

Lyn (pTyr508) Polyclonal Antibody

Store at -20°C

Catalog Number: A16552

Pub. No. MAN0009252 Rev. 1.00

Clonality: Polyclonal
Reactivity: Human Lyn

Amount: 100 µL
Species Reactivity: Human, Mouse

Host/Class: Rabbit

Product Description

Lyn, one of the Src family members, is predominantly expressed in hematopoietic cells (1). Two tyrosine residues have been reported to play a crucial role in the regulation of protein tyrosine kinases of the Src family. Autophosphorylation of Tyr396 (equivalent to Tyr416 of Src), located in the catalytic domain, correlates with enzyme activation. Csk-mediated phosphorylation of the carboxy-terminal Tyr507 (equivalent to Tyr527 of Src) inactivates the kinase. Tyrosine phosphorylation and activation of Lyn occurs upon association with cell surface receptors such as the B cell Ag receptor (BCR) and CD40 (2-4). Studies using knockout mice have shown that the net effect of Lyn deficiency is to render B cells hypersensitive to BCR stimulation (5-7), suggesting that the most critical role for Lyn in vivo is in the down-regulation of B cell responses. Lyn is also involved in controlling the migration and development of specific B cell populations (8).

Product Specifications

Immunogen:	A synthetic phosphopeptide corresponding to residues surrounding Tyr507 of human Lyn
Predicted Reactivity:	Rat
Apparent MW:	53, 56 kDa
Gene ID:	4067
Protein Accession No.:	P07948
Lot:	See product label

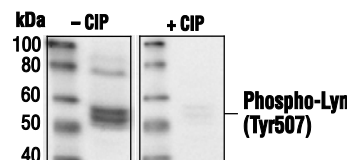


Figure 1 Western blot analysis of extracts from Ramos cells treated with anti-IgM antibody (12 µg/mL for 2 minutes), using Lyn (pTyr508) Polyclonal Antibody. The phospho-specificity of the antibody was characterized by treating the membrane without or with calf intestinal alkaline phosphatase (CIP) after western transfer.

Product Applications

Applications reported for this antibody include western blot. Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for each application.

Application	Recommended Dilution
Western blotting	1:1000

Storage and Handling

Store reagents at -20°C. Avoid repeated freezing and thawing. Do not aliquot the antibody.

Stability

When stored as instructed, expires one year from date of receipt unless otherwise indicated on product label.

Storage Buffer

10 mM sodium HEPES (pH 7.5), 150 mM NaCl, 100 µg/mL BSA, 50% glycerol.

References

1. Yamanashi, Y. et al. (1989) Proc. Natl. Acad. Sci. USA 86, 6538-6542.
2. Yamanashi, Y. et al. (1991) Science 251, 192-194.
3. Burkhardt, A.L. et al. (1991) Proc. Natl. Acad. Sci. USA 88, 7410-7414.
4. Ren, C.L. et al. (1994) J. Exp. Med. 179, 673-680.
5. Wang, J. et al. (1996) J. Exp. Med. 184, 831-838.
6. Chan, V.W. et al. (1997) Immunity 7, 69-81.
7. Hibbs, M. L. et al. (1995) Cell 83, 301-311.
8. Seo, S.J. et al. (2001) J. Immunol. 166, 3710-3723.








Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit www.lifetechnologies.com/support.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Explanation of symbols

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		

DISCLAIMER: LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

©2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

28 August 2013

