

Product Information

CHAK1(TRPM7)

Product Number : 10-111

Product description

Human CHAK1, catalytic domain [1158-1865 amino acids of accession number NP_060142.3] was expressed as N-terminal GST-fusion protein (107 kDa) using baculovirus expression system. GST-CHAK1 was purified by using glutathione sepharose chromatography.

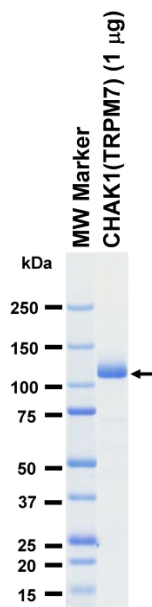
Storage buffer:

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.

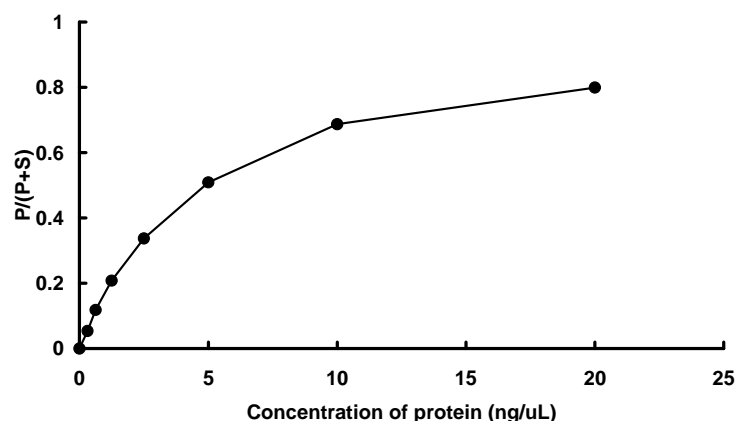
SDS-PAGE



Purity: 99 %

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 : EEF2Ktide

ATP : 1000 µM