

PPP2R1A Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # P162-30G

Lot # U1692-6

Product Description

Recombinant full-length human PPP2R1A was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PPP2R1A gene accession number is [NM_014225](#).

Gene Aliases

MRD36; PP2A-Aalpha; PP2AAALPHA; PR65A

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

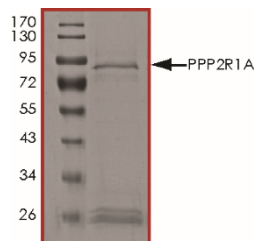
Scientific Background

PPP2R1A or the PR65 subunit of protein phosphatase 2A (PP2A) acts as a scaffolding protein to coordinate the assembly of the catalytic subunit and a variable regulatory B subunit, together forming the core of heterodimeric PP2A. Required for proper chromosome segregation and for centromeric localization of SGOL1 in mitosis.

References

1. Tang Z, et al: PP2A is required for centromeric localization of Sgo1 and proper chromosome segregation. Dev Cell. 10(5):575-85, 2006.
2. <http://www.uniprot.org/uniprot/P30153>.

Purity



The purity of PPP2R1A was determined to be **>70%** by densitometry, approx. MW **92 kDa**.

PPP2R1A Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	P162-30G
Lot #	U1692-6
Purity	>70%
Concentration	0.05 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.