

## Product Information

### TYRO3

Product Number : **08-109**

#### Product description

Human TYRO3, cytoplasmic domain of [453-890(end) amino acids of accession number NP\_006284.2] was expressed as N-terminal GST fusion protein (76 kDa) using baculovirus expression system. GST-TYRO3 was purified by using glutathione sepharose chromatography

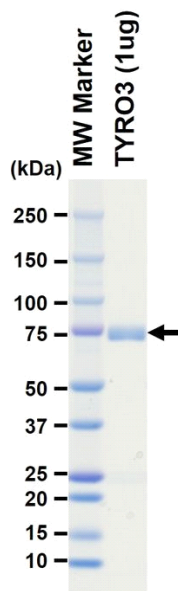
#### Storage buffer:

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

#### Storage and Handling:

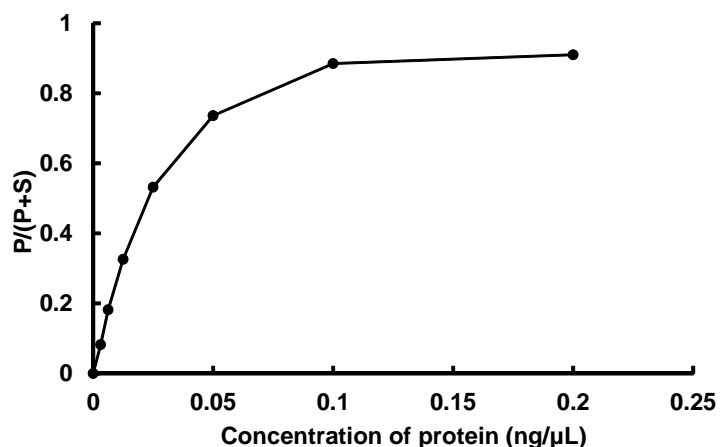
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 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 : CSKtide

ATP : 100 μM