

Datasheet: IRAK4

Lot: SP001

Description: Human IRAK4 Fragment c-DNA (GenBank accession No. AF445802.1), amino acids A₁₀₄-S₄₆₀, N-terminally fused to GST-HIS₆-Thrombin cleavage site

Theoretical MW_{Fusionprotein}: 69,839

Expression: Baculovirus infected Sf9 cells

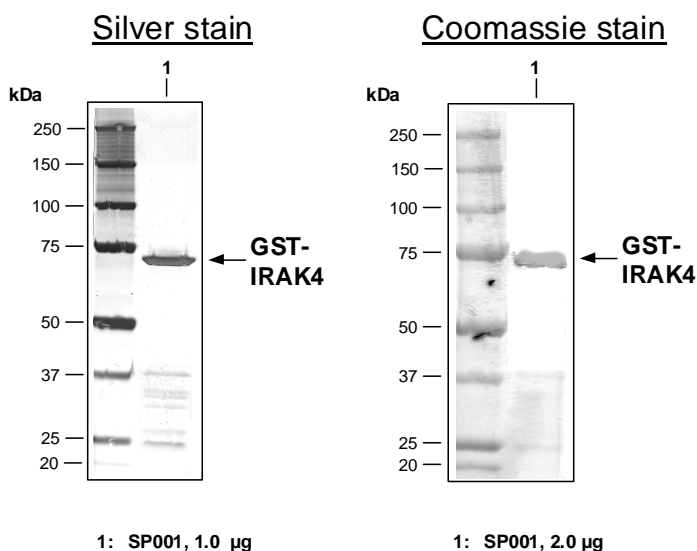
Purification: One-step affinity purification using GSH-agarose

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

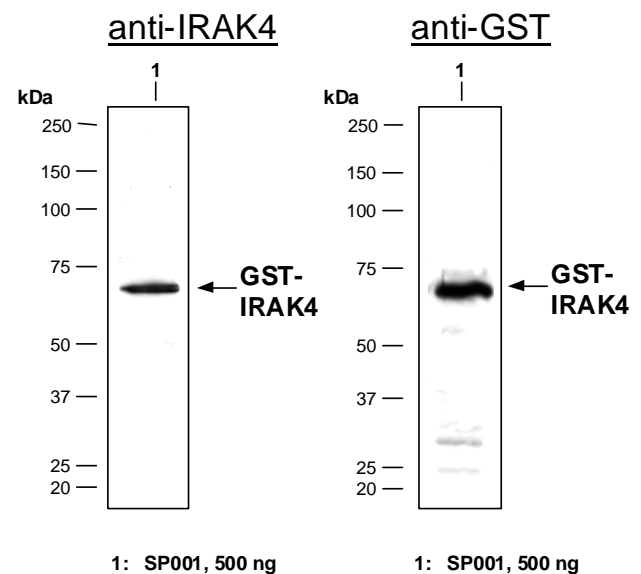
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.195 µg/µl (Bradford method using BSA Sigma A-7638, lot 79H7641 as standard protein)

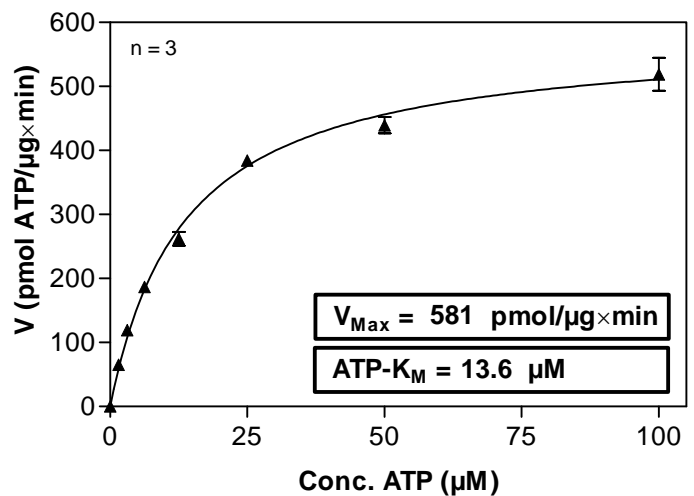
Purified protein (SDS-PAGE):



Purified protein (Western blot):



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: Histone H2B, 5 µg / 50 µl
- Recombinant IRAK4: 20 ng / 50 µl

Assay: MAFC (Millipore) filter assay

Specific activity: 581 pmol/µg min