

WWP2 (WW Domain Containing Protein 2)

Cat. # UB316H

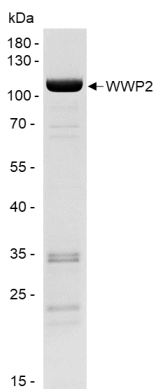
Background

WWP2, or WW domain containing E3 ubiquitin protein ligase 2, is a ubiquitin ligase enzyme that 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.

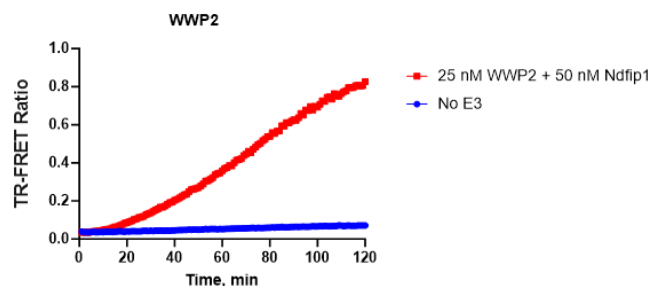
Product Information

Purity	≥ 90% by SDS-PAGE
Molecular Weight	112 kDa
Quantity	25 µg
Physical State	Liquid, 50 mM Tris-HCl pH 8.0, 150 mM NaCl, 10% glycerol, 1 mM DTT
Species	Human
Source	<i>E. coli</i>
Tag	His6
Activity	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
Storage	-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 was tested in a TR-FRET assay for 120 minutes and shows robust S/B.

References

1. Zhang, N., et al., Cell Death Differ., 2020. 27(9):2605-2619.
2. Zhu W, et al., J Biol Chem. 2017;292(27):11178-88.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic use.

Copyright © 2009 LifeSensors, Inc. All Rights Reserved