

## Product Information

### DDR1(SRC treated)

 Product Number : **08-132**

#### Product description

Human DDR1, cytoplasmic domain [444-876(end) amino acids of accession number NP\_001945.3] was co-expressed as N-terminal GST-fusion protein (75 kDa) with human His-tagged SRC [1-536(end) amino acids of accession number NP\_005408.1] using baculovirus expression system. GST-DDR1 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

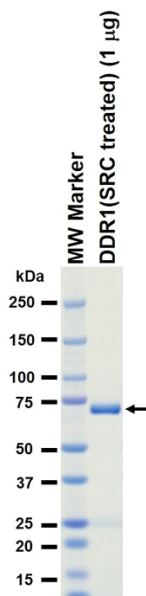
#### Storage buffer:

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.

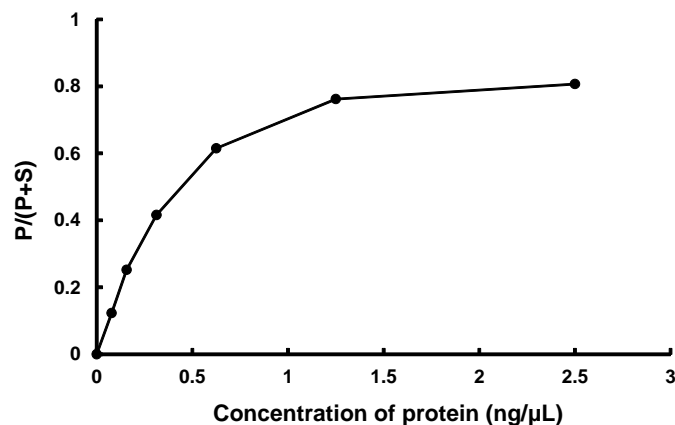
#### SDS-PAGE



Purity: 90 %

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 : IRS1

ATP : 100 μM