

PRPF4B Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P87-35G

Lot # A1266-4

Product Description

Recombinant human PRPF4B (551-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC034969](#).

Gene Aliases

PRP4, PRP4H, PRP4K, PR4H, KIAA0536, dJ1013A10.1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

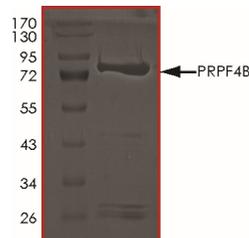
Scientific Background

PRPF4B or pre-mRNA processing factor 4B belongs to a kinase family that includes serine/arginine-rich protein-specific kinases and cyclin-dependent kinases (CDKs) which is regarded as a CDK-like kinase (Clk) with homology to mitogen-activated protein kinases (MAPKs). PRPF4B is involved in pre-mRNA splicing and isolation of a putative mammalian homologue (1).

References

- Gross, T.et.al: Functional analysis of the fission yeast Prp4 protein kinase involved in pre-mRNA splicing and isolation of a putative mammalian homologue. Nucleic Acids Res. 25: 1028-1035, 1997.

Purity



The purity of PRPF4B protein was determined to be **>90%** by densitometry. Approx. MW **82 kDa**.

PRPF4B Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P87-35G
Lot #	A1266-4
Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.