

Product Information

LTK[L650F]

 Product Number : **08-597**

Product description

Human LTK, catalytic domain [498-796 amino acids and L650F of accession number NP_002335.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-LTK[L650F] was purified by using glutathione sepharose chromatography.

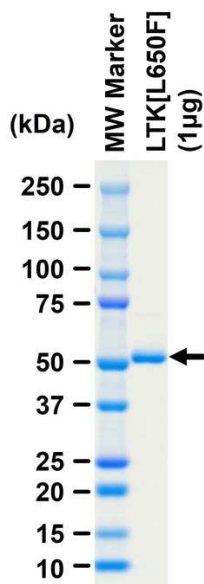
Storage buffer:

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

Storage and Handling:

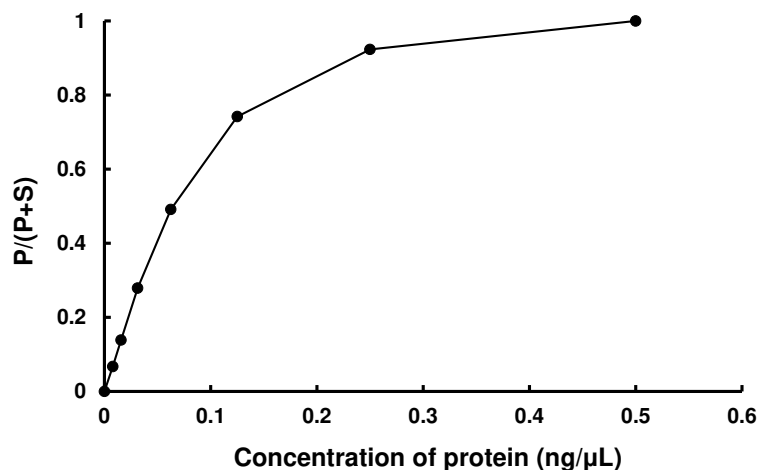
Store at -80C.
 Avoid repeating freeze-thaws.

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(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Srctide

ATP : 100 µM