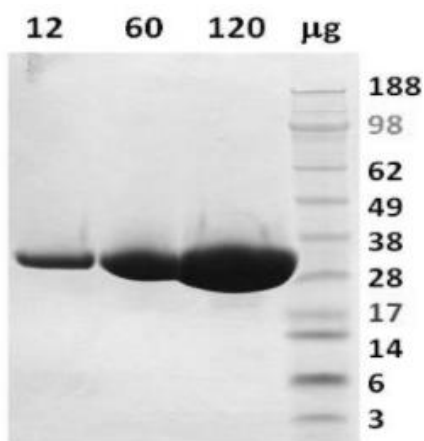


Product Information

Construct:	PAK5 (425-719)
Concentration:	1.00 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 150 mM NaCl, 10% glycerol, 5 mM DTT.
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_177990
MW:	34 kDa
Specific Activity:	4,199 pmole/min/μg
Assay Conditions:	Analysis of enzymatic activity was performed according to the Zlyte assay protocol (Invitrogen): Different concentrations of PAK5 were incubated in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl ₂ , 1 mM EGTA, 200 mM ATP, 0.01% Brij-35, and 2 μM substrate (SER/THR 14, Invitrogen) at room temperature for 1 hour. Developer solution was added to the reaction and the reaction was stopped after 1 hour of incubation at room temperature. Fluorescence was detected using $\lambda_{exc}=460\pm40$ nm and $\lambda_{em}=528\pm20$ nm filters.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

SDS-Page Analysis



Specific Activity

