

## CDK5P35 Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # C34-30G**

Lot # A1435-2

### Product Description

Recombinant full-length human CDK5P35 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_003885](#).

### Gene Aliases

p35, CDK5R1, CDK5R, NCK5A, p35nck5a

### 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  $-70^{\circ}\text{C}$ . 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

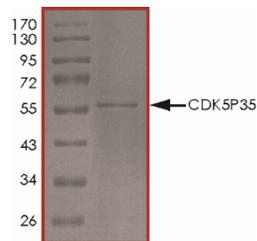
CDK5P35 is a neuron-specific activator of cyclin-dependent kinase 5 (CDK5) that bind to cyclin to form active holoenzymes that play a pivotal role in the regulation of the eukaryotic cell cycle. The activation of CDK5 is required for proper development of the central nervous system. p35 is a regulatory subunit for CDK5 which is proteolytically cleaved by calpain and generating a p25 form that involved in the pathogenesis of cytoskeletal abnormalities and neuronal death in neurodegenerative diseases (1). Cdk/p35 plays a role in primary afferent nociceptive signaling and is required for cortical plate neurons to migrate past preexisting neurons and take up superficial positions to constitute the inside-outside layering order of cortical lamination(2).

### References

1. Patrick, G. N.et.al: Conversion of p35 to p25 deregulates Cdk5 activity and promotes neurodegeneration. Nature 402: 615-622, 1999.
2. Chae, T. et.al: Mice lacking p35, a neuronal specific activator of Cdk5, display cortical lamination defects, seizures, and adult lethality. Neuron 18: 29-42, 1997.

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

### Purity



The purity was determined to be **>80%** by densitometry. Approx. MW **60 kDa**.

## CDK5P35 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	C34-30G
Lot #	A1435-2
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.