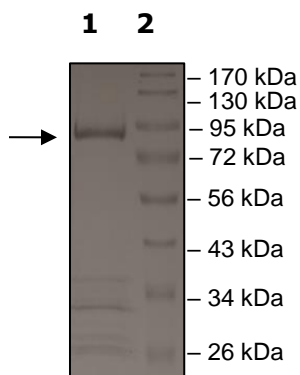


Product Information

Construct:	WNK4 (GST-1-441)
Concentration:	0.10 mg/ml
Species:	Mouse
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_175638
MW:	86 kDa
Purity:	85%
Specific Activity:	1.5 pmol/min/μg
Assay Conditions:	10 μl of varying amounts of WNK4 was mixed with 10 μl of 0.5 mg/ml PKA protein substrate. Reaction was initiated by adding 5 μl of 250 μM [33P]-ATP and incubating for 15 min at 30°C followed by spotting the reaction on phosphocellulose paper, fixing in 1% phosphoric acid, and assaying with a scintillation counter.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

