

## Product Information

<b>Construct:</b>	CLPP(His-15-207(end))
<b>Concentration:</b>	1.58 mg/ml
<b>Species:</b>	<i>E. coli</i>
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 200 mM imidazole
<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>	NP_414971
<b>MW:</b>	23 kDa
<b>Purity:</b>	80%

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining

