

# ARK5

aa1-661

**Product No.:** 0300-0000-1

**Lot:** 002

**Description:** Human ARK5

Amino acids M<sub>1</sub>-N<sub>661</sub> (as in GenBank entry NM\_014840)\*, N-terminally fused to GST-HIS<sub>6</sub>-Thrombin cleavage site

\*Sequence may contain documented polymorphisms  
Detailed sequence on request

**Product identity:** ARK5, Lot 002, was confirmed as human ARK5 by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

**Theoretical MW**<sub>Fusion Protein</sub>: 104,220 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, 4 mM reduced glutathione, 20% glycerol

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

**Protein concentration:** 0.176 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Method for determination of K<sub>m</sub> value & Specific activity:**

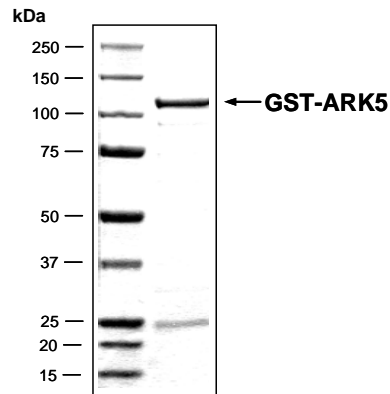
• Assay conditions:

60 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 µM Na-orthovanadate  
1.2 mM DTT  
50 µg / ml PEG<sub>20,000</sub>  
ATP (variable)  
Substrate: CHKtide-peptide (KKKVSRSGLYRSPSPENLNRPR), 50 µg / ml  
Recombinant ARK5: 4 µg / ml

• Filter binding assay  
MSPH membrane (Millipore)

**Specific activity:** 32 pmol/µg×min

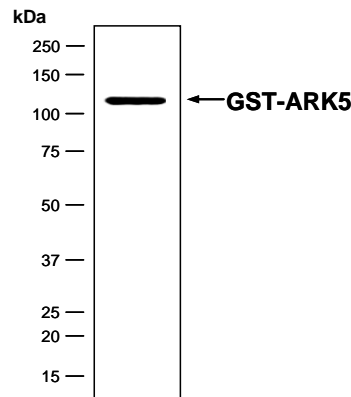
**Coomassie stain:**



2.0 µg GST-ARK5

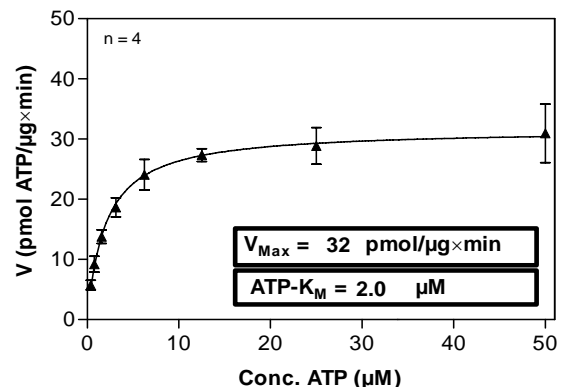
**Western blot analysis:**

anti-GST



500 ng GST-ARK5

**Determination of K<sub>m</sub> value for ATP:**



This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.