

mTOR (human FRB Domain) polyclonal antibody

mTOR (FKBP12-rapamycin associated protein (FRAP)) is one of a family of proteins involved in cell cycle progression, DNA recombination, and DNA damage detection. It has been shown to associate with the immunophilin FKBP12 in a rapamycin-dependent fashion. FKBP12-rapamycin-binding (FRB) domain is a conserved 11kDa region necessary for FKBP12-rapamycin binding. FRB is one of several functional domains of mTOR: HEAT (Huntington elongation factor 1A-protein phosphatase 2A-A subunit-TOR); FAT (FRAP, ATM, TTRAP2); PIKK (PI3-kinase-related kinase); RD (regulatory domain); and FATC (FAT, C terminal).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 22

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-215-065-1	1Vial
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

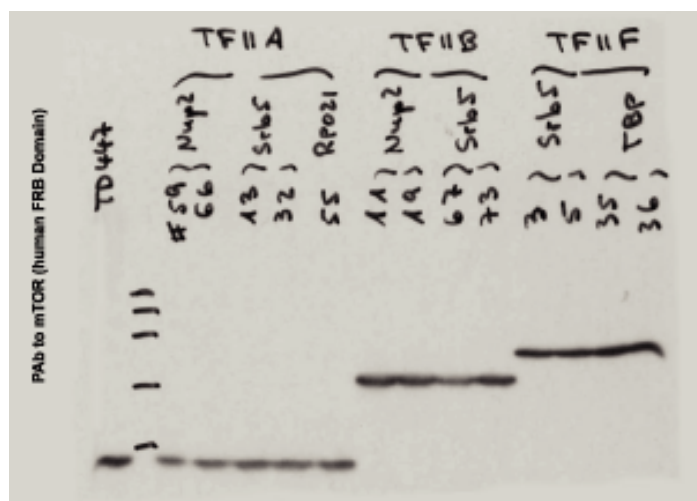


Figure: WB using PAb to mTOR (human FRB Domain) (Prod. No. [ALX-215-065](#)) on total yeast extracts derived from cells expressing FRB-tagged genes coding for three different transcription factors. Separation SDS-PAGE. "", mTOR (human FRB Domain) polyclonal antibody WB, mTOR (human FRB Domain) polyclonal antibody Western Blot, yes ALX-215-065, 215-

065_2_WB.gif, <http://static.enzolifesciences.com/fileadmin/fi> 065_2_WB.gif, **Figure 2:** Western Blot of mTOR (human FRB Domain)

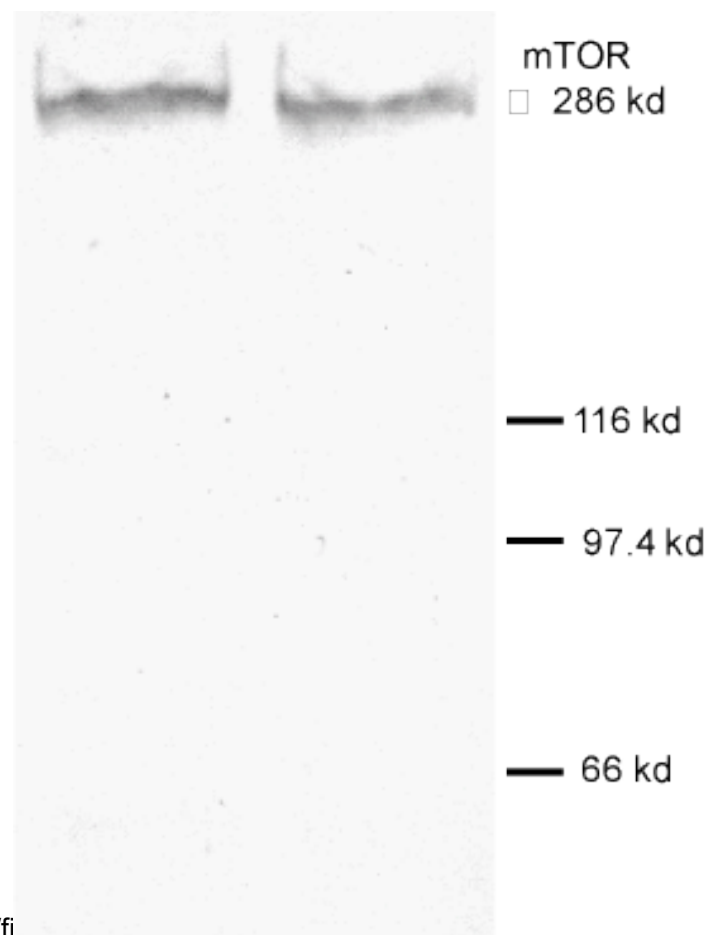


Figure 2: Western Blot of mTOR (human FRB Domain), pAb (Prod. No. ALX-215-065) with HeLa cells.

Method: Two ml of an exponential HeLa suspension culture ($2-4 \times 10^5$ cells) was washed twice with PBS quickly without resuspension of the pellet by spinning in a Eppendorf tube and finally with 25% TCA. TCA was carefully removed and pellet resuspended 100 μ l FSB + β ME (final sample buffer) and the pH adjusted with about a few μ l of 1M Tris base (≈ 10 μ l). The sample was heated at 100°C for 5 min. and sonicated. SDS PAGE was at 7.5%, 20 & 25 μ l loaded in lane 1 and 2, respectively. The TCA step is performed to prevent proteolytic breakdown of the very large mTOR.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name mTOR (human FKBP12-rapamycin-binding domain)

Application ChIP, ICC, WB

Application Notes Among other applications this antibody can be used to control FRB-tagging of genes and the level of expression thereof in yeast cells (see H. Haruki, et al. 2008).

Formulation Lyophilized from neat serum containing sodium azide.

Host Rabbit

Immunogen Recombinant human mTOR (FKBP12-rapamycin-binding (FRB) domain).

Recommendation Dilutions/Conditions Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Reconstitution Reconstitute with 50µl water.

Species Reactivity Human

Specificity Recognizes the FRB domain of fusion proteins expressed in *Saccharomyces cerevisiae* (see H. Haruki, et al. 2008) and the human FRB domain of mTOR.

UniProt ID P42345

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: January 23, 2025

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.