

Product Information

RET [G691S]

 Product Number : **08-522**

Product description

Human RET, cytoplasmic domain [658-1114(end) amino acids and G691S of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[G691S] was purified by using glutathione sepharose chromatography.

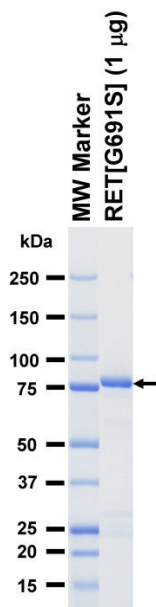
Storage buffer:

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

Storage and Handling:

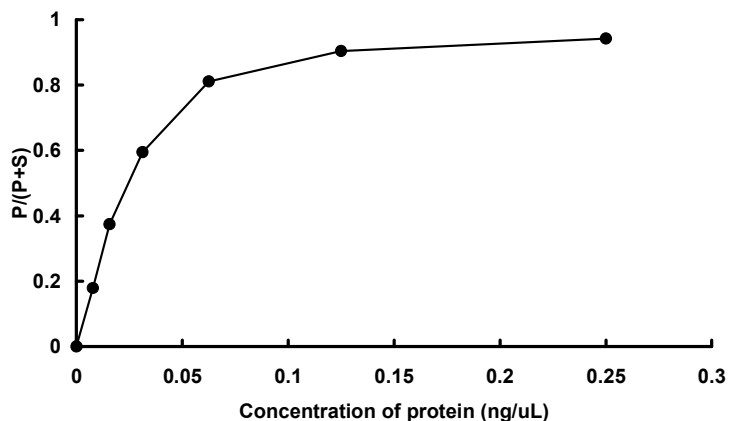
Store at -80C.
 Avoid repeating freeze-thaws.
 Use within 12 months from the date of shipment.

SDS-PAGE



Purity: 86 %
 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 : CSKtide

ATP : 100 µM