

## Product Information

<b>Construct:</b>	Sestrin 2 (FLAG-2-480(end))
<b>Concentration:</b>	0.70 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 100 µg/ml FLAG peptide, and 10% glycerol.
<b>Expression System:</b>	Sf9
<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_031459
<b>MW:</b>	55 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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

