

Recombinant Spt16 / FACT p140 protein

Catalog No: 31344

Source: Human

Quantity: 10 µg

Expressed In: Sf9 cells

Buffer: 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 µg recombinant Spt16 / FACT p140 protein, 0.5 µg/µl.

Protein Details:

Spt16 (p140) is a component of FACT. Spt16 was expressed in Sf9 cells (accession number Q9Y5B9) with an N-terminal FLAG tag. The molecular weight of the protein is ~140 kDa.

Dilution of Protein:

Recombinant Spt16 is supplied in solution and should be kept as a solution in order to maintain full activity.

Application:

Recombinant Spt16 (p140) protein is suitable for use in in vitro transcription assays. We recommend to start with a range of protein concentrations (e.g. 5 ng, 50 ng, and 500 ng) to determine the optimal amount needed for each in vitro transcription assay.

Storage:

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

Guarantee:

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.