

Data Sheet

PAK6

Human, recombinant

Catalog #: 40081

Lot#: 110520

Conc.: 1 mg/ml

Formulated in: 25 mM Tris-HCl, pH8, 150 mM NaCl, 10% glycerol, 5 mM DTT.

Stability: >6 months at -80°C

References:

1. Kaur, R., et al., Prostate 68 (14), 1510-1516 (2008)
2. Hosgood, H.D. et al., Carcinogenesis 29 (10), 1938-1943 (2008)

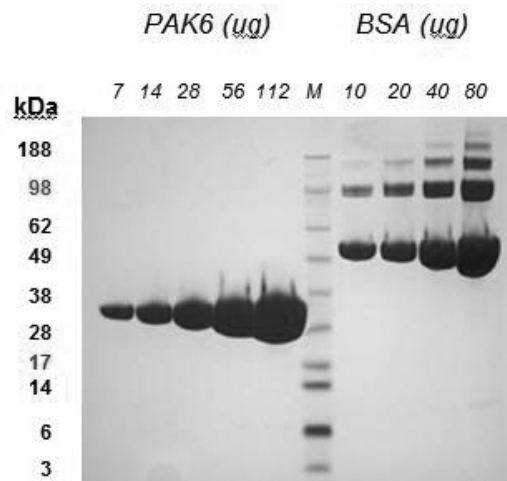
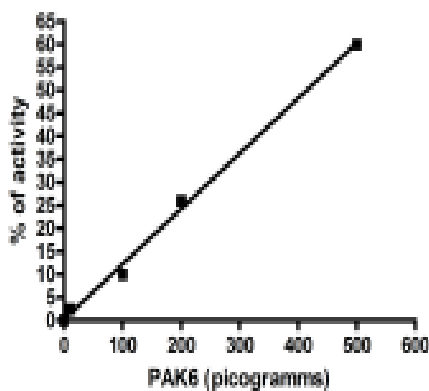
Description: Human PAK6 (GenBank Accession No. NM_020168), amino acids 385-680, MW=34.2 kDa, expressed in E.coli.

Specific Activity: 2.1 nmoles/min/μg. Analysis of enzymatic activity was performed according to Zlyte assay protocol (Invitrogen): 1. Different concentration of PAK6 were incubated in a buffer containing 50 mM HEPES (pH7.5), 10 mM MgCl₂, 1 mM EGTA, 200 μM ATP, 0.01% Brij-35, and 2 μM substrate (SER/THR 20, Invitrogen) at room temperature for 1 hour.

2. Developer solution was added to reaction and reaction was stopped after 1h of incubation at RT. 3. Fluorescence was detected using λ_{exc}=460±40 nm and λ_{em}=528±20 nm filters.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com