

WWP2 (WW Domain Containing Protein 2)

Cat. # UB316H

Background

The E3 ubiquitin ligase WW domain containing protein 2 (WWP2) plays a crucial role in regulating protein degradation within cells. WWP2 contains WW domains, which interact with specific protein motifs, and a HECT (Homologous to E6-AP Carboxyl Terminus) domain responsible for transferring ubiquitin molecules onto target proteins. WWP2 has been implicated in various cellular processes, including the regulation of protein stability, cell growth, and signal transduction pathways. It is particularly known for its role in ubiquitinating and targeting specific substrates for degradation, thereby influencing cellular homeostasis and physiological responses.

Alternate Names

AIP2, Atrophin-1-Interacting Protein 2, HECT-Type E3 Ubiquitin Transferase WWP2

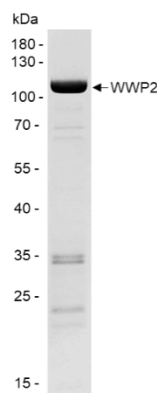
Application(s)

Ubiquitin ligation reactions

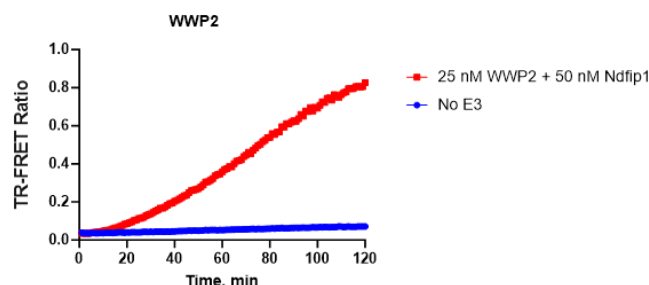
Product Specifications

Tag	His ₆
Purity	≥ 90% by SDS-PAGE
Molecular Weight	112 kDa
Quantity	25 µg
Species	Human
Expression System	<i>E. coli</i>
Physical State	Liquid
Buffer	50 mM Tris-HCl pH 8.0, 150 mM NaCl, 10% glycerol, 1 mM DTT
Solubility	> 3 mg/mL
Activity	A typical enzyme concentration of 10-200 nM is used for in vitro conjugation, depending on assay conditions.
Stability & Storage	Over 1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page analysis of purified WWP2. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of WWP2. 25 nM WWP2 + 50 nM Ndfip1 were tested in a TR-FRET assay for 120 minutes and showed a robust Signal to Background ratio.

References

1. Yang Y., et al., Proc Natl Acad Sci U S A. 2013;110(13):5115-20.
2. Chantray A. Cell Cycle. 2011;10(15):2437-9.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved