

## Anti-LIMK1

Rabbit Polyclonal Antibody

**Catalog # L04-63R**

Lot # B3216-55

### Cited Applications

ELISA, IF, WB

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

### Specificity

Recognizes the human LIM Kinase protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 630-647 of human LIM Kinase protein

### Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

### Stability

1yr at -20°C from date of shipment

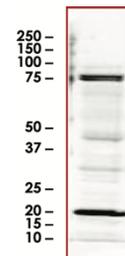
### Scientific Background

LIMK1 or LIM domain kinase 1 contains a unique combination of 2 N-terminal LIM domain and a C-terminal protein kinase domain. LIM domains are highly conserved cysteine-rich structures containing 2 zinc fingers that can bind to DNA/RNA as well as mediating protein-protein interactions (1). LIMK1 is thought to be a component of an intracellular signaling pathway that may be involved in brain development especially development of nerve cells. LIMK1 may play an important role in areas of the brain that are responsible for processing visual-spatial information (visuospatial constructive cognition). LIMK1 can regulate aspects of the cytoskeleton, the structural framework that helps to determine cell shape, size, and movement (2).

### References:

1. Davila, M. et al: LIM kinase 1 is essential for the invasive growth of prostate epithelial cells: implications in prostate cancer. *J Biol Chem.* 2003;19;278(38):36868-75.
2. Davila, M. et al: Expression of LIM kinase 1 is associated with reversible G1/S phase arrest, chromosomal instability and prostate cancer. *Mol Cancer.* 2007;8;6:40.

### Sample Data



Western blot using Anti-LIMK1 (1:500 dilution) in 18 µg of lysates from mouse brain. The doublet band at ~75 kDa may represent phosphorylated and non-phosphorylated forms of the protein.

## Anti-LIMK1

Rabbit Polyclonal Antibody

Catalog #

L04-63R

Lot #

B3216-55

Purification

Immunoaffinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**