

Anti-Raptor

Rabbit Polyclonal Antibody

Catalog # R47-363BR

Lot # O2361-49

Cited Applications

E, WB, ICC, IF

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

Specificity

Recognizes the Raptor protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Raptor polyclonal antibody was raised against a 13 amino acid synthetic peptide from near the amino-terminus of human Raptor

Formulation

PBS + 0.02% sodium azide.

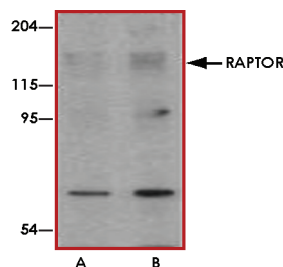
Stability

1yr at -20°C from date of shipment.

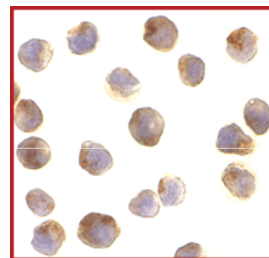
References

1. Shamji AF et al: Integration of growth factor and nutrient signaling: implications for cancer biology. *Mol. Cell* 2003; 12:271-80.
2. 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.
3. Yonezawa K et al: Raptor, a binding partner of target of rapamycin. *Biochem. Biophys. Res. Commun.* 2004; 313:437-441.
4. Hara K et al: Amino acid sufficiency and mTOR regulate p70 S6 kinase and eIF-4E BP1 through a common effector mechanism. *J. Biol. Chem.* 1998; 273:14484-94.

Sample Data



Western blot analysis of Raptor in L1210 cell lysate with Raptor antibody at (A) 2 and (B) 4 µg/mL.



Immunocytochemistry of Raptor in L1210 cells with Raptor antibody at 10 µg/mL.

Anti-Raptor

Rabbit Polyclonal Antibody

Catalog Number

R47-363BR

Specific Lot Number

O2361-49

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.