

Anti-mTOR

Rabbit Polyclonal Antibody

Catalog # F17-63R-100

Lot # H257-1

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the mTOR protein

Cross Reactivity

- Western blot of mTOR from human cells

mTOR from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 2446-2450 of human mTOR

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

Scientific Background

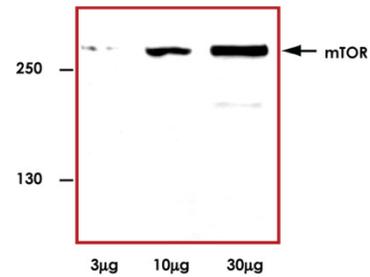
The mammalian target of rapamycin, mTOR, also known as FRAP or RAFT, is a Ser/Thr protein kinase. mTOR acts as a sensor for ATP and amino acids, balancing the availability of nutrients and cell growth (1). When sufficient nutrients are available, mTOR responds to a phosphatidic acid-mediated signal, transmits a positive signal to p70 S6 kinase and participates in the inactivation of the eIF4E inhibitor, 4E-BP1 (2).

References

1. Sabers CJ et al: Isolation of a protein target of the FKBP12-rapamycin complex in mammalian cells. *J Biol Chem.* 270(2): 815-22, 1995.
2. Fang Y. et al: Phosphatidic acid-mediated mitogenic activation of mTOR signaling. *Science.* 294(5548): 1942-5, 2001.

Sample Data

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



Representative western blot with Anti-mTOR (1:1000) using 3 µg, 10 µg, and 30 µg of HeLa cell lysate.

Anti-mTOR

Rabbit Polyclonal Antibody

Catalog Number	F17-63R-100
Quantity	100 µg
Specific Lot Number	H257-1
Purification	Affinity chromatography
Format	100 µg in 100 µl
Concentration	1.0 µg/µL
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.

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