

Recombinant SSRP1 / FACT p80 protein

Catalog No: 81094, 81894

Lot No: 32217001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: Recombinant SSRP1 / FACT p80 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.

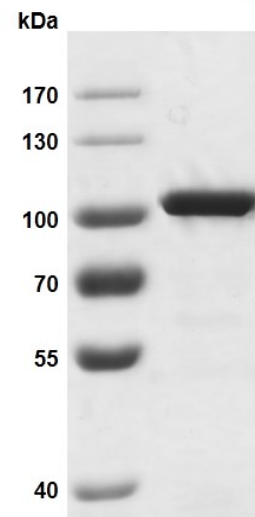
Background: SSRP1 (Structure Specific Recognition Protein 1), also known as FACT p80 (Facilitates Chromatin Transcription Complex 80 KDa Subunit), is a component of the FACT complex, a general chromatin factor that acts to reorganize nucleosomes. The FACT complex is involved in multiple processes that require DNA as a template such as mRNA elongation, DNA replication and DNA repair. During transcription elongation the FACT complex acts as a histone chaperone that both destabilizes and restores nucleosomal structure. It facilitates the passage of RNA polymerase II and transcription by promoting the dissociation of one histone H2A-H2B dimer from the nucleosome, then subsequently promotes the reestablishment of the nucleosome following the passage of RNA polymerase II. The FACT complex is probably also involved in phosphorylation of Ser-392 of p53/TP53 via its association with CK2 (casein kinase II).

Protein Details: Recombinant human SSRP1 / FACT p80 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_003137.1) with an N-terminal FLAG tag. The molecular weight of the protein is 82.2 kDa.

Application Notes: This protein is suitable for use in protein-protein interaction, *in vitro* transcription assay, binding assay.

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.

SSRP1 / FACT p80



Recombinant SSRP1 / FACT p80 protein gel

7.5% SDS-PAGE Coomassie staining
MW: 82.2 kDa
Purity: >95%