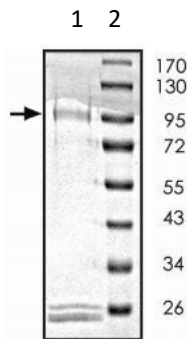


Product Information

<b>Construct:</b>	ULK2 (GST-1-478)
<b>Concentration:</b>	0.1 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris-HCL, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 25% 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_014683
<b>MW:</b>	98 kDa
<b>Purity:</b>	80%
<b>Specific Activity</b>	102 pmol/min/μg

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

