

p33ING1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # I32-30H

Lot # S168-4

Product Description

Recombinant full-length human p33ING1 was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_198219](#).

Gene Aliases

ING1, p33, p47, p24ING1c, p33ING1b, p47ING1a

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.2mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

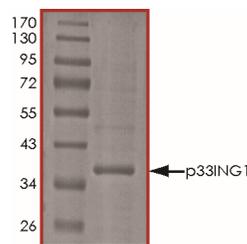
Scientific Background

p33ING1 displays characteristics of a tumour suppressor protein and can induce cell growth arrest and apoptosis. p33ING1 is a nuclear protein and acute expression of p33ING1 inhibits cell growth. p33ING1 physically interacts with the tumor suppressor protein TP53 and is a component of the p53 signaling pathway. Mutation of p33ING1 gene occurs in neuroblastoma cells and reduced expression is observed in some breast cancer cell lines (1). It has been proposed that p33ING1 can act as a potent growth regulator in normal and in established cells and plays a role as a candidate tumour suppressor gene whose inactivation may contribute to the development of cancers (2).

References

1. Garkavtsev, I. Et al: Suppression of the novel growth inhibitor p33ING1 promotes neoplastic transformation. Nat Genet, 1999; 23(3):373.
2. Goeman, F. Et al: Growth Inhibition by the Tumor Suppressor p33ING1 in Immortalized and Primary Cells: Involvement of Two Silencing Domains and Effect of Ras.

Purity



The purity of p33ING1 was determined to be **>95%** by densitometry. Approx. MW **38kDa**.

p33ING1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #	I32-30H
Lot #	S168-4
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.