

Product Information

EGFR[d746-750/T790M/C797S]

 Product Number : **08-565**

Product description

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M/C797S of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa/T790M/C797S] was purified by using glutathione sepharose chromatography.

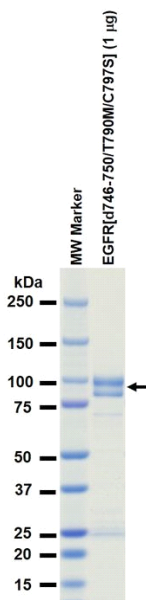
Storage buffer:

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.

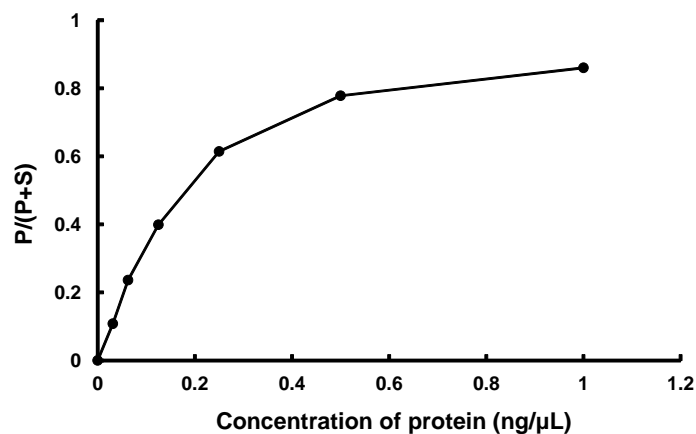
SDS-PAGE



Purity: 78 %

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 : Srcptide

ATP : 100 µM