

Data Sheet

PAK1

human, recombinant

Catalog #: 40072

Lot#110413

Conc. 1 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0,
 150 mM NaCl, 10% glycerol, and 5 mM DTT.

Stability: >6 months at -80°C

References:

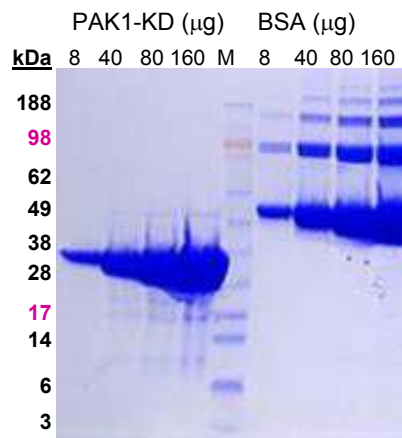
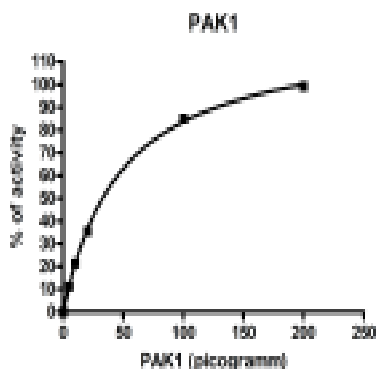
1. Sheehan, K.A., et al., Am. J. Physiol., Cell Physiol. 296 (1), C47-C58 (2009)
2. Reddy, S.D., et al., Cancer Res. 68 (20), 8195-8200 (2008)

Description: Human PAK1 (GenBank Accession No. NM_002576.4), amino acids 248-545, MW=33.6 kDa, expressed in E.coli.

Specific Activity: 32638 pmoles/min/μg. Analysis of enzymatic activity was performed according to Zlyte assay protocol (Invitrogen): 1. Different concentration of PAK1 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 14, 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 then 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