

## Product Information

### BTK[C481S]

 Product Number : **08-547**

#### Product description

Full-length human BTK [2-659(end) amino acids and C481S of accession number NP\_000052.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-BTK[C481S] was purified by using glutathione sepharose chromatography.

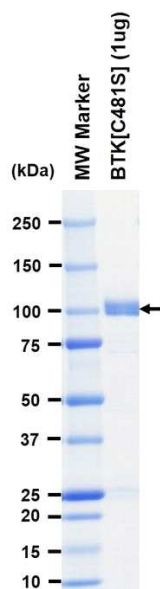
#### Storage buffer:

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

#### Storage and Handling:

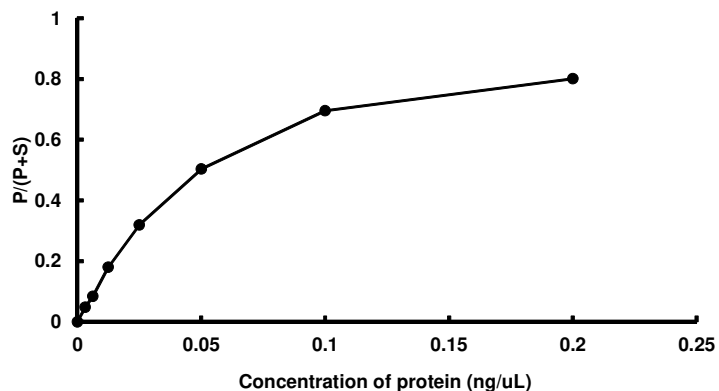
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 85 %  
 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 : Srcptide

ATP : 100  $\mu$ M