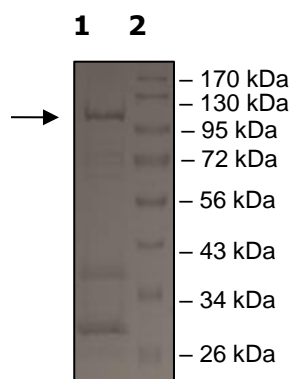


Product Information

| | |
|---------------------------|---|
| Construct: | WNK2 (GST-1-489) |
| 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_006648 |
| MW: | 100 kDa |
| Purity: | 70% |
| Specific Activity: | 2.0 pmol/min/μg |
| Assay Conditions: | 10 μl of varying amounts of WNK2 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

