

## Product Information

### EGFR [d746-750/T790M]

Product Number : **08-528**

#### Product description

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR [d746-750aa/T790M] was purified by using glutathione sepharose chromatography.

#### Storage buffer:

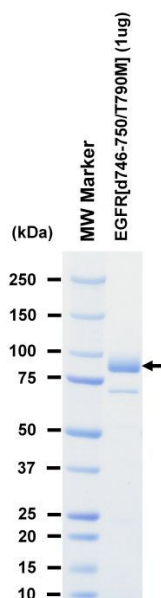
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

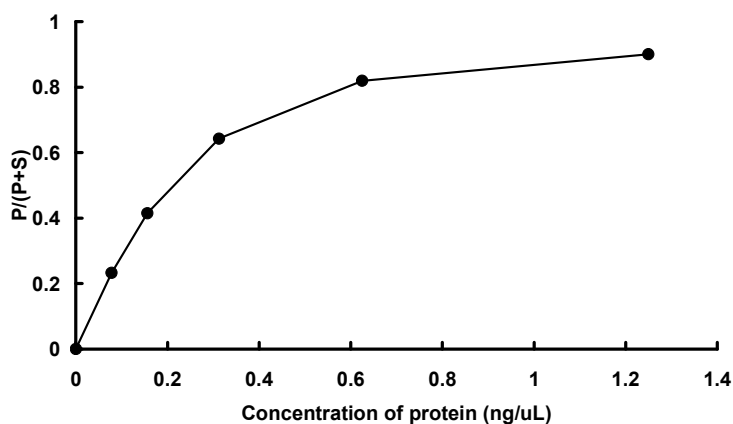
#### SDS-PAGE



Purity: 84 %

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 : Src tide

ATP : 100  $\mu$ M