

## Product Information

<b>Construct:</b>	ERK4 (GST-Full length)
<b>Concentration:</b>	0.05 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 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_002747
<b>MW:</b>	110 kDa
<b>Purity:</b>	85%

## Quality Control Data

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

