

Anti-phospho-STK10-α (Ser482 Ser486 Ser490)

Rabbit Polyclonal Antibody

Catalog # S29-165R

Lot # J1274-31

Cited Applications

WB

Suggested Dilutions:

WB 1:1,000

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the STK10-α protein phosphorylated at serine 482/486/490

Cross Reactivity

Human, Mouse, Rat and Zebrafish

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser482/486/490 conjugated to KLH

Formulation

100 µl in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

Scientific Background

Polo-like kinase (PLK) plays an important role in cell cycle regulation. PLK is also required for bipolar spindle formation, activation of the anaphase-promoting complex/cyclosome, and cytokinesis. Deregulation of PLKs often results in abnormalities in centrosome duplication, maturation and/or microtubule dynamics (1). STK10-α, which is also known as PLK1 is likely responsible for activating PLK. STK10-α itself is activated by phosphorylation at Ser482, Ser486 and Ser490 (2).

References

1. Nigg, EA. et al: Dynamic changes in nuclear architecture during mitosis: on the role of protein phosphorylation in spindle assembly and chromosome segregation. *Exp Cell Res.* 1996 Dec 15;229.
2. Erikson E et al. A feedback loop in the polo-like kinase activation pathway. *J Biol Chem.* 2004, 279:32219-32224.

Anti-phospho-STK10-α (Ser482 Ser486 Ser490)

Rabbit Polyclonal Antibody

Catalog Number

S29-165R

Specific Lot Number

J1274-31

Purification
Stability
Storage & Shipping

Affinity chromatography
1yr at -20°C from date of shipment
Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.