

## Anti-PIAS3

Rabbit Polyclonal Antibody

**Catalog # P282-63R**

Lot # O2361-39

### Cited Applications

E, WB, IHC, IF

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

### Specificity

Recognizes the PIAS3 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

PIAS3 antibody was raised against a 12 amino acid synthetic peptide near the carboxy terminus of human PIAS3

### Formulation

PBS + 0.02% sodium azide.

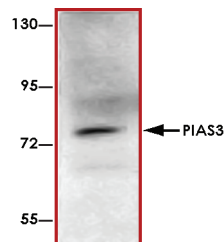
### Stability

1yr at -20°C from date of shipment.

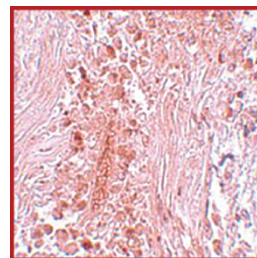
### References

1. Liu B et al: Inhibition of Stat1-mediated gene activation by PIAS1. Proc. Natl. Acad. Sci. USA 1998; 95: 10626-31.
2. Shuai K et al: Regulation of gene-activation pathways by PIAS proteins in the immune system. Nat. Rev. Immunol. 2005; 5:593-605.
3. Chung CD et al: Specific inhibition of Stat3 signal transduction by PIAS3. Science 1997; 278:1803-5.
4. Wang L B et al: Differential PIAS3 expression in human malignancy. Oncol. Rep. 2004; 11:1319-24.
5. Zhang Q et al: Multilevel dysregulation of STAT3 activation in anaplastic lymphoma kinase-positive T/null-cell lymphoma. J. Immunol. 2002; 168:466-7.

### Sample Data



Western blot analysis of PIAS3 in K562 cell lysate with PIAS3 antibody at 1 µg/mL.



Immunohistochemistry of PIAS3 in human breast carcinoma tissue with PIAS3 antibody at 5 µg/mL.

## Anti-PIAS3

Rabbit Polyclonal Antibody

Catalog Number

**P282-63R**

Specific Lot Number

**O2361-39**

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.**