

ULK4 Protein

Recombinant protein expressed in Sf9 cells

Catalog # U04-35G

Lot # 1027-2

Product Description

Recombinant human ULK4 (1-616) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_017886](#).

Gene Aliases

DKFZp434E1822; FAM7C1; FLJ20574; REC01035

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.

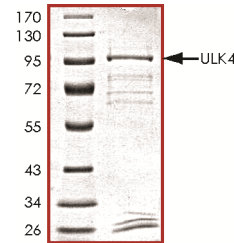
Scientific Background

ULK4, or Unc-51-like kinase 4, is a Ser/Thr protein kinase that is a member of the APG1/unc-51/ULK1 subfamily. ULK4 contains five HEAT repeats and a catalytic kinase domain. The ULK4 gene is conserved in many species including chimpanzee, dog, cow, mouse, rat, chicken, zebrafish, *A.thaliana*, and rice (1). ULK4 has been implicated as a gene involved in the development of hydrocephalus (2). This neurologic disorder in animal models is being used to elucidate factors responsible for the excessive accumulation of cerebrospinal fluid in hydrocephalic humans.

References

1. Krupnova T et al: Microtubule-associated kinase-like protein RUNKEL needed [corrected] for cell plate expansion in Arabidopsis cytokinesis. *Curr Biol.* 2009 Mar 24;19(6):518-23.
2. Vogel, P. et al: Congenital Hydrocephalus in Genetically Engineered Mice. *Vet Pathol.* 2011 Jul 11.

Purity



The purity was determined to be **>75%** by densitometry. Approx. MW **97 kDa**.

ULK4 Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	U04-35G
Specific Lot Number	1027-2
Purity	>75%
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.