

## Product Information

### JAK2(JH1 JH2) [V617F]

Product Number : 08-515

#### Product description

Human JAK2, catalytic/pseudo-kinase domain [532-1132(end) amino acids and V617F of accession number NP\_004963.1] was expressed as N-terminal GST-fusion protein (97 kDa) using baculovirus expression system. GST-JAK2(JH1 JH2)[V617F] was purified by using glutathione sepharose chromatography and gel filtration chromatography.

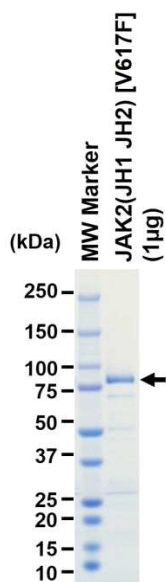
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.  
Avoid repeating freeze-thaws.

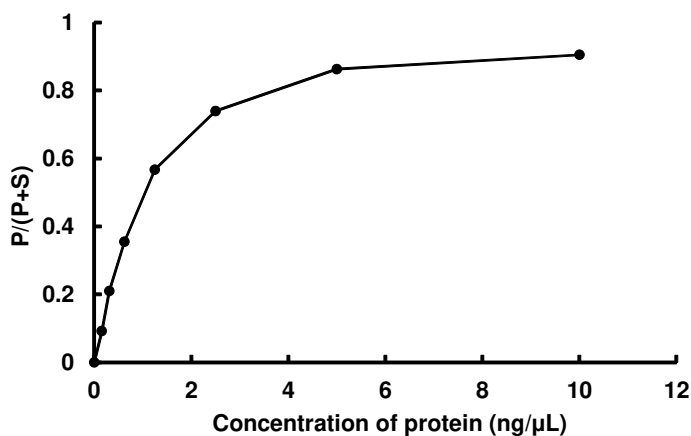
#### SDS-PAGE



Purity: 70 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Src tide

ATP : 100 µM