

## Recombinant LXR $\alpha$ protein

---

**Catalog No:** 31122

**Expressed In:** *E. coli*

**Quantity:** 10  $\mu$ g

**Concentration:** 0.2  $\mu$ g/ $\mu$ l

**Source:** Human

**Buffer Contents:** 10  $\mu$ g of Recombinant LXR $\alpha$  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 $\alpha$  is isolated from an *E. coli* strain that carries the coding sequence of the human LXR $\alpha$  under the control of a T7 promoter (accession number NM 005693). 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 $\alpha$  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 ~56 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.