

## Product Information

### FGFR3[G697C]

 Product Number : **08-542**

#### Product description

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and G697C of accession number NP\_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3[G697C] was purified by using glutathione sepharose chromatography.

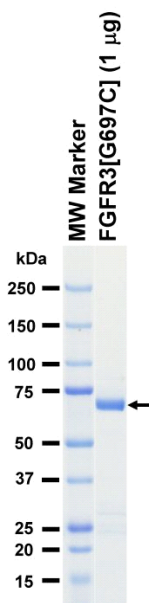
#### Storage buffer:

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.

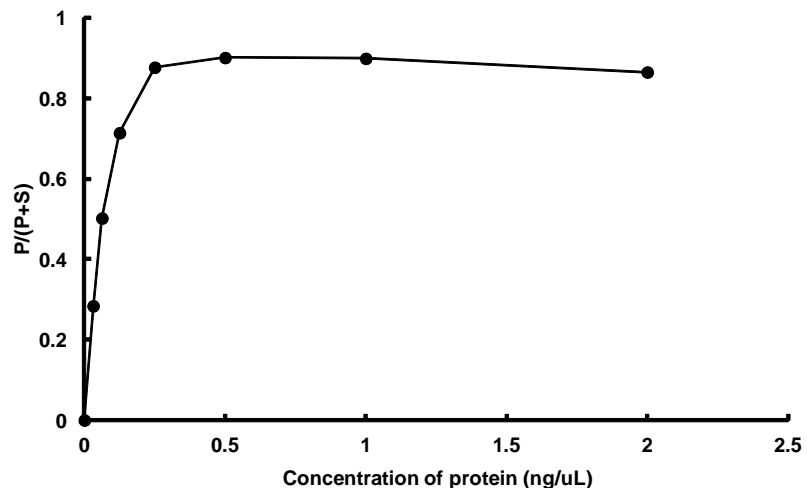
#### SDS-PAGE



Purity: 91 %

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 : CSKtide

 ATP : 100  $\mu$ M