

CDK5P25 Protein

Recombinant protein expressed in Sf9 cells

Catalog # C34-31G

Lot # N092-3

Product Description

Recombinant human CDK5P25 (CDK5R1) (98-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_003885](#).

Gene Aliases

p25, CDK5R1; CDK5R, NCK5A, p23

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

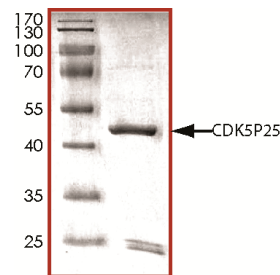
CDK5P25 (cyclin-dependent kinase 5, regulatory subunit 1) is a neuron-specific activator of CDK5 and is required for proper development of the central nervous system (1). The p35 form of this protein is proteolytically cleaved by calpain, generating a p25 form of the protein which relocates from the cell periphery to nuclear and perinuclear regions. CDK5P25 regulates CDK5 activity by prolonging its activation and changing its cellular location and this may be important in pain signaling (2). The CDK5P25 accumulates in the brain neurons of patients with Alzheimer's disease and this accumulation correlates with an increase in CDK5 kinase activity which may result in aberrantly phosphorylated forms of the microtubule-associated protein tau.

References

1. Tsai, L.-H.et.al: p35 is a neural-specific regulatory subunit of cyclin-dependent kinase 5. Nature 371: 419-423, 1994.
2. Pareek, T. K.et.al: Cyclin-dependent kinase 5 activity regulates pain signaling. Proc. Nat. Acad. Sci. 103: 791-796, 2006.

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

Purity



The purity of CDK5P25 was determined to be **>85%** by densitometry. Approx. MW **49 kDa**.

CDK5P25 Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	C34-31G
Specific Lot Number	N092-3
Purity	>85%
Concentration	0.2 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.