

Anti-p28ING5

Goat Polyclonal Antibody

Catalog # **I32-363EG**

Lot # B3216-63

Cited Applications

ELISA, WB

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

Specificity

Recognizes the p85 ING5 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Goat, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 127-140 of human p28 ING5 protein (inhibitor of growth family, member 5)

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

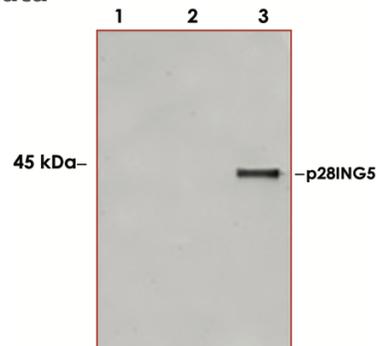
Scientific Background

p28 ING5 is a tumor suppressor protein similar to ING1 that can interact with TP53, inhibit cell growth, and induce apoptosis. This protein contains a PHD-finger, which is a common motif in proteins involved in chromatin remodeling. This protein can bind TP53 and EP300/p300, a component of the histone acetyl transferase complex, suggesting its involvement in TP53-dependent regulatory pathway. Multiple alternatively spliced transcript variants have been observed. The accession number listed below is for variant (1) encodes the longest isoform

References:

1. Shiseki, M., et al.: p29ING4 and p28ING5 bind to p53 and p300, and enhance p53 activity. *Cancer Res.* 2003; 63 (10): 2373-2378.

Sample Data



Western blot analysis is shown using Anti-p28ING5 (1:500 dilution) to detect over expressed Human ING5 present in HeLa cell nuclear extracts. This western blot shows reactivity with purified recombinant TAP tagged human ING5 protein (lane 3) and does not recognize TAP tagged ING4 on the same membrane (lane 2). A mock purification is shown in lane 1

Anti-p28ING5

Goat Polyclonal Antibody

Catalog # **I32-363EG**

Lot # **B3216-63**

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.