

STK9 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # C46-31G

Lot # Q2488-5

Product Description

Recombinant human STK9 (CDKL5) (1-498) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_003159](#).

Gene Aliases

CDKL5; EIEE2; ISSX

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.

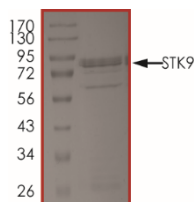
Scientific Background

STK9 or CDKL5 is a member of Ser/Thr protein kinase family which has protein kinase activity and is associated with X-linked infantile spasm syndrome (ISSX), also known as X-linked West syndrome, and Rett syndrome (RTT). STK9 can control nuclear speckle morphology by regulating the phosphorylation state of splicing regulatory proteins, and may be involved indirectly in pre-mRNA processing, by controlling splicing factor dynamics (1). The mutations in the STK9 is a major cause of infantile spasms and early epileptic seizures in female patients, and of a later intractable seizure disorder (2).

References

1. Ricciardi, S. et.al: CDKL5 influences RNA splicing activity by its association to the nuclear speckle molecular machinery. Hum. Molec. Genet. 18: 4590-4602, 2009.
2. Archer, H. L. et.al: CDKL5 mutations cause infantile spasms, early onset seizures, and severe mental retardation in female patients. J. Med. Genet. 43: 729-734, 2006.

Purity



The purity of STK9 was determined to be **>85%** by densitometry, approx. MW **84 kDa**.

STK9 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	C46-31G
Lot #	Q2488-5
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.