

AXIN1 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog # A71-30G

Lot # G1474-1

Product Description

Full-length recombinant human AXIN1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC044648](#).

Gene Aliases

AXIN; MGC52315

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

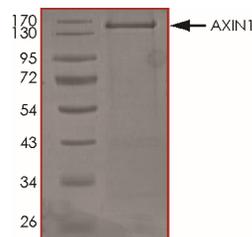
AXIN 1 encodes a cytoplasmic protein which contains a regulation of G-protein signaling (RGS) domain and a dishevelled and axin (DIX) domain that interacts with adenomatous polyposis coli, catenin beta-1, glycogen synthase kinase 3 beta, protein phosphate 2, and itself. AXIN1 has both positive and negative regulatory roles in Wnt-beta-catenin signaling. AXIN1 is a core component of a 'destruction complex' that promotes phosphorylation and polyubiquitination of cytoplasmic beta-catenin, resulting in beta-catenin proteasomal degradation in the absence of Wnt signaling. Nuclear accumulation of AXIN1 can positively influence beta-catenin-mediated transcription during Wnt signalling and can induce apoptosis (1). AXIN1 may be an effective therapeutic molecule for suppressing growth of hepatocellular and colorectal cancers (2).

References

1. Lui, T. T. H.et.al: The ubiquitin-specific protease USP34 regulates axin stability and Wnt/beta-catenin signaling. *Molec. Cell. Biol.* 31: 2053-2065, 2011.
2. Satoh, S.et.al: AXIN1 mutations in hepatocellular carcinomas, and growth suppression in cancer cells by virus-mediated transfer of AXIN1. *Nature Genet.* 24: 245-250, 2000.

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

Purity



The purity of AXIN1 protein was determined to be **>75%** by densitometry.
Approx. MW **135 kDa**.

AXIN1 Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	A71-30G
Lot #	G1474-1
Purity	>75%
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.