

Recombinant LXR β protein

Catalog No: 31123

Expressed In: *E. coli*

Quantity: 10 μ g

Concentration: 0.5 μ g/ μ l

Source: Human

Buffer Contents: 10 μ g of Recombinant LXR β 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 LXR β is isolated from an *E. coli* strain that carries the coding sequence of the human LXR β under the control of a T7 promoter (accession number NM 007121). The purified recombinant protein has an N-terminal His-Tag and is greater than 95% homogeneous and contains no detectable protease, DNase and RNase activity.

Application Notes: Recombinant LXR β is suitable for DNA and protein-protein interaction assays. 20 ng is sufficient for gelshift assays and 100 ng is sufficient for protein-protein interaction studies. The molecular weight of the protein is ~62 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.