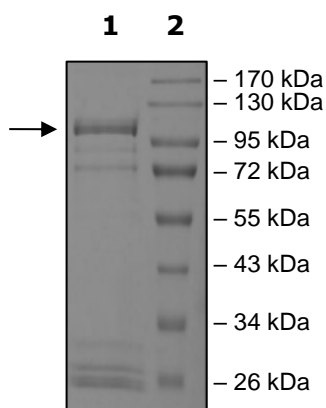


Product Information

Construct:	NUAK2 (GST-Full Length)
Concentration:	0.10 mg/ml
Species:	Human
Formulated In:	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.
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_030952
MW:	110 kDa
Purity:	70%
Specific Activity:	10 pmol/min/μg
Assay Conditions:	Kinase activity was measured using ADP-Glo™ Kinase Assay Kit (Promega; Cat# V9101) which quantifies the amount of ADP produced by the NUAK2 reaction. The ADP-Glo™ Reagent is added to terminate the reaction and quench the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly converted ATP using a luciferase/luciferin reaction. Assay: Kinase activity was measured using substrate CHKtide (KKKVSRSGLYRSPSPENLNRPR) (diluted in distilled water to a final concentration of 1 mg/ml. Increasing amounts of NUAK2 kinase were incubated with a final concentration of 0.2 mg/ml substrate in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl ₂ , 0.1 mg/ml BSA, and 50 μM fresh DTT. The reaction was initiated by adding a final concentration of 25 μM ATP and incubating for 40 minutes at room temperature. The reaction was terminated by the addition of 5 μl of ADP-Glo™ Reagent and a subsequent incubation at room temperature for 40 minutes. Luminescence was measured by the addition of 10 μl Kinase Detection Reagent followed by incubation at room temperature for an additional 30 minutes. The blank was determined from a “no kinase” sample.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

