

RXR- γ antibody (mAb)

Catalog No: 61033

RRID: AB_2793481

Clone: 12RX-2D3

Isotype: IgG

Application(s): EMSA, IHC, WB

Reactivity: Human

Quantity: 100 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 51 kDa

Background: RXR- γ (retinoid X receptor gamma, NR2B3) is a member of the nuclear receptors superfamily, specifically those involved in mediating the effects of retinoic acid (RA). This receptor 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-gamma antibody was raised against a peptide within the D-region of mouse RXR-gamma.

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:

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.