

Anti-p53

Rabbit Polyclonal Antibody

Catalog # P05-63R

Lot # L020-3

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the p53 protein

Cross Reactivity

- Western blot p53 from human cells

p53 from other species may also be detectable

Host/Isotype/Clone#

Rabbit

Immunogen

Human p53 peptide (15-40aa)-KLH conjugates

Formulation

TBS pH 7.2, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

Scientific Background

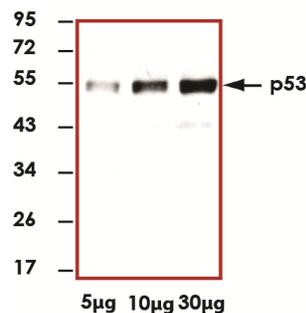
p53 is a key regulator of cell growth and acts as a tumor suppressor gene. Wild-type p53 gene can suppress transformation of rat embryo fibroblasts in cell culture by other oncogenes such as the adenovirus E1A and Ras (1). Mutational inactivation of the p53 gene is detected in more than 50% of human cancers. Mutation of p53 renders cancer cells more resistant to current cancer therapies due to lack of p53-mediated apoptosis (2).

References

- 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.
- Sun, Y. p53 and its downstream proteins as molecular targets of cancer. *Mol Carcinog*. 2006 Jun;45(6):409-15.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Sample Data



Representative western blot with Anti-p53 (1:1000) using 5 µg, 10 µg, and 30 µg of HEK 293 cell lysate.

Anti-p53

Rabbit Polyclonal Antibody

Catalog Number

P05-63R

Specific Lot Number

L020-3

Purification

Affinity purified by p53 peptide antigen

Format

100 µg in 400 µl

Concentration

0.25 µg/µl

Stability

1 yr 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.