

Product Information

FGFR1[V561M]

 Product Number : **08-536**

Product description

Human FGFR1, cytoplasmic domain [398-822(end) amino acids and V561M of accession number NP_075598.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR1[V561M] was purified by using glutathione sepharose chromatography.

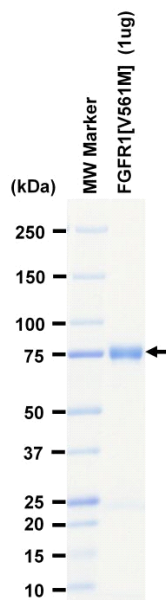
Storage buffer:

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

Storage and Handling:

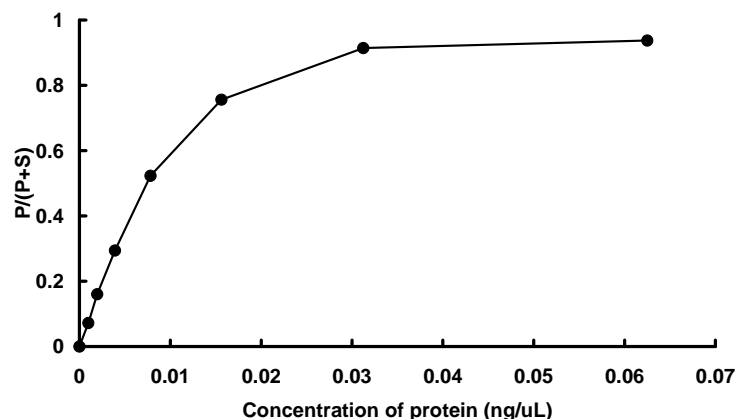
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 94 %
 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 μM