

RAR- β antibody (mAb)

Catalog Nos: 61027, 61028

RRID: AB_2793478

Clone: 8B-10B2

Isotype: IgG1

Application(s): EMSA, IHC, WB

Reactivity: Human

Quantities: 100 μ g, 10 μ g

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 μ g/ μ l

Molecular Weight: 50 kDa

Background: RAR- β (HAP, NR1B2) is a member of the thyroid-steroid hormone receptor superfamily of nuclear transcriptional regulators. RAR- β localizes to the cytoplasm and to subnuclear compartments. RAR- β binds retinoic acid, the biologically active form of vitamin A which mediates cellular signaling in embryonic morphogenesis, cell growth and differentiation. It is thought that RAR- β limits growth of many cell types by regulating gene expression.

Immunogen: This RAR-beta antibody was raised against amino acids 413-448 of human RAR-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:

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.