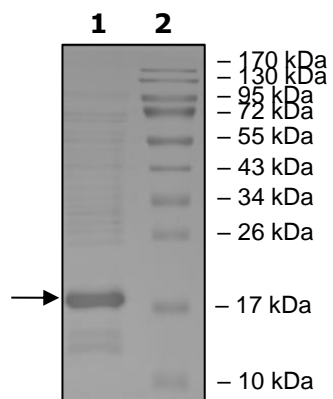


## Product Information

<b>Description:</b>	Recombinant full length human UBE2A (ubiquitin-conjugating enzyme E2). This construct contains an N-terminal His-tag. The recombinant protein was affinity purified and is active.
<b>Species:</b>	Human
<b>Construct:</b>	UBE2A (His-Full Length)
<b>Concentration:</b>	0.10 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM Imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol
<b>MW:</b>	17 kDa
<b>Genbank Accession:</b>	NM_003336
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Specific Activity:</b>	30 pmol/min/μg
<b>Assay Conditions:</b>	<p>UBE2A activity was measured by using the wild type ubiquitin protein diluted in Ubiquitination Buffer [(40 mM Tris pH 7.5), 20 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA and 0.5 mM DTT]) to a working concentration of 170 ng/μl (20 μM), in AMP-Glo™ based assay (Promega #V5011). Reaction was initiated by mixing 5 μl of a solution of 2x the final concentrations of UBE2A with 5 μl of a 2x Reaction Cocktail (1 μl of stock solution of ubiquitin, 15 ng/μl UBA1 + + 50 μM ATP).</p> <p>The reaction was initiated by mixing 5 μl of 2x UBE2A protein dilutions with 5 μl of 2x Reaction Cocktail for 60 minutes at 37°C. After the plate was equilibrated to Room Temperature (RT), 10 μl of AMP-Glo™ Reagent I was added, mixed for 1-2 minutes and incubated at RT for 60 minutes. AMP-Glo™ Reagent II was diluted 100-fold with Kinase-Glo™ One Solution, and 20 μl were added to each well, mixed for 1-2 minutes and incubated at RT for 30 minutes. The plate was read in a GloMax plate reader (Promega # E7031). The Enzyme Specific Activity was calculated as follows: [ (AMP concentration in μM) * (Reaction Volume in μl)] / Reaction time in min) *(Enzyme amount in mg)] x 10<sup>-3</sup>. A blank can be determined from a “no substrate” sample by replacing the enzyme solution with an equal volume of Ubiquitination Buffer.</p>
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining



### Specific Activity

