

## Anti-QIK

Rabbit Polyclonal Antibody

**Catalog # S15-63R**

Lot # O2121-44

### Cited Applications

WB, ELISA

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

### Specificity

Recognizes the QIK protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

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

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

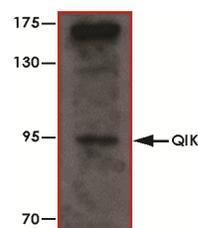
### Scientific Background

QIK is a serine/threonine protein kinase that contains an N-terminal kinase domain, a central domain with ubiquitin-associate motif, and a C-terminal PKA phosphorylation site. QIK can phosphorylate IRS1 and overexpression of QIK in adipocyte elevates the phosphorylation of IRS1 (1). The QIK-mediated phosphorylation of IRS1 may modulate the efficiency of insulin signal transduction and could be responsible for insulin resistance associated with diabetes (1). Insulin disrupts TORC2 activity by induction of QIK which then stimulates the phosphorylation and cytoplasmic translocation of TORC2. Phosphorylated TORC2 is subsequently degraded by the 26S proteasome (2).

### References

1. Horike N, et al: Adipose-specific expression, phosphorylation of ser794 in insulin receptor substrate-1, and activation in diabetic animals of salt-inducible kinase-2. J. Biol. Chem. 278: 18440-18447, 2003.
2. Dentin R, et al: Insulin modulates gluconeogenesis by inhibition of the coactivator TORC2. Nature 449: 366-369, 2007.

### Sample Data



Western blot analysis of QIK in SW480 cell lysate with QIK antibody at 1 µg/ml.

## Anti-QIK

Rabbit Polyclonal Antibody

Catalog Number

S15-63R

Specific Lot Number

O2121-44

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.**