

PP2B Protein

Recombinant full-length protein expressed in Sf9 cells

Catalog # P18-34G

Lot # S334-5

Product Description

Recombinant full-length human PP2B was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PP2B gene accession number is [NM_001130691](#).

Gene Aliases

PPP2B, PPP3CA, CALN, CCN1, CNA1, CALNA, CALNA1

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.

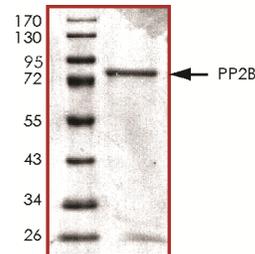
Scientific Background

PP2B is a key regulator of the anaphase-promoting factor which is an excellent substrate for protein phosphatase. PP2B is involved in bovine ocular tissues and is postulated to be involved in the immunologic privilege of the cornea, in retinal signal transduction, and in the toxic effects of immunosuppressants on the eye. PP2B acts as an inhibitory constraint on L-LTP that is relieved by PKA, and that this inhibitory constraint acts as a gate to regulate the synaptic induction of L-LTP (1). PP2B plays a role in the transition from short- to long-term memory and also regulates bone formation through an effect on osteoblast differentiation (2).

References

1. Winder. et.al: Genetic and pharmacological evidence for a novel, intermediate phase of long-term potentiation suppressed by calcineurin. *Cell* 92: 25-37, 1998.
2. Sun, L. et.al: Calcineurin regulates bone formation by the osteoblast. *Proc. Nat. Acad. Sci.* 102: 17130-17135, 2005.

Purity



The purity of PP2B was determined to be **>85%** by densitometry. Approx. MW **83 kDa**.

PP2B Protein

Recombinant full-length protein expressed in Sf9 cells

Catalog Number	P18-34G
Specific Lot Number	S334-5
Purity	>85%
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.