

Recombinant Estrogen Receptor α protein

Catalog No: 81276, 81976

Expressed In: Baculovirus

Quantity: 20, 1000 μ g

Concentration: 0.2 μ g/ μ l

Source: Human

Buffer Contents: Recombinant Estrogen Receptor-alpha protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

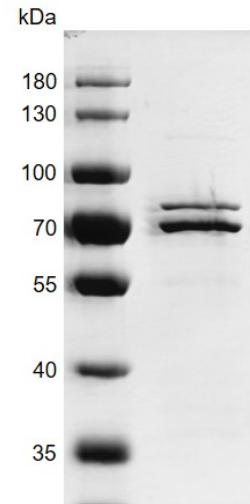
Background: Estrogen Receptor-alpha protein, also called as ESRA or ER-alpha, is a nuclear hormone receptor. The steroid hormones and their receptors are involved in the regulation of eukaryotic gene expression and affect cellular proliferation and differentiation in target tissues. Ligand-dependent nuclear transactivation involves either direct homodimer binding to a palindromic estrogen response element (ERE) sequence or association with other DNA-binding transcription factors, such as AP-1/c-Jun, c-Fos, ATF-2, Sp1 and Sp3, to mediate ERE-independent signaling. Ligand binding induces a conformational change allowing subsequent or combinatorial association with multiprotein coactivator complexes through LXXLL motifs of their respective components. Mutual transrepression occurs between the estrogen receptor (ER) and NF-kappa-B in a cell-type specific manner. ER-alpha can also act synergistically with NF-kappa-B to activate transcription involving respective recruitment adjacent response elements. It also mediates membrane-initiated estrogen signaling involving various kinase cascades.

Protein Details: Recombinant Estrogen Receptor-alpha protein was expressed in baculovirus system as the full length protein (accession number NP_000116.2) with an N-terminal FLAG tag. The molecular weight of the protein is 67.5 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

Estrogen Receptor-alpha



Recombinant Estrogen Receptor-alpha protein gel

9% SDS-PAGE gel with Coomassie blue staining

MW: 67.5 kDa

Purity: >80%