

## NHERF1 Polyclonal Antibody

Store at -20°C

Catalog Number: A16581

Pub. No. MAN0009299 Rev. 1.00

**Clonality:** Polyclonal  
**Reactivity:** Human NHERF1

**Amount:** 100 µL  
**Species Reactivity:** Human

**Host/Class:** Rabbit

### Product Description

NHERF1 (Na<sup>+</sup>/H<sup>+</sup> exchanger regulatory factor) is one of several related PDZ domain-containing proteins (1). It was first identified as a necessary co-factor for cyclic AMP-associated inhibition of NHE3 (2). NHERF1 is a multifunctional adaptor protein that interacts with receptors and ion transporters via its PDZ domains, and with the ERM family of proteins, including merlin, via its carboxy-terminus (2,3). NHERF1 has been found to promote the formation of a ternary complex containing PTEN, NHERF1, and PDGFR, and may be involved in inhibition of PI3K/Akt signaling by recruiting PTEN to PDGFR (4). Other proposed functions of NHERF1 include testicular differentiation, endosomal recycling, membrane targeting, protein sorting, and trafficking (6).

### Product Specifications

**Immunogen:** A synthetic peptide surrounding Ala310 of human NHERF1

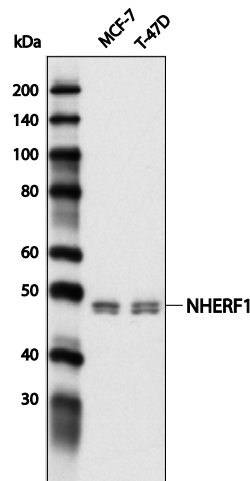
**Alternate Names:** EBP-50

**Apparent MW:** 50 kDa

**Gene ID:** 9368

**Protein Accession No.:** O14745

**Lot:** See product label



**Figure 1** Western blot analysis of extracts from MCF-7 and T-47D cells using NHERF1 Polyclonal Antibody.

### Product Applications

Applications reported for this antibody include western blot, and immunofluorescence (immunocytochemistry).

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
Immunofluorescence (immunocytochemistry)	1:50

### 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

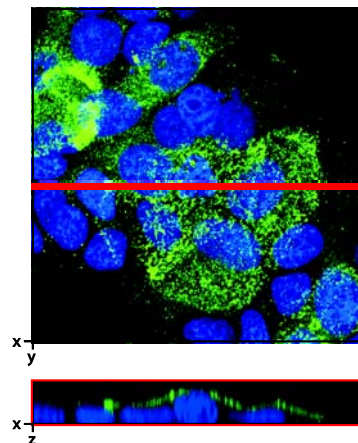
1. Donowitz, M. et al. (2005) J Physiol 567, 3-11.
2. Voltz, J.W. et al. (2001) Oncogene 20, 6309-14.
3. Stemmer-Rachamimov, A.O. et al. (2001) Am J Pathol 158, 57-62.
4. Takahashi, Y. et al. (2006) EMBO J 25, 910-20.
5. Weinman, E.J. et al. (2000) Am J Physiol Renal Physiol 279, F393-9.

## Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit [www.lifetechnologies.com/support](http://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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).



**Figure 2** Projected confocal z-stack of MCF-7 cells showing staining at the apical membrane when labeled with NHERF1 Polyclonal Antibody (green). A fluorescent DNA dye was used to produce the blue pseudocolor.

## 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](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

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

28 August 2013

*life*  
technologies™