

## Product Information

<b>Construct:</b>	CBL-B (His-Avi-39-426)
<b>Concentration:</b>	1.04 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM HEPES, pH 7.4, 120 mM NaCl, and 20% glycerol.
<b>Expression System:</b>	<i>E. coli</i>
<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_170662
<b>MW:</b>	48 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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

