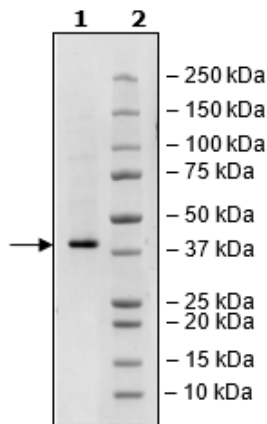


## Product Information

<b>Construct:</b>	WDR5 (Avi-His-2-334(end))-(Biotin)
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
<b>Concentration:</b>	0.48 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 mM Imidazole, and 20% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_017588
<b>MW:</b>	39 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

