

RAR- α antibody (mAb)

Catalog No: 39971

RRID: AB_2793414

Clone: 9 α -9A6

Isotype: IgG1, k

Application(s): EMSA, IHC, WB

Reactivity: Human, Mouse

Quantity: 100 μ g

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 μ g/ μ 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- α , RAR- β and RAR- γ . RAR- α (Nuclear receptor subfamily 1 group B member 1, NR1B1) interacts with NCoA3, NCoA6, NCoA7 and MLL to activate transcription of target genes. Chromosomal aberrations of RAR- α are sometimes involved in Promyelocytic leukemia.

Immunogen: This RAR- α antibody was raised against a recombinant protein corresponding to amino acids 420-443 (the F region) of human RAR α .

Buffer: Purified IgG in 60 mM Sodium Citrate, 105 mM Tris-HCl, 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.