

Recombinant RAR- α protein

Catalog No: 31130

Expressed In: Baculovirus

Quantity: 5 μ g

Concentration: 0.5 μ g/ μ l

Source: Human

Buffer Contents: 5 μ g of Recombinant RAR- α protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 8), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

Protein Details: Recombinant RAR- α protein was expressed in baculovirus (accession number NM 000964) with an N-terminal His-Tag and purified by an affinity column in combination with FPLC chromatography. The purified recombinant protein is greater than 90% homogeneous and contains no detectable protease, DNase and RNase activity.

Application Notes: Recombinant RAR- α is suitable for DNA-protein, *in vitro* transcription and protein-protein interaction assays. 20 ng is sufficient for DNA-protein assays, 20-100 ng is sufficient for *in vitro* transcription assays and 100 ng is sufficient for protein-protein interaction studies. The molecular weight of the protein is ~55 kDa.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. 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.