

## Lck, Rabbit Monoclonal Antibody

Publication Part no. MAN0004913

Rev. Date: 21 July 2011

Catalog Number: A13995

Store at -20°C

**Clonality:** Monoclonal  
**Volume:** 100 µL

**Host/Class:** Rabbit IgG  
**Reactivity:** Human Lck  
**Predicted Reactivity:** Human

### Product Description

Lck is a member of the Src family of protein tyrosine kinases, and is important in the regulation of growth and differentiation of eukaryotic cells. Src activity is regulated by tyrosine phosphorylation at two sites, but with opposing effects. Phosphorylation of Tyr<sup>416</sup> in the activation loop of the kinase domain upregulates enzyme activity. Phosphorylation of Tyr<sup>527</sup> in the carboxy-terminal tail by Csk renders the enzyme less active. p56lck (lymphocyte cellular kinase) has been shown to be a pivotal enzyme to both maturation of thymocytes and activation and proliferation of peripheral lymphocytes. The p56lck sequence appeared highly homologous to that of the oncogene p60c-src as did its exon-intron organisation.

### Product Specifications

**Immunogen:** Synthetic peptide corresponding to residues surrounding Tyr<sup>505</sup> of human Lck

**Alternate Names:** Lymphocyte-specific protein tyrosine kinase ,T-lymphocyte specific protein tyrosine kinase

**Apparent MW:** 56 kDa

**Gene ID:** 3932

**Protein Accession No.:** P06239

**Sequence Identity:** Human

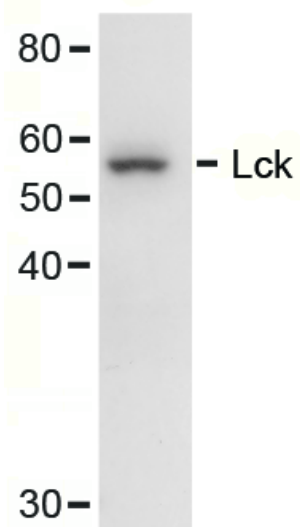
**Clone/PAD:** 73A5

**Lot:** See product label

### Product Applications

Application	Species	Test Material	Concentration
Western blotting	Human	Jurkat cells	1:1000
Immuno fluorescence (IF-IC)	Human	Jurkat cells	1:400
Flow Cytometry	Human	Jurkat cells	1:400

Additional applications for which this antibody has been used include: Immunoprecipitation (1:50 dilution).



**Figure 1** Western blot analysis of Jurkat cell extracts using Lck Rabbit Monoclonal Antibody.

**For research use only. Not for human or animal therapeutic or diagnostic use.**

Manufacturing Site • 7335 Executive Way • Frederick • MD 21704 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

## Storage and Handling

Store the antibody at  $-20^{\circ}\text{C}$ . Do not make aliquots of the antibody.

## Stability

Expires one year from date of receipt when stored as instructed.

## Storage Buffer

10 mM sodium HEPES (pH 7.5), 150 mM NaCl, 100  $\mu\text{g}/\text{mL}$  BSA and 50% glycerol.

## Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at [www.invitrogen.com/sds](http://www.invitrogen.com/sds).




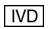



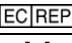
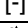
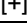


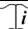
## Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to [www.invitrogen.com/support](http://www.invitrogen.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the box.

## Related Products

Product Name	Quantity	Catalog No.
iBlot <sup>®</sup> Dry Blotting System	1 unit	IB1001
WesternBreeze <sup>™</sup> Chromogenic Kit Anti-Rabbit	1 kit	WB7105
WesternBreeze <sup>™</sup> Chemiluminescent Kit, Anti-Rabbit	1 kit	WB7106
Goat anti-mouse (H+L), HRP conj.	1 mg	G21040
Goat anti-rabbit (H+L), HRP conj.	1 mg	G21234
Goat anti-mouse (H+L), AP conj.	1 mg	G21060
Goat anti-rabbit (H+L), AP conj.	1 mg	G21079
Nitrocellulose, 0.2 $\mu\text{m}$	20/pack	LC2000

### Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

## Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [outlicensing@lifetech.com](mailto:outlicensing@lifetech.com) or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit [www.invitrogen.com/support](http://www.invitrogen.com/support) or email [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)

*life*  
technologies<sup>™</sup>