

Product Information

PDGFR α (PDGFRA) [D842V]

 Product Number : **08-506**

Product description

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and D842V of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFR α [D842V] was purified by using glutathione sepharose chromatography.

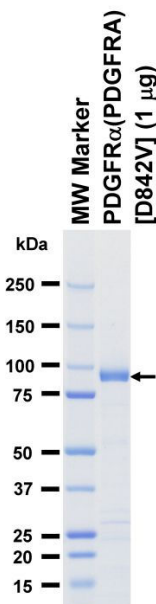
Storage buffer:

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.

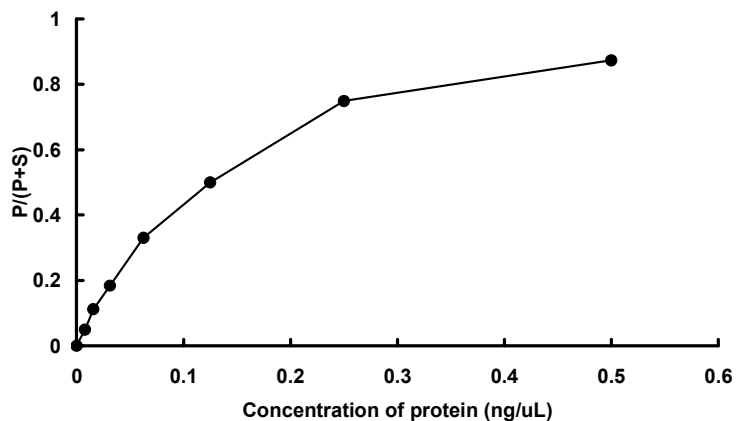
SDS-PAGE



Purity: 83 %

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

Activity data



The activity was measured by off-chip mobility shift assay. 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 μ M