

p33ING2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # 132-30BG

Lot # S250-4

Product Description

Recombinant full-length human p33ING2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001564](#).

Gene Aliases

ING1L, ING2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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.

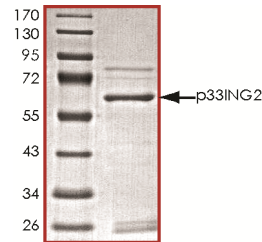
Scientific Background

p33ING2 is a member of the inhibitor of growth (ING) family that is associated with and modulates the activity of histone acetyltransferase (HAT) and histone deacetylase (HDAC) complexes and function in DNA repair and apoptosis. P33ING2 is a DNA damage-inducible gene that negatively regulates cell proliferation through activation of p53 by enhancing its acetylation (1). P33ING2 is required for initial DNA damage sensing and chromatin remodeling in the nucleotide excision repair process (2).

References

1. Nagashima, M.et.al: DNA damage-inducible gene p33ING2 negatively regulates cell proliferation through acetylation of p53. Proc. Nat. Acad. Sci. 98: 9671-9676, 2001.
2. Wang, J.et.al: The novel tumor suppressor p33ING2 enhances nucleotide excision repair via induction of histone H4 acetylation and chromatin relaxation. Cancer Res. 66: 1906-1911, 2006.

Purity



The purity of p33ING2 was determined to be **>75%** by densitometry. Approx. MW **62kDa**.

p33ING2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	132-30BG
Specific Lot Number	S250-4
Purity	>75%
Concentration	0.2 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.