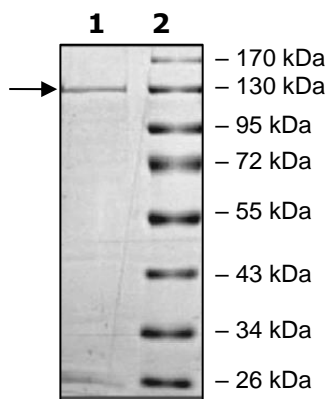


## Product Information

<b>Description:</b>	Recombinant full-length human EEF2K was expressed in Sf9 insect cells using an N-terminal GST-tag.
<b>Species:</b>	Human
<b>Construct:</b>	EEF2K (GST-Full Length)
<b>Concentration:</b>	0.10 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥70%
<b>Format:</b>	Aqueous buffer solution.
<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>MW:</b>	125 kDa
<b>Genbank Accession:</b>	NM_013302
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Specific Activity:</b>	42 pmol/min/μg
<b>Assay Conditions:</b>	EEF2K activity was measured by using the Ef2tide synthetic peptide substrate (RKKFGESEKTKTEFL), diluted in distilled water to a final concentration of 1 mg/ml, in a [33P]-ATP based assay. Reaction was initiated by mixing increasing amounts of the EEF2K with 1250 pmoles of [33P]-ATP in 25 mM MOPS, pH 7.2, 12.5 mM β-glycerol-phosphate, 25 mM MgCl <sub>2</sub> , 5 mM EGTA, 2 mM EDTA prepared with 250 μM DTT, co-factor calcium/calmodulin and substrate at a final concentration of 250 μg/ml. The reaction was initiated by addition of [33P]-ATP Assay Cocktail, followed by a 15-minute incubation at 30°C. The reaction was terminated by spotting the reaction mixture on phosphocellulose P81 paper, followed by three 10-minute washes with 1% phosphoric acid solution. Radioactivity was measured in a scintillation counter. The corrected activity (RLU) was calculated by removing the blank value for each sample. The Kinase Specific Activity was calculated as follows: $RLU / [(specific\ activity\ of\ [33P]-ATP\ in\ cpm/pmol) * (Reaction\ time\ in\ min) * (Enzyme\ amount\ in\ \mu g\ or\ mg)] * [(Reaction\ Volume) / (Spot\ Volume)]$ . The blank was determined from a “no substrate” sample by replacing the substrate solution with an equal volume of distilled water.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

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



### Specific Activity

