

## Product Information

<b>Construct:</b>	NUAK2 (K81R) (GST-Full Length)
<b>Mutation:</b>	K81R
<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_030952
<b>MW:</b>	110 kDa
<b>Purity:</b>	70%
<b>Assay Conditions:</b>	Kinase activity was measured using a CHKtide peptide substrate (KKKVSRSGLYRSPSPENLNRPR) diluted in distilled water to a final concentration of 1 mg/ml. Increasing amounts of kinase were mixed with CHKtide peptide substrate with a final concentration of 200 µg/ml in a buffer containing 5 mM MOPS, pH 7.2, 2.5 mM β-glycerol-phosphate, 5 mM MgCl <sub>2</sub> , 1 mM EGTA, 0.4 mM EDTA and 0.05 mM fresh DTT to a final volume of 20 µl. The reaction was initiated by addition of 5 µl of [33P]-ATP diluted in kinase buffer: 6 ml kinase buffer containing 1 mCi [33P]-ATP, 0.25 mM ATP, 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. After incubating for 30°C for 15 minutes, the reaction was terminated by spotting 20 µl of the mixture onto phosphocellulose paper strips that were fixed in 1% phosphoric acid and washed three times. Radioactivity was determined using a scintillation counter. The blank was determined from a “no substrate” sample.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

