LIMK1 monoclonal antibody (1B11)

The LIM kinases 1 and 2 (LIMK1 and LIMK2) are activated downstream of the Rho-ROCK, Rac-PAK, and Cdc42-MRCK pathways, thereby acting as network nodes in the signaling pathways that regulate the actin cytoskeleton. LIMK is required for matrix remodeling activities for path generation by leading cells in collective invasion. LIMK1 phosphorylates and inactivates the actin binding/depolymerizing factor cofilin, thereby stabilizing the actin cytoskeleton.

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

Ordering Information

Order Online »

ALX-803-343-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -20°C.

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name LIM domain kinase 1

Application ICC, IHC, IP, WB

Clone 1B11

Crossreactivity Does not cross-react with LIMK2.

Formulation Liquid. In PBS containing 10% glycerol and 0.02% sodium

azide.

Host Rat

Immunogen Recombinant LIM + PDZ domain of LIMK1.

lsotype lgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (20ng/ml)Immunohistochemistry

(paraffin sections 10μg/ml)Immunocytochemistry (10μg/ml)Western Blot (200ng/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Source Purified from concentrated hybridoma tissue culture

supernatant.

Species Reactivity

Species independent

Specificity

Recognizes full length GFP-LIMK1 and GFP-PDZ domain but not the GFP-LIM or GFP-kinase regions.

UniProt ID

P53667 (human)

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-