

## Anti-IRS1

Rabbit Polyclonal Antibody

**Catalog # I40-363R**

Lot # O2109-42

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the IRS1 protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

IRS1 antibody was raised against a 16 amino acid synthetic peptide from near the center of human IRS1

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

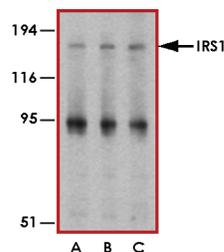
### Scientific Background

IRS1 is the substrate for the insulin tyrosine kinase receptor and is found in a variety of insulin-responsive cells and tissues. IRS1 protein has no intrinsic enzymatic activity but acts as a docking protein, via the SH2 domains, for mediating the insulin downstream signaling events. IRS1 has been shown to associate with the 14-3-3 family of proteins and this could play a role in the regulation of insulin sensitivity by interrupting the association between the insulin receptor and IRS1 (1). IRS1 may be associated with colorectal cancer and diet and related factors may affect the risk by modifying plasma insulin levels. Thus, the inter-individual variation in insulin signaling mediated by IRS1 may play a plausible role in the development of colorectal cancer (2).

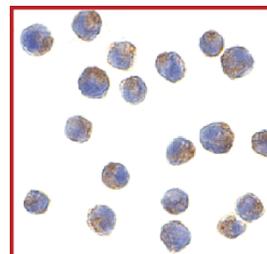
### References

- Ogihara, T. et al: 14-3-3 protein binds to insulin receptor substrate-1, one of the binding sites of which is in the phosphotyrosine binding domain. *J. Biol. Chem.* 272: 25267-25274, 1997.
- Slattery, M.L. et al: Genetic variation in IGF1, IGFBP3, IRS1, IRS2 and risk of breast cancer in women living in Southwestern United States. *Breast Cancer Res Treat.* 2007.

### Sample Data



Western blot analysis of IRS1 in PC-3 cell lysate with IRS1 antibody at (A) 1, (B) 2 and (C) 4 µg/ml.



Immunocytochemistry of IRS1 in P815 cells with IRS1 antibody at 2 µg/ml.

## Anti-IRS1

Rabbit Polyclonal Antibody

Catalog Number

**I40-363R**

Specific Lot Number

**O2109-42**

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.**