

Datasheet: PAK2

Lot: SP001

Description: Human PAK2 Full Length c-DNA (GenBank accession No. NM_002577), amino acids M₁-R₅₃₄, N-terminally fused to GST

Theoretical MW_{Fusionprotein}: 87,938 Da

Expression: Baculovirus infected Sf9 cells

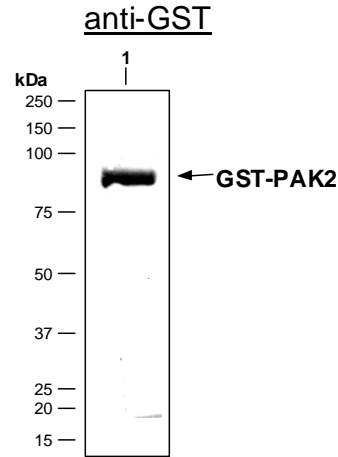
Purification: One-step affinity purification using GSH-agarose

Storage buffer: 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathion, 20% glycerol

Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

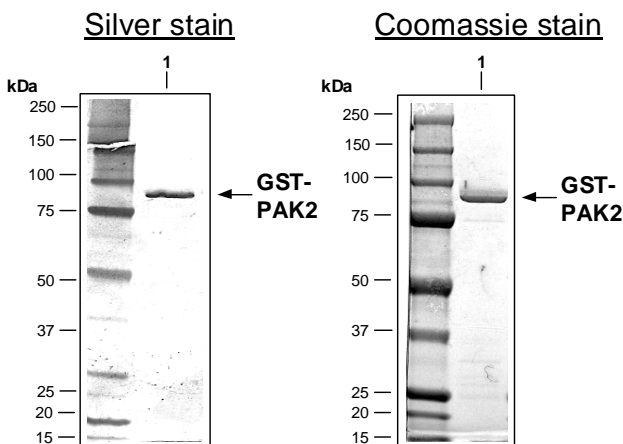
Protein concentration: 0.403 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Purified protein (Western blot):



1: SP001, 500 ng

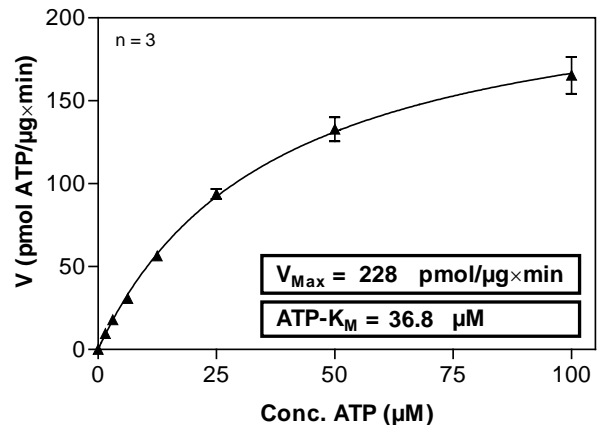
Purified protein (SDS-PAGE):



1: SP001, 1.0 µg

1: SP001, 2.0 µg

Kinase activity:



Final assay concentrations:

- 60 mM HEPES-NaOH, pH 7.5
- 3 mM MgCl₂
- 3 mM MnCl₂
- 3 µM Na-orthovanadate
- 1.2 mM DTT
- ATP (variable)
- 2.5 µg / 50 µl PEG_{20,000}
- Substrate: Tetra(LRRWSLG), 5 µg / 50 µl
- Recombinant PAK2: 50 ng / 50 µl

Assay: MAFC (Millipore) filter assay

Specific activity: 228 pmol/µg min