

Product Information

BTN-FGFR2[C491A][non-activated]

 Product Number : **08-903-23N**

Product description

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and C491A of accession number NP_000132.3] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (52 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

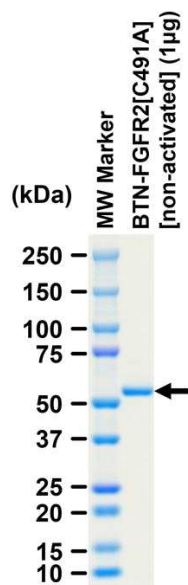
Storage buffer:

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.

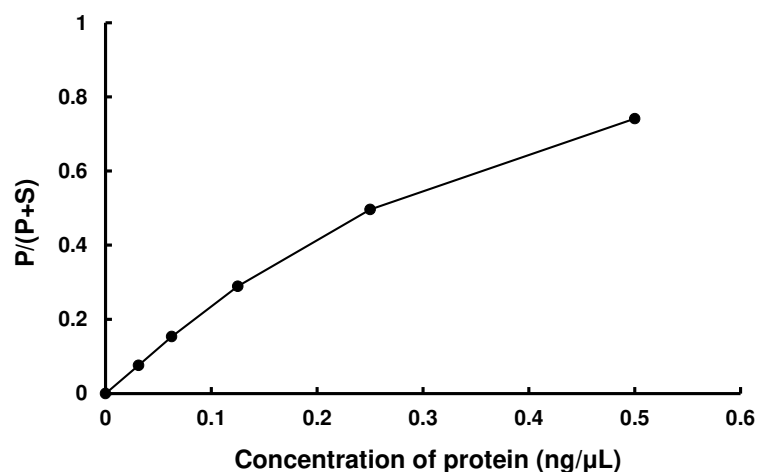
SDS-PAGE



Purity: 99 %

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