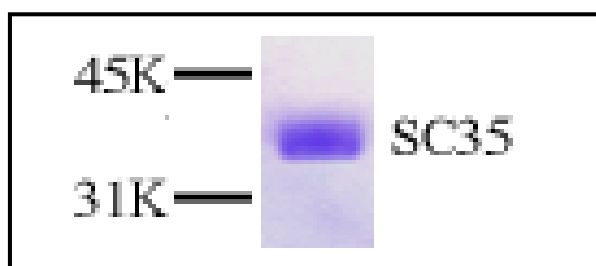


SC35 (SFRS2 or SRp30b)

Splicing Component, 35 kDa, Pre-mRNA Splicing Factor

human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-776	5 µg



For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 100 mM KCl, 0.2 mM EDTA, 1 mM DTT, 20% glycerol.

Activity

1 ng is the amount sufficient for a gel mobility shift assay in a 20 µl reaction, 50 and 100 ng are sufficient for a protein-protein interaction assay.

Application

Recombinant SC35 protein can be used 1) for *in vitro* function studies including pre-mRNA splicing, cross linking and other RNA binding assays, 2) for protein-protein interaction assay, and 3) for cell growth and proliferation assays.

Molecular Weight

35 kDa

Purity

> 95% by SDS-PAGE

Description

SC35, a member of SR protein family, is an essential pre-mRNA splicing factor. Phosphorylation on serine residues located within the SR domain directly regulates SC35 activity and compartmentalization of other SR splicing factors. In addition to interacting with RNA and other splicing factors, SC35 has been shown to interact either directly or indirectly with the C-terminal domain (CTD) of the largest subunit of RNA polymerase II, thereby suggesting a potential role of SC35 in coordinating of transcription and pre-mRNA splicing.

The human SC35 wild type protein was expressed in baculovirus system and purified by an affinity column in combination with FPLC chromatography.

Selected References:

- Fu *et al.* (1992) Isolation of a complementary DNA that encodes the mammalian splicing factor SC35. *Science* **256**:535.
Zahler *et al.* (1992) SR proteins: a conserved family of pre-mRNA splicing factors. *Genes & Dev.* **6**:837.
Gui *et al.* (1994) Purification and characterization of a kinase specific for the serine- and arginine-rich pre-mRNA splicing factors. *Proc. Natl. Acad. Sci. USA* **91**:10824.
Colwill *et al.* (1996) The Clk/Sty protein kinase phosphorylates SR splicing factors and regulates their intranuclear distribution. *EMBO J.* **15**:265.
Misteli *et al.* (1997) The dynamics of a pre-mRNA splicing factor in living cells. *Nature* **387**:523.