

Recombinant STK10 / LOK (1-348) protein

Catalog No: 81378, 81678

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.3 µg/µl

Source: Human

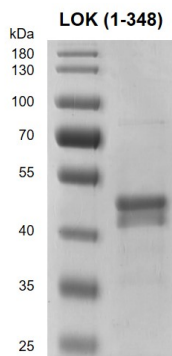
Buffer Contents: Recombinant STK10 / LOK (1-348) protein is supplied in 25 mM HEPESNaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: STK10 (Serine/Threonine Kinase 10) is a serine/threonine-protein kinase involved in regulation of lymphocyte migration. STK10 phosphorylates MSN and involved in regulation of lymphocyte migration by mediating phosphorylation of ERM proteins such as MSN. It acts as a negative regulator of MAP3K1/MEKK1 and may also act as a cell cycle regulator by acting as a polo kinase kinase. Diseases associated with STK10 include Testicular Germ Cell Tumor.

Protein Details: Recombinant STK10 / LOK (1-348) protein that includes amino acids 1-348 of human LOK protein (accession number NP_005981.3) was expressed in a baculovirus expression system, and contains an N-terminal FLAG tag. The molecular weight of the protein is 40.4 kDa

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.



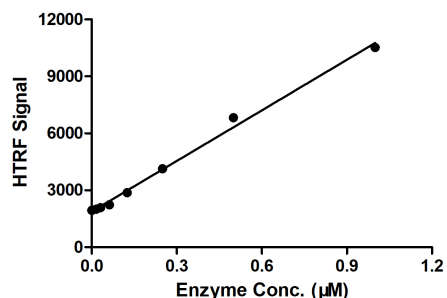
Recombinant STK10 / LOK (1-348) protein

10% SDS-PAGE with Coomassie staining

MW: 40.4

kDa Purity: >80%

STK10 / LOK (1-348) Titration



HTRF assay for STK10 / LOK (1-348) activity

1 μM STK S3 substrate was incubated with different concentrations of STK10 / LOK (1-348) protein in a 10 μl reaction system containing 1×Enzymatic Buffer, 5 mM MgCl₂, 1 mM DTT and 100 μM ATP for 1 hour. Then 10 μl detection reagents containing STK antibody (1:2) and SA-XL665 (1:100) diluted with 1× Detection Buffer were added and incubated with the reactions for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.