

LATS2, Unactive

Recombinant protein expressed in Sf9 cells

Catalog # L02-35G

Lot # C104-2

Product Description

Recombinant human LATS2 (480-1088) was expressed by baculovirus in Sf9 cells using an N-terminal GST tag. The gene accession number is [NM_014572](#).

Gene Aliases

KPM, FLJ13161

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

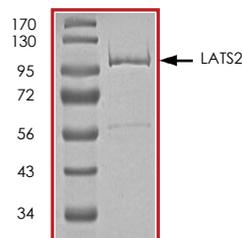
LATS2 is a serine/threonine protein kinase belonging to the LATS tumor suppressor family (1). The kinase activity and two conserved domains within LATS2 are responsible for the suppression of tumorigenicity and inhibition of cell growth. LATS2 negatively regulates the cell cycle by controlling G1/S and/or G2/M transition. LATS2 interacts with the centrosomal proteins AURORA A and Ajuba and is required for accumulation of γ -tubulin and spindle formation at the onset of mitosis (2). LATS2 also interacts with a negative regulator of p53 and may function in a positive feedback loop with p53 that responds to cytoskeleton damage. LATS2 can also function as a co-repressor of androgen-responsive gene expression.

References

- Li, Y. et al: Lats2, a putative tumor suppressor, inhibits G1/S transition. *Oncogene*. 2003; 22(28):4398-405.
- Yabuta, N. et al: Structure, expression, and chromosome mapping of LATS2, a mammalian homologue of the Drosophila tumor suppressor gene lats/warts. *Genomics*. 2000 63(2):263-70.

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

Purity



The purity of LATS2 was determined to be **>80%** by densitometry. Approx. MW **98kDa**.

LATS2, Unactive

Recombinant protein expressed in Sf9 cells

Catalog Number	L02-35G
Specific Lot Number	C104-2
Purity	>80%
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.