

PRPF19 Protein

Recombinant protein expressed in insect cells

Catalog # P270-30G

Lot # O988-2

Product Description

Recombinant human PRPF19 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_014502](#).

Gene Aliases

PSO4; SNEV; PRP19; UBOX4; hPSO4; NMP200

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.

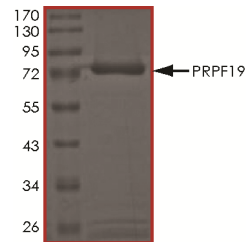
Scientific Background

PRPF19 or pre-mRNA-processing factor 19 is the human homolog of yeast Pso4. PRPF19 is a gene essential for cell survival, DNA repair and association with terminal deoxynucleotidyl transferase (1). PRPF19 forms a stable complex with SETMAR (Metnase) which is a SET and transposase fusion protein in humans and plays a positive role in double-strand break (DSB) repair. PRPF19 mediates association of SETMAR with sites of DNA damage (2). SNEV is an orthologue of PRPF19 and is an evolutionarily conserved splicing factor whose oligomerization is necessary for spliceosome assembly.

References

1. Mahajan, K. N. et.al: Role of human Pso4 in mammalian DNA repair and association with terminal deoxynucleotidyl transferase. Proc. Nat. Acad. Sci. 100: 10746-10751, 2003.
2. Beck, B. D. et.al: Human Pso4 is a Metnase (SETMAR)-binding partner that regulates Metnase function in DNA repair. J. Biol. Chem. 283: 9023-9030, 2008.

Purity



The purity of PRPF19 protein was determined to be **>90%** by densitometry. Approx. MW **74 kDa**.

PRPF19 Protein

Recombinant protein expressed in insect cells

Catalog Number	P270-30G
Specific Lot Number	O988-2
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.