

LKB1(STK11) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # L15-30G

Lot # H2668-9

Product Description

Recombinant full-length human LKB1 (STK11) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_000455](#).

Gene Aliases

STK11, LKB1, PJS

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.

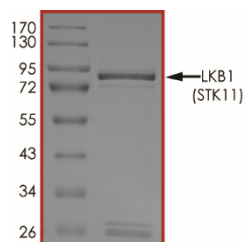
Scientific Background

LKB1 (STK11) is a member of the serine/threonine kinase family which regulates cell polarity and functions as a tumor suppressor. STK11 is a critical barrier to pulmonary tumorigenesis which controlling initiation, differentiation, and metastasis. STK11 plays a central role in restricting HSC entry into cell cycle and in broadly maintaining energy homeostasis in hematopoietic cells through a novel metabolic checkpoint (1). STK11 serves as an essential regulator of HSCs and hematopoiesis, and more generally, points to the critical importance of coupling energy metabolism and stem cell homeostasis (2).

References

1. Gurumurthy, S. et.al: The Lkb1 metabolic sensor maintains haematopoietic stem cell survival. Nature 468: 659-663, 2010. Note: Erratum: Nature 476: 114 only, 2011.
2. Gan, B.et.al: Lkb1 regulates quiescence and metabolic homeostasis of haematopoietic stem cells. Nature 468: 701-704, 2010.

Purity



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

Upstream Active Kinases

SRC, Active
MEK1, Active

Cat # S19-10G
Cat # M02-13G

LKB1 (STK11) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	L15-30G
Lot #	H2668-9
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.

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.