

Product Information

CLK2

Product Number : **04-127**

Product description

Full-length human CLK2 [1-499(end) amino acids of accession number AAH53603.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-CLK2 was purified by using glutathione sepharose chromatography.

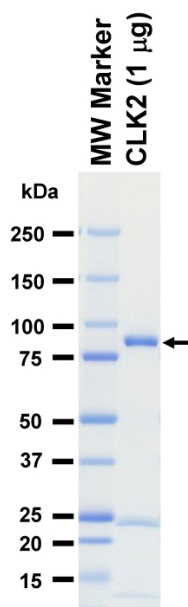
Storage buffer:

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

Storage and Handling:

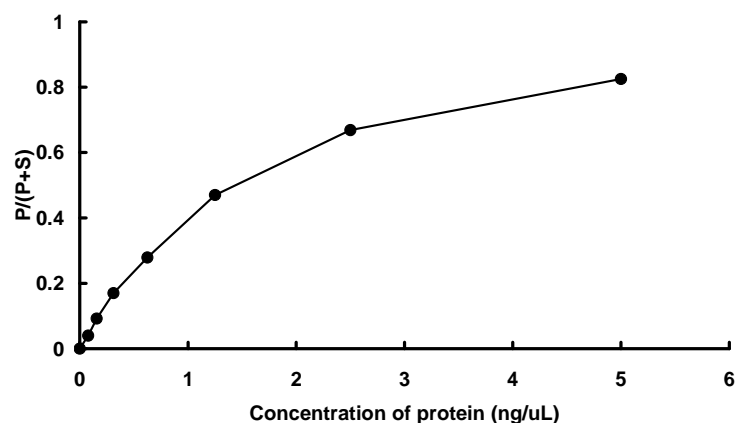
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 69 %
 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 : DYRKtide-F

ATP : 100 µM