

## Anti-GAB1

Rabbit Polyclonal Antibody

**Catalog # G03-63R**

Lot # M132-5

### Cited Applications

- Western blot (1:1000)

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

### Specificity

Recognizes the GAB1 protein

### Cross Reactivity

- Western blot human and mouse GAB1

*GAB1 from other species may also be detectable*

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 625-629 of human GAB1

### Formulation

PBS, pH7.4, 0.02%NaN<sub>3</sub>, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) 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

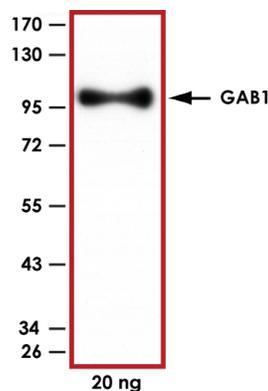
GAB1 is a member of the IRS1-like multisubstrate docking protein family and is a direct substrate of the epidermal growth factor receptor and the insulin receptor (1). GAB1 is tyrosine phosphorylated upon stimulation of cells with various cytokines, growth factors, and antigen receptor. Tyrosine phosphorylation of GAB1 mediates interaction with several proteins that contain SH2 domains such as SHP2 and phosphatidylinositol 3-kinase. GAB1 is an important mediator of branching tubulogenesis and plays a central role in cellular growth response, transformation and apoptosis (2).

### References

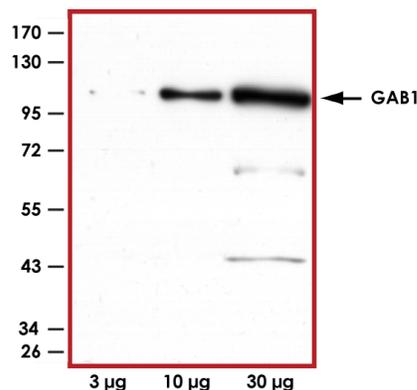
- Holgado-Madruga M, et al: A Grb2-associated docking protein in EGF- and insulin-receptor signalling. *Nature* 379: 560-564, 1996.
- Kiyatkin A, et al: Scaffolding protein Grb2-associated binder 1 sustains epidermal growth factor-induced mitogenic and survival signaling by multiple positive feedback loops. *J Biol Chem.* 2006 Jul 21;281(29):19925-38.

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)

### Sample Data



Representative Western blot with Anti-GAB1 (1:1000) using 20 ng of His-tagged human recombinant GAB1.



Representative Western blot with Anti-GAB1 (1:1000) using 3 µg, 10 µg and 30 µg of Jurkat cell lysate.

## Anti-GAB1

Rabbit Polyclonal Antibody

Catalog Number

G03-63R

Specific Lot Number

M132-5

Purification

Affinity chromatography

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.