

## Product Information

<b>Construct:</b>	MSP1D (His-TEV-79-267(end))
<b>Concentration:</b>	14.79 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	20 mM Tris-HCl, pH 7.4, 100 mM NaCl, 0.5 mM EDTA
<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_000039
<b>MW:</b>	25 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining

