

Product Information

KIT [D816E]

Product Number : **08-541**

Product description

Human KIT, cytoplasmic domain [544-976(end) amino acids and D816E of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[D816E] was purified by using glutathione sepharose chromatography.

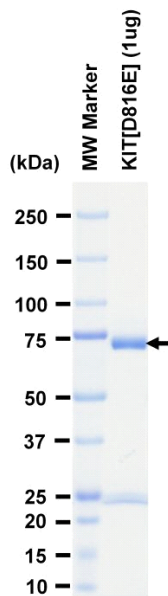
Storage buffer:

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

Storage and Handling:

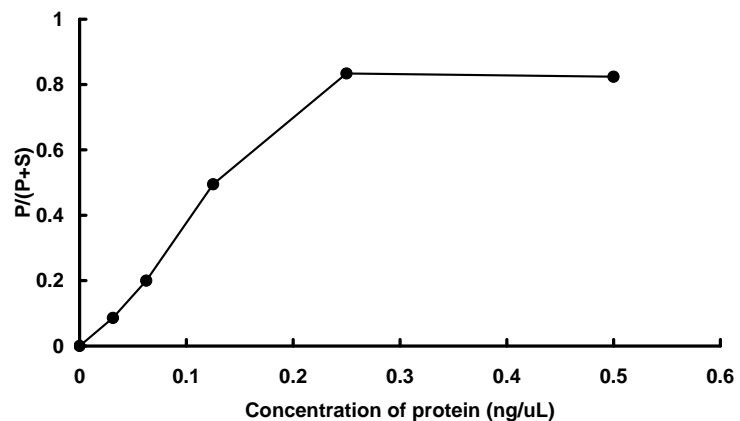
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 75 %
 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 : Srctide

ATP : 100 μ M