

Product Information

BTN-BTK[C481S]

Product Number : **08-417-20N**

Product description

Full-length human BTK [2-659(end) amino acids and C481S of accession number NP_000052.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (80 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with ATP. Activated protein was purified by using gel filtration.

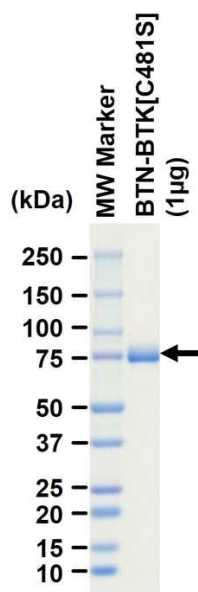
Storage buffer:

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.

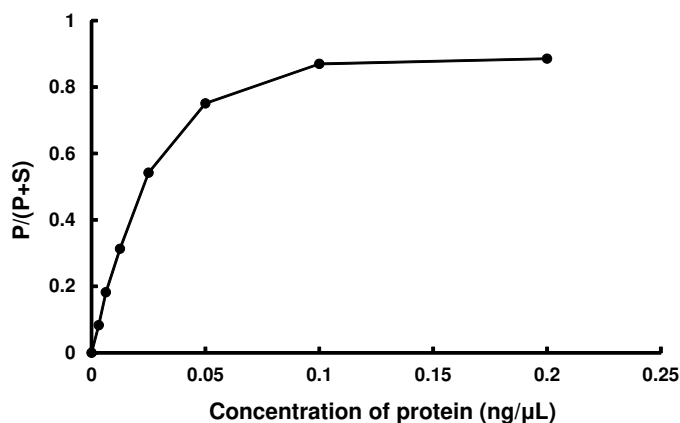
SDS-PAGE



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