

## Anti-phospho-p53 (Ser392)

Rabbit Polyclonal Antibody

**Catalog # P05-363R**

Lot # J1274-26

### Cited Applications

WB

Suggested Dilutions:  
WB 1:1,000

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

### Specificity

Recognizes the p53 protein phosphorylated at serine 392

### Cross Reactivity

Human and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser392 conjugated toKLH

### Formulation

100 µl in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

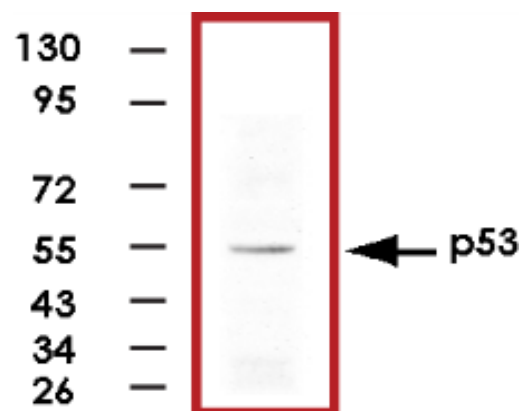
### Scientific Background

p53 is a key regulator of cell growth and acts as a tumor suppressor gene. Wild-type p53 gene can suppress transformation of rat embryo fibroblasts in cell culture by other oncogenes such as the adenovirus E1A and Ras (1). DNA tumor viruses such as SV40 large T antigen and the adenovirus E1A plus E1B-55Kd proteins bind to p53 and inactivate its tumor suppressor activities leading to cellular transformation. Mutational inactivation of the p53 gene is detected in more than 50% of human cancers. Mutation of p53 renders cancer cells more resistant to current cancer therapies due to lack of p53-mediated apoptosis (2).

### References

1. Levine, A J. et al: Tumor suppressor genes: the p53 and retinoblastoma sensitivity genes and gene products. *Biochim Biophys Acta*. 1990 Jun 1;1032(1):119-36.
2. Sun, Y. p53 and its downstream proteins as molecular targets of cancer. *Mol Carcinog*. 2006 Jun;45(6):409-15.

### Sample Data



Western blot of rat brain nuclear fraction lysate showing specific immunolabeling of the ~53kDa p53 phosphorylated at Ser392 (Control). The phosphospecificity of this labeling is shown in the second lane (lambda-phosphatase: lambda-Ptase). The blot is identical to the control except that it was incubated in lambda-Ptase (1200 units for 30 min) before being exposed to the Anti-phospho-p53 (Ser392)antibody. The immunolabeling is completely eliminated by treatment with lambda-Ptase.

## Anti-phospho-p53 (Ser392)

Rabbit Polyclonal Antibody

Catalog Number

P05-363R

Specific Lot Number

J1274-26

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.**