

# 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.
Isotype	IgG2a
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

<b>Specificity</b>	Recognizes full length GFP-LIMK1 and GFP-PDZ domain but not the GFP-LIM or GFP-kinase regions.
<b>UniProt ID</b>	P53667 (human)
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)