

## Product Information

<b>Construct:</b>	Tim1 (21-293-Avi-His)
<b>Concentration:</b>	1.37 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	HEK293
<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_012206
<b>MW:</b>	32 kDa
<b>Purity:</b>	≥90%
<b>Notes:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

## Quality Control Data

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

