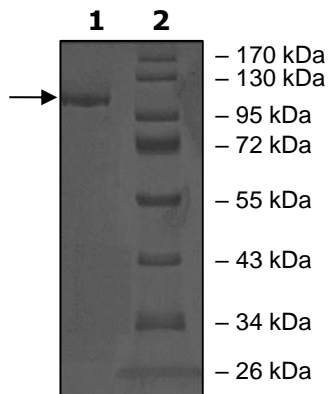


## Product Information

<b>Description:</b>	Recombinant full-length human RSK1 was expressed in Sf9 insect cells using an N-terminal GST-tag.
<b>Species:</b>	Human
<b>Construct:</b>	RSK1 (GST-Full Length)
<b>Concentration:</b>	0.10 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥90%
<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>	108 kDa
<b>Genbank Accession:</b>	NM_002953
<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>	210 pmol/min/μg
<b>Assay Conditions:</b>	<p>Kinase activity was measured using an S6K synthetic peptide substrate (KRRRLASLR) diluted in distilled water at 1.0 mg/ml.</p> <p>The protein kinase was diluted in a buffer containing 5 mM MOPS, pH 7.2, 2.5 mM β-glycero-phosphate, 5 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.4 mM EDTA, 50 ng/μl BSA (Bovine Serum Albumin) and 0.05 mM fresh DTT.</p> <p>Samples were prepared with 10 μl of diluted protein, 5 μl of substrate (1 mg/ml), and 5 μl of distilled water. The blank was determined from a “no substrate” sample. The reaction was initiated by addition of 5 μl of an ATP mix containing 1 mCi [33P]-ATP and 0.25 mM ATP, prepared in 6 ml of buffer: 25 mM MOPS, pH 7.2, 12.5 mM β-glycero-phosphate, 25 mM MgCl<sub>2</sub>, 5 mM EGTA, 2 mM EDTA and 0.25 mM fresh DTT.</p> <p>Samples were incubated for 15 minutes at 30°C, and the reaction was terminated by spotting 20 μl of the samples onto P81 phosphocellulose paper strips that were fixed in 1% phosphoric acid and washed three times. Radioactivity was determined using a scintillation counter.</p>
<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

