclear receptors superfamily, specifically those involved in mediating the effects of retinoic acid (RA). RXR- β forms heterodimers with the RAR (retinoic acid receptor), TR (thyroid hormone receptor) and the VDR (vitamin D receptor), increasing both DNA binding and transcriptional function on their respective response elements.

Immunogen: This RXR-beta antibody was raised against a peptide within the N-terminal A region of mouse RXR-beta.

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:

IHC

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.