

Product Information

BTN-BTK[M437R][non-activated]

Product Number : 08-705-23N

Product description

Full-length human BTK [2-659(end) amino acids and M437R 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.

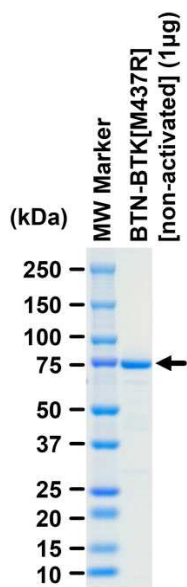
Storage buffer:

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.

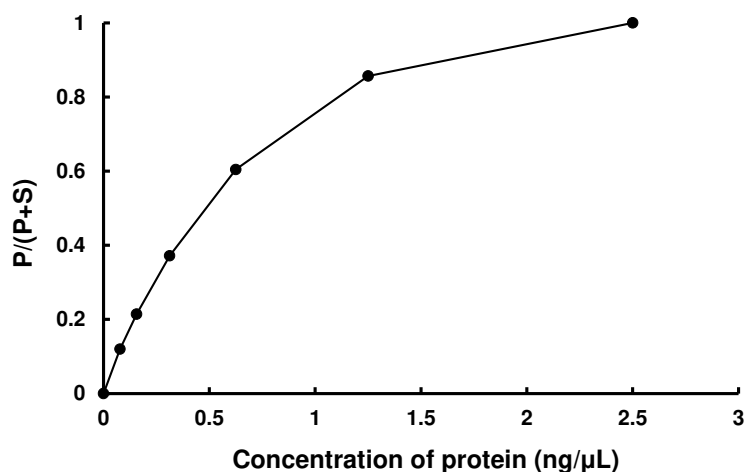
SDS-PAGE



Purity: 92 %

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