

BCKDK Protein

Recombinant protein expressed in Sf9 cells

Catalog # B15-31G

Lot # O845-3

Product Description

Recombinant human BCKDK (30-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The BCKDK gene accession number is [NM_005881](#).

Gene Aliases

BCKDKD; BDK

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.

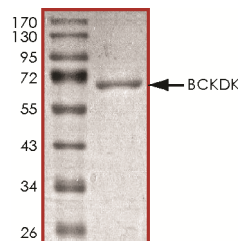
Scientific Background

BCKDK or branched chain alpha-keto acid dehydrogenase kinase is an important regulator of the valine, leucine, and isoleucine catabolic pathways. BCKDK is found in the mitochondrion where it phosphorylates and inactivates BCKD. The branched-chain alpha-keto dehydrogenase complex (BCKDH) catalyzes the rate-limiting step in the oxidation of branched-chain amino acids and BCKDK inactivates the BCKDH complex by catalyzing serine phosphorylation of the E1-alpha subunit (1). BCKDH kinase catalyzes the serine phosphorylation of the E1- alpha subunit of BCKDH (2).

References

1. Joshi, M. A. et.al: growth and neurological abnormalities in branched-chain alpha-keto acid dehydrogenase kinase-deficient mice. *Biochem. J.* 400: 153-162, 2006.
2. Popov, K. M. et.al: A. Branched-chain alpha-ketoacid dehydrogenase kinase: molecular cloning, expression, and sequence similarity with histidine protein kinases. *J. Biol. Chem.* 267: 13127-13130, 1992.

Purity



The purity of BCKDK was determined to be **>95%** by densitometry, approx. MW **69kDa**.

BCKDK Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	B15-31G
Lot #	O845-3
Purity	>95%
Concentration	0.1 µg/µl
Stability	1 yr 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.