

Product Information

Description:	Recombinant human WWP1 (WW Domain Containing E3 Ubiquitin Protein Ligase 1), full length, encompassing amino acids 2-end. This construct contains an N-terminal FLAG-tag. This protein was affinity purified.
Background:	WW Domain Containing E3 Ubiquitin Protein Ligase 1 (WWP1) is a HECT-type E3 Ub ligase belonging to the NEDD4 family of E3 ligases. WWP1 interacts with and ubiquitinates many substrates, including transcription factor Δ Np63 as well as phosphatase and tensin homolog (PTEN), and regulates the degradation of sodium channels and membrane receptors. Genetic mutations or regulatory defects involving WWP1 are associated with neurological disorders and cancer. Aberrant expression of WWP1 in gastric, prostate, and breast cancer, for instance, is an area of high interest, and therefore, WWP1 represents an excellent candidate for therapeutic targeting in multiple cancer types.
Species	Human
Construct:	WWP1 (FLAG-2-922(end))
Concentration:	0.57 mg/ml
Expression System:	Sf9
Purity:	$\geq 90\%$
Format:	Aqueous buffer solution.
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 10% glycerol, 3 mM DTT, 100 μ g/ml FLAG peptide
MW:	106 kDa
Genbank Accession:	NM_007013
Stability:	At least 6 months at -80°C .
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

