

Anti-mTOR

Rabbit Polyclonal Antibody

Catalog # F17-63BR

Lot # O2361-33

Cited Applications

ELISA, WB, ICC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the mTOR protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

mTOR antibody was raised against a 15 amino acid synthetic peptide from near the amino terminus of human mTOR

Formulation

PBS + 0.02% sodium azide.

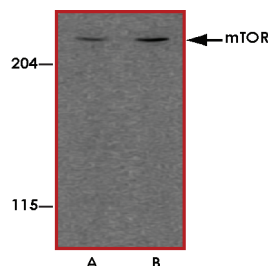
Stability

1yr at -20°C from date of shipment.

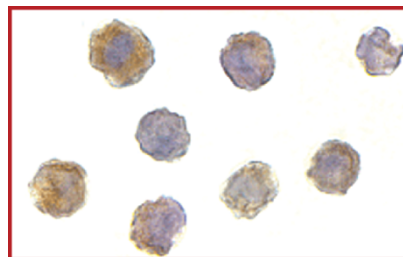
References

1. Shamji AF et al: Integration of growth factor and nutrient signaling: implications for cancer biology. *Mol. Cell* 2003; 12:271-80.
2. Sabatini DM et al: RAFT1: a mammalian protein that binds to FKPI2 in a rapamycin-dependent fashion and is homologous to yeast TORs. *Cell* 1994; 78:35-43.
3. Cardenas ME et al: Molecular mechanisms of immunosuppression by cyclosporine, FK506, and rapamycin. *Curr. Opin. Nephrol. Hypertens.* 1995; 4:472-7.
4. Fingar DC et al: Target of rapamycin (TOR): an integrator of nutrient and growth factor signals and coordinator of cell growth and cell cycle progression. *Oncogene* 2004; 23:3151-71.

Sample Data



Western blot analysis of mTOR in L1210 cell lysate with mTOR antibody at (A) 1 and (B) 2 ug/mL.



Immunocytochemistry of mTOR in L1210 cells with mTOR antibody at 2 ug/mL.

Anti-mTOR

Rabbit Polyclonal Antibody

Catalog Number

F17-63BR

Specific Lot Number

O2361-33

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.