

## RAR- $\gamma$ antibody (mAb)

---

**Catalog Nos:** 39973, 39974

**RRID:** AB\_2614944

**Clone:** 4 $\gamma$ -7A11

**Isotype:** IgG1, k

**Application(s):** EMSA, IHC, WB

**Reactivity:** Human

**Quantities:** 100  $\mu$ g, 10  $\mu$ g

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1  $\mu$ g/ $\mu$ l

**Molecular Weight:** 52 kDa

---

**Background:** Retinoic acid receptors (RAR) are nuclear receptors that are activated by both all-trans retinoic acid and 9-cis retinoic acid. There are three retinoic acid receptors (RAR), RAR- $\alpha$ , RAR- $\beta$  and RAR- $\gamma$ . RAR- $\gamma$  has two splice isoforms. RAR- $\gamma$  heterodimerizes with RXR and in the absence of ligand, the RAR/RXR dimer binds to hormone response elements complexed with co-repressor protein. Binding of agonist ligands to RAR results in dissociation of co-repressor and recruitment of coactivator protein which in turn promotes transcription of the downstream target gene into mRNA and eventually protein.

**Immunogen:** This RAR- $\gamma$  antibody was raised against a recombinant protein corresponding to amino acids 428-441 (F-region) of human RAR $\gamma$ .

**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.