

## Anti-DRAK2 (STK17B)

Rabbit Polyclonal Antibody

Catalog # S34-63R

Lot # O2109-13

### Cited Applications

WB, ELISA, ICC

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

### Specificity

Recognizes the DRAK2 (STK17B) protein and has no cross responses to DAP or ZIP kinases.

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

DRAK2 (STK17B) antibody was raised against a peptide corresponding to amino acids 351 to 365 of human DRAK2 (STK17B)

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

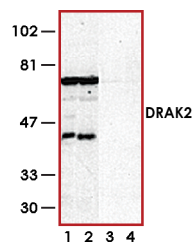
### Scientific Background

DRAK2 is a member of the serine/threonine kinase family and is related to death-associated protein kinase that triggers apoptosis (1). DRAK2 is selectively important for T-cell survival and inhibition of DRAK2 has therapeutic potential for autoimmune disease (2). T-cell survival depends on a balance of T-cell receptor and co-stimulatory signals and deficiency of DRAK2 can affect autoimmune disease susceptibility without generalized suppression of the immune system (3).

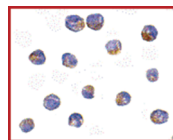
### References

1. Sanjo, H. et.al: DRAKs, novel serine/threonine kinases related to death-associated protein kinase that trigger apoptosis. J. Biol. Chem. 273: 29066-29071, 1998.
2. Ramos, S. J. et.al: Enhanced T cell apoptosis within Drak2-deficient mice promotes resistance to autoimmunity. J. Immun. 181: 7606-7616, 2008.
3. McGargill, M. A. et.al: Drak2 regulates the survival of activated T cells and is required for organ-specific autoimmune disease. J. Immun. 181: 7593-7605, 2008.

### Sample Data



Western blot analysis of DRAK2 (STK17B) in Jurkat (1,3) and Raji (2,4) cell lysate in the absence (1,2) or presence (3,4) of blocking peptide with DRAK2 (STK17B) antibody at 1:500 dilution.



Immunocytochemistry of DRAK2 (STK17B) in Jurkat cells with DRAK2 (STK17B) antibody at 10 ug/ml.

## Anti-DRAK2 (STK17B)

Rabbit Polyclonal Antibody

Catalog Number

S34-63R

Specific Lot Number

O2109-13

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.**