

## Product Information

<b>Construct:</b>	ACHE (32-614(end)-His)
<b>Concentration:</b>	0.16 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 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_000665
<b>MW:</b>	66 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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

