

Afadin (pS1718) Polyclonal Antibody

Store at -20°C

Catalog Number: A16692

Pub. No. MAN0009172 Rev. 1.00

Clonality: Polyclonal

Amount: 100 µL

Host/Class: Rabbit

Reactivity: Human Afadin

Species Reactivity: Human

Product Description

Intercellular junctions play essential roles in tissue integrity and maintenance of cell polarity. Tight junctions (TJs) form a continuous barrier to fluids (1), while adherens junctions (AJs) link cells in a continuous sheet (2). Afadin is an actin filament-binding protein that binds to nectin, forming a connection to the actin cytoskeleton (3). AJs are formed when nectin assembles cadherin at cell-cell adhesion sites, which are then involved in the formation and maintenance of TJs (4,5). Afadin has two splice variants, l- and s-afadin. l-afadin is ubiquitously expressed, while the shorter s-afadin, lacking one of the three proline-rich regions and a carboxyl-terminal F-actin binding region, is expressed predominantly in neural tissue (6). Afadin is also involved in controlling the directionality of cell movement when it is localized at the leading edge of moving cells (7,8).

Product Specifications

Immunogen:	A synthetic phosphopeptide corresponding to residues surrounding Ser1718 of human afadin
Predicted Reactivity:	Mouse, Rat, Monkey, Dog
Apparent MW:	205 kDa
Gene ID:	4301
Protein Accession No.:	P55196
Lot:	See product label

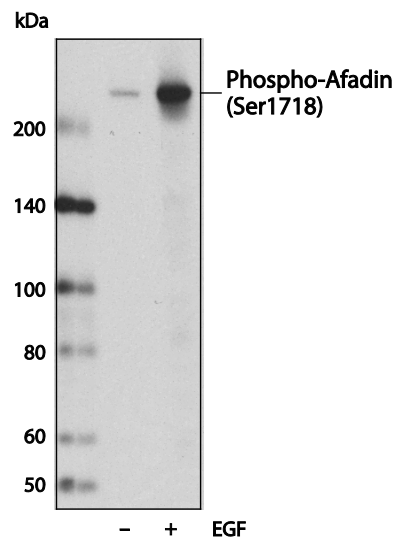


Figure 1 Western blot analysis of extracts from A431 cells, untreated or EGF-treated (100 ng/mL 5 minutes), using Afadin (pS1718) Polyclonal Antibody.

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.

For research use only. Not for use in diagnostic procedures

Manufacturing Site • 7335 Executive Way • Frederick • MD 21704 • E-mail: techsupport@lifetech.com

References

1. Shin, K. et al. (2006) Annu Rev Cell Dev Biol 22, 207-35.
2. Harris, T.J. and Tepass, U. (2010) Nat Rev Mol Cell Biol 11, 502-14.
3. Ikeda, W. et al. (1999) J Cell Biol 146, 1117-32.
4. Sato, T. et al. (2006) J Biol Chem 281, 5288-99.
5. Ooshio, T. et al. (2007) J Cell Sci 120, 2352-65.
6. Mandai, K. et al. (1997) J Cell Biol 139, 517-28.
7. Miyata, M. et al. (2009) J Cell Sci 122, 4319-29.
8. Miyata, M. et al. (2009) J Biol Chem 284, 24595-609.








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

