

Product Information

RET[V804M]

Product Number : **08-560**

Product description

Human RET, cytoplasmic domain [658-1114(end) amino acids and V804M of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[V804M] was purified by using glutathione sepharose chromatography.

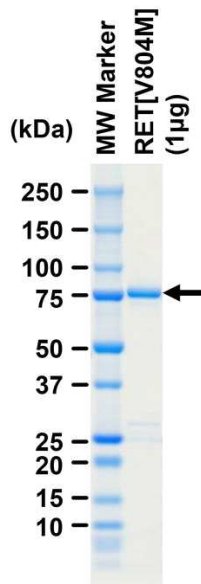
Storage buffer:

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

Storage and Handling:

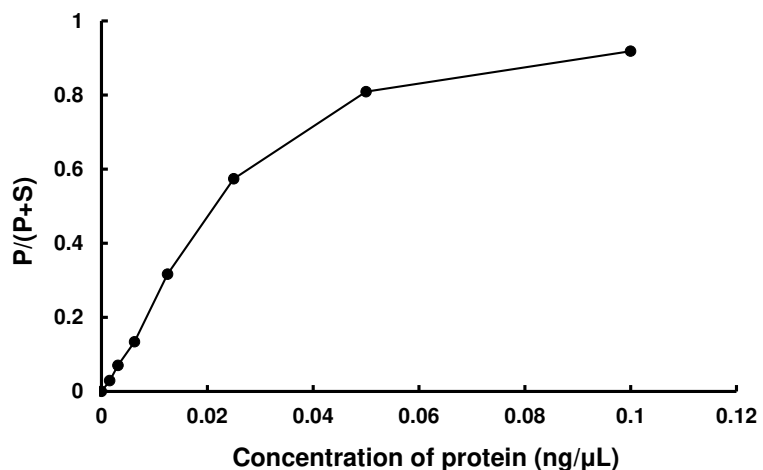
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 90 %
 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 : CSKtide

ATP : 100 μM