

AXIN2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # A72-30G

Lot # C2030-5

Product Description

Full-length recombinant human AXIN2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC101533](#).

Gene Aliases

AXIL

Formulation

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

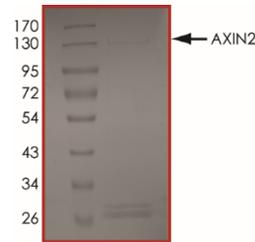
Scientific Background

AXIN2 plays an important role in the regulation of the stability of beta-catenin in the Wnt signaling pathway and Axin2-expressing cells themselves produce Wnt signals as well as long-range secreted Wnt inhibitors, suggesting an autocrine mechanism of stem cell self-renewal (1). AXIN2 is highly expressed in thymus, prostate, testis, ovary, and small intestine. The AXIN2 gene has been mapped to 17q23-q24, a region that shows frequent loss of heterozygosity in breast cancer, neuroblastoma, and other tumors. AXIN2 as a gene that is responsible for hereditary colorectal cancer and tooth agenesis that may be an indicator of cancer susceptibility (2).

References

1. Lim, X.et.al: Interfollicular epidermal stem cells self-renew via autocrine Wnt signaling. Science 342: 1226-1230, 2013.
2. Lammi, L.et.al: Mutations in AXIN2 cause familial tooth agenesis and predispose to colorectal cancer. Am. J. Hum. Genet. 74: 1043-1050, 2004.

Purity



The purity of AXIN2 protein was determined to be **>70%** by densitometry. Approx. MW **130 kDa**.

AXIN2 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #

A72-30G

Lot #

C2030-5

Purity

>70%

Concentration

0.05 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.