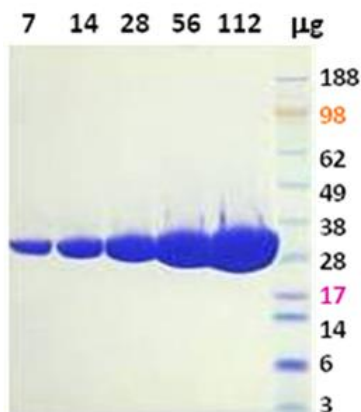


Product Information

Description:	PAK4 is a recombinant wild type human PAK4 kinase domain of full-length human PAK4 expressed in <i>E. coli</i> and purified to homogeneity with full activity.
Construct:	PAK4 (300-591)
Concentration:	1.0 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
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Genbank Accession:	NM_005884
MW:	33.3 kDa
Purity:	≥95%
Specific Activity:	5,595 pmol/min/μg
Assay Conditions:	Analysis of enzymatic activity was performed according to the Zlyte assay protocol (Invitrogen): 1. Different concentrations of PAK4 were incubated in a buffer containing 50 mM HEPES pH 7.5, 10 mM MgCl ₂ , 1 mM EGTA, 200 μM ATP, 0.01% Brij-35, and 2 μM substrate (SER/THR 14, Invitrogen) at room temperature for 1 hour. 2. Developer solution was added to the reaction and the reaction was stopped after 1 hour of incubation at room temperature. 3. Fluorescence was then detected using λ _{exc} =460±40 nm and λ _{em} =528±20 nm filters.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

Coomassie Staining



Specific Activity

