

## Product Information

### IRAK4

 Product Number : **09-145**

#### Product description

Full-length human IRAK4 [1-460(end) amino acids of accession number NP\_057207.2] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-IRAK4 was purified by using glutathione sepharose chromatography.

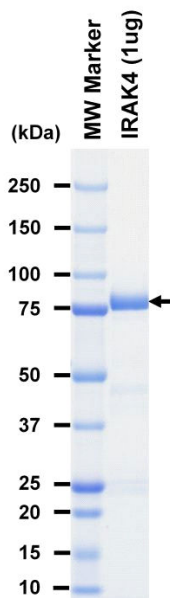
#### Storage buffer:

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

#### Storage and Handling:

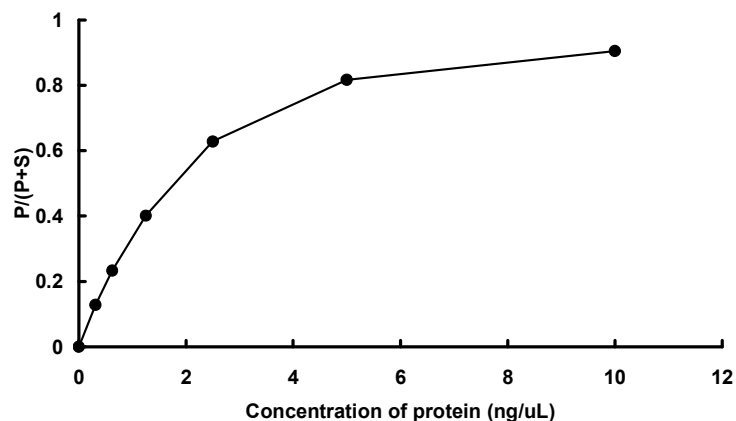
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 88 %  
 The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : IRAK1 peptide

ATP : 1000  $\mu$ M