

Product Information

RET [S891A]

Product Number : **08-523**

Product description

Human RET, cytoplasmic domain [658-1114(end) amino acids and S891A of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[S891A] 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

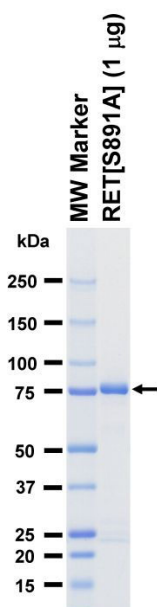
Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

Use within 12 months from the date of shipment.

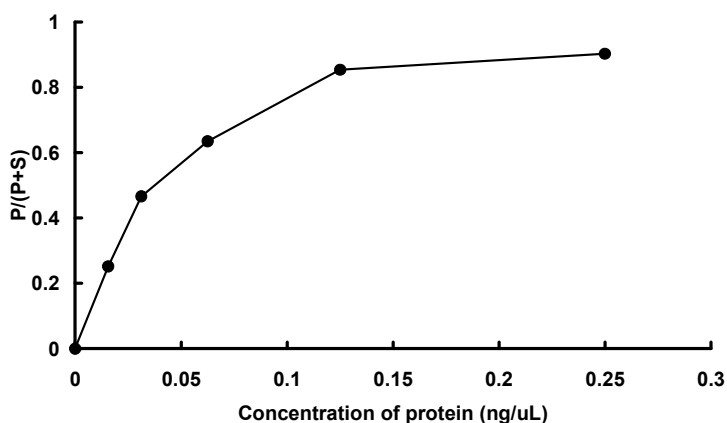
SDS-PAGE



Purity: 88 %

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