

Product Information

BTN-EGFR[d746-750/C797S]

 Product Number : **08-764-20N**

Product description

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and C797S of accession number NP_005219.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (65 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose

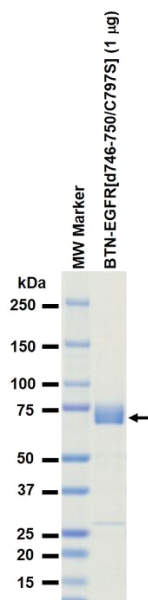
Storage buffer:

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.

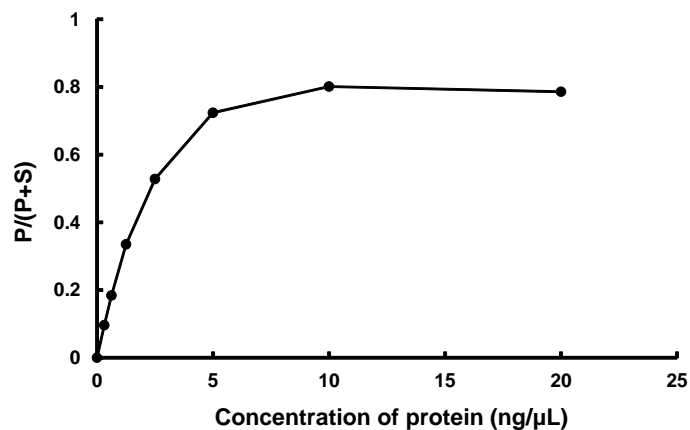
SDS-PAGE



Purity: 91 %

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 : Srctide

ATP : 100 µM