

## Product Information

### BTN-ABL(ABL1)[T315I]

Product Number : **08-493-20N**

#### Product description

Full-length human ABL [2-1130(end) amino acids and T315I of accession number NP\_005148.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (127 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

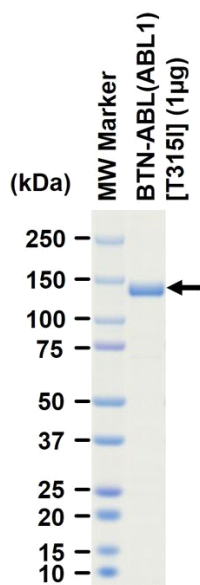
#### Storage buffer:

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.

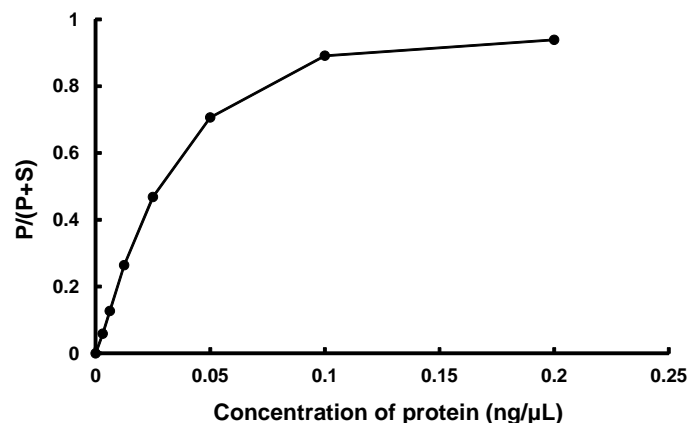
#### SDS-PAGE



Purity: 96 %

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 : ABLtide

ATP : 100 µM