

Product Information

ICK

Product Number : **04-156**

Product description

Full-length human ICK [1-632(end) amino acids of accession number NP_057597.2] was expressed as N-terminal GST-fusion protein (98 kDa) using baculovirus expression system. GST-ICK 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

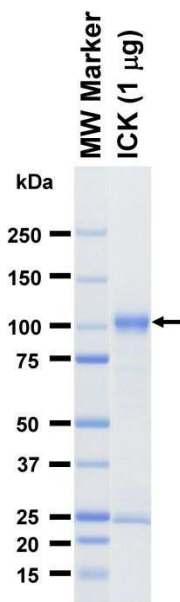
Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

Use within 12 months from the date of shipment.

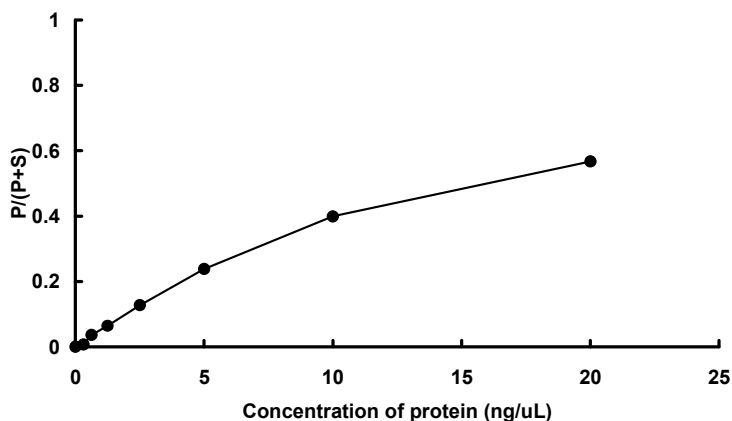
SDS-PAGE



Purity: 74 %

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 : Modified Erktide

ATP : 1000 µM