

WWP1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # W296-30G

Lot # Y791-2

Product Description

Recombinant full-length human WWP1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_007013](#).

Gene Aliases

AIP5, Tiu1, hSDRP1

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.

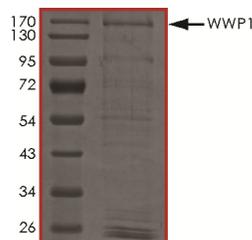
Scientific Background

WWP1 or WW domain containing E3 ubiquitin protein ligase 1 belongs to a family of NEDD4-like proteins, which are E3 ubiquitin-ligase molecules and regulate key trafficking decisions, including targeting of proteins to proteasomes or lysosomes (1). WWP1 are found in all eukaryotes and play an important role in the regulation of a wide variety of cellular functions such as protein degradation, transcription, and RNA splicing. WWP1 negatively regulates TGF-beta signaling by directing the degradation of active components of the TGF-beta network. Wwp1 and Ubc18 function to ubiquitinate substrates that regulate longevity induced by diet restriction (2).

References

1. Flasza, M.et.al: Alternative splicing determines the domain structure of WWP1, a Nedd4 family protein. Biochem. Biophys. Res. Commun. 290: 431-437, 2002.
2. Carrano, A.et.al: A conserved ubiquitination pathway determines longevity in response to diet restriction. Nature 460: 396-399, 2009.

Purity



The purity of WWP1 protein was determined to be **>70%** by densitometry. Approx. MW **160kDa**.

WWP1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog #

W296-30G

Lot #

Y791-2

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.