

Product Information

EGFR[T790M/L858R]

Product Number : 08-510

Product description

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M, L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[T790M/L858R] was purified by using glutathione sepharose chromatography.

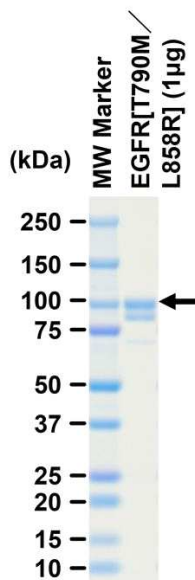
Storage buffer:

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

Storage and Handling:

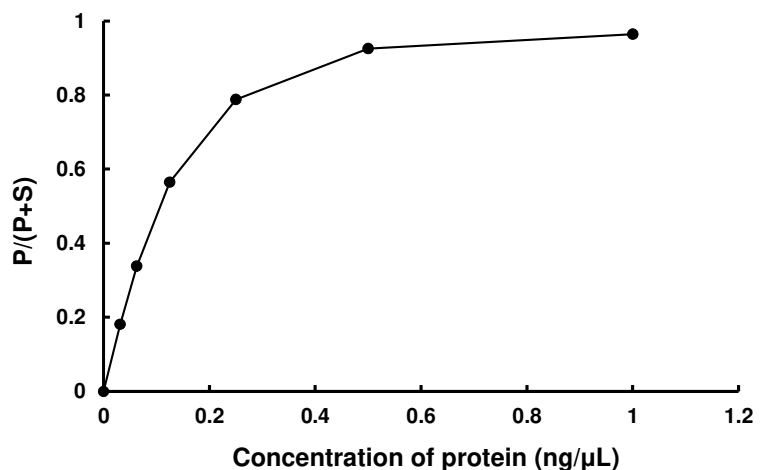
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 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