

## Product Information

### EGFR[D770\_N771insNPG]

Product Number : 08-553

#### Product description

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and D770\_N771insNPG of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[D770\_N771insNPG] was purified by using glutathione sepharose chromatography.

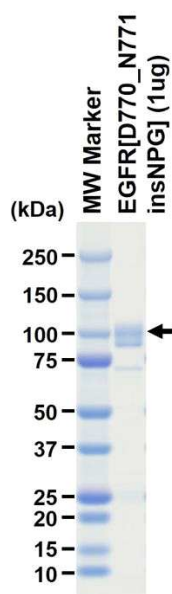
#### Storage buffer:

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.

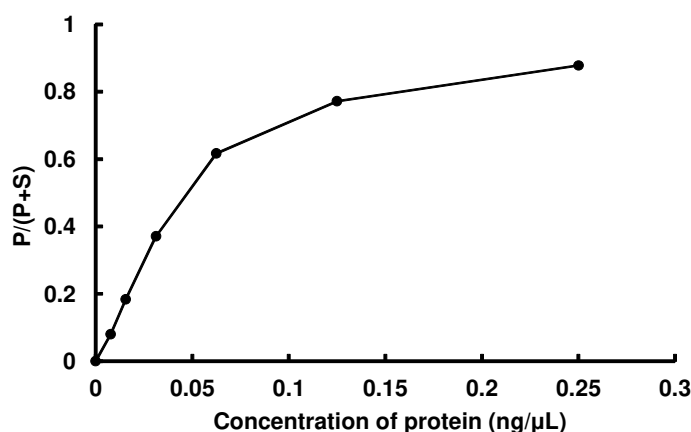
#### SDS-PAGE



Purity: 83 %

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 : Src tide

ATP : 100 μM

---

## Product Information

### EGFR[insNPG]

Product Number : **08-553**

#### Product description

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and D770\_N771insNPG of accession number NP\_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[D770\_N771insNPG] 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

#### Activity data

Data is not registered

Purity: 88 %

The purity was assessed  
by SDS-PAGE/CBB  
staining.

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 : Src tide

ATP : 100  $\mu$ M