

DDX3 Polyclonal Antibody

Store at -20°C

Catalog Number: A16543

Pub. No. MAN0009242 Rev. 1.00

Clonality: Polyclonal

Amount: 100 µL

Host/Class: Rabbit

Reactivity: Human DDX3

Species Reactivity: Human, Mouse, Rat, Hamster, Monkey

Product Description

The DEAD box family of RNA helicases is characterized in part by a common D-E-A-D amino acid motif. The family is composed of a growing number of proteins found in a wide range of organisms from bacteria to mammals. DEAD helicases have distinct biological functions in RNA metabolism and ribonucleoprotein (RNP) processing (1,2). DDX3 is a DEAD box family RNA helicase with diverse cellular functions including export of viral transcripts (3), and proliferation (4,5). DDX3 is phosphorylated at Thr204 and Thr323 by the mitotic cyclin dependent kinase, cyclin B/cdc2. This phosphorylation is thought to cause a loss of DDX3 function and a concomitant repression of ribosome biogenesis and translation in mitosis (6).

Product Specifications

Immunogen:	A synthetic peptide corresponding to the amino-terminal sequence of human DDX3
Apparent MW:	75 kDa
Gene ID:	1654
Protein Accession No.:	O00571
Lot:	See product label

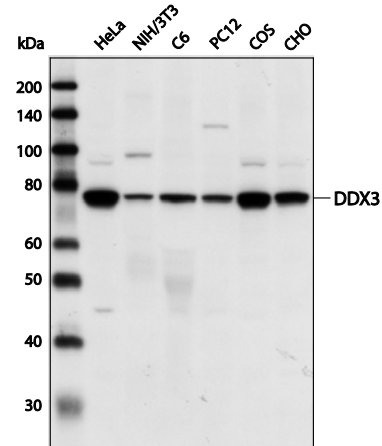


Figure 1

Western blot analysis of extracts from various cell types using DDX3 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.

References

1. Rocak, S. and Linder, P. (2004) Nat Rev Mol Cell Biol 5, 232-41.
2. Linder, P. (2006) Nucleic Acids Res 34, 4168-80.
3. Yedavalli, V.S. et al. (2004) Cell 119, 381-92.
4. Chang, P.C. et al. (2006) Oncogene 25, 1991-2003.
5. Chao, C.H. et al. (2006) Cancer Res 66, 6579-88.
6. Sekiguchi, T. et al. (2007) Biochem Biophys Res Commun 356, 668-73.








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

