

Product Information

BTN-FGFR2[N549H]

 Product Number : **08-901-20N**

Product description

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and N549H 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 and activated with ATP. Activated protein was purified by using gel filtration.

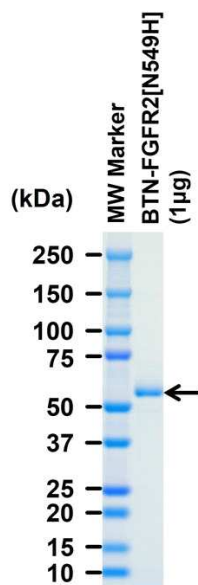
Storage buffer:

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.

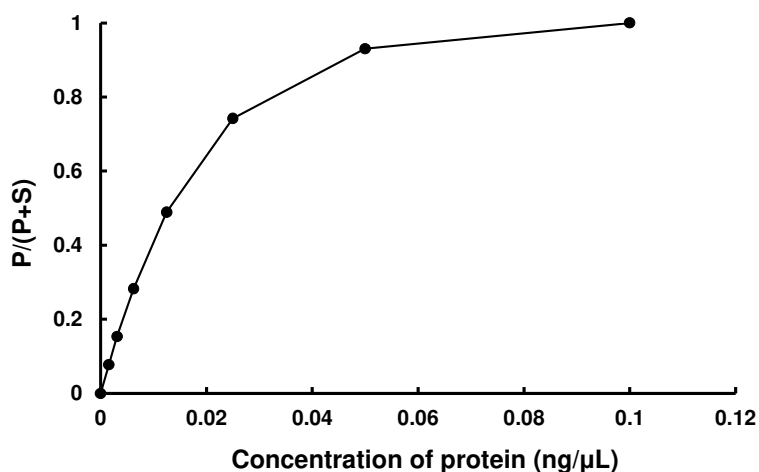
SDS-PAGE



Purity: 97 %

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