

Anti-p33ING1

Goat Polyclonal Antibody

Catalog # **I32-363G**

Lot # B3216-65

Cited Applications

ELISA, WB

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

Specificity

Recognizes the p33 ING1 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Goat, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 285-296 of Human p33 ING1 protein (inhibitor of growth family, member 1)

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

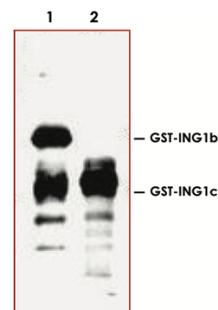
Scientific Background

p33 ING1 is a tumor suppressor protein that can induce cell growth arrest and apoptosis. The encoded protein is a nuclear protein that physically interacts with the tumor suppressor protein TP53 and is a component of the p53-signaling pathway. Reduced expression and rearrangement of this gene have been detected in various cancers. Multiple alternatively spliced transcript variants encoding distinct isoforms have been reported. The accession number listed below is for variant (4) that encodes the longest isoform (D). Other synonyms for p33 ING1 include p33, p47, p33ING1, p24ING1c, p33ING1b, p47ING1a, growth inhibitor ING1, inhibitor of growth 1, tumor suppressor ING1 and growth inhibitory protein ING1.

References:

1. Campos, E.I., et al.: Biological functions of the ING family tumor suppressors. *Cell. Mol. Life Sci.* 2004; 61 (19-20): 2597-2613.
2. Cheung, K.J., et al.: p33(ING1) enhances UVB-induced apoptosis in melanoma cells. *Exp. Cell Res.* 2002; 279 (2): 291-298.
3. Vieyra, D., et al.: Human ING1 proteins differentially regulate histone acetylation. *J. Biol. Chem.* 2002; 277 (33): 29832-29839.

Sample Data



Western blot analysis using anti-p33ING1 (1:1,000 dilution) to detect over expressed Human ING1 present in cell nuclear extracts. This western blot shows reactivity with purified recombinant GST tagged human ING1b (lane 1) and ING1c protein (lane 2).

Anti-p33ING1

Goat Polyclonal Antibody

Catalog #	I32-363G
Lot #	B3216-65
Purification	Immunoaffinity 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
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.