

## Anti-phospho-mTOR (Ser2448)

Rabbit Polyclonal Antibody

**Catalog # F17-65R**

Lot # Z2014-25

### Cited Applications

IHC

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

### Specificity

Recognizes the mTOR protein phosphorylated at serine 2448

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

The antiserum was produced against synthesized phosphopeptide derived from human mTOR around the phosphorylation site of serine 2448 (T-D-SP-Y-S).

### Formulation

Phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

### Stability

1yr at -20°C from date of shipment

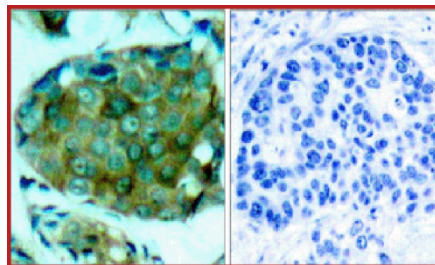
### Scientific Background

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.
3. Holz MK, et.al. (2005) J Biol Chem ;280:26089-26093.
4. Chiang GG, et.al. (2005) J Biol Chem ;280: 25485-25490.
5. Mothe-Satney I, et.al. (2004) J Biol Chem ;279: 42628-42637.
6. Bolster DR, et.al. (2003) J Physiol ;553:213-220.

### Sample Data



Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using Anti-phospho-mTOR (Ser2448).

## Anti-phospho-mTOR (Ser2448)

Rabbit Polyclonal Antibody

Catalog # F17-65R

Lot # Z2014-25

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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