

## Product Information

### BTN-IKK $\alpha$ (CHUK)[inactive mutant]

 Product Number : **05-412-11-20N**

#### Product description

Full-length human IKK $\alpha$  [1-745(end) amino acids and K44A of accession number NP\_001269.3] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (88 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

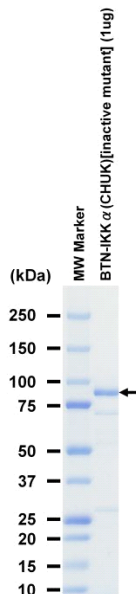
#### Storage buffer:

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

#### Storage and Handling:

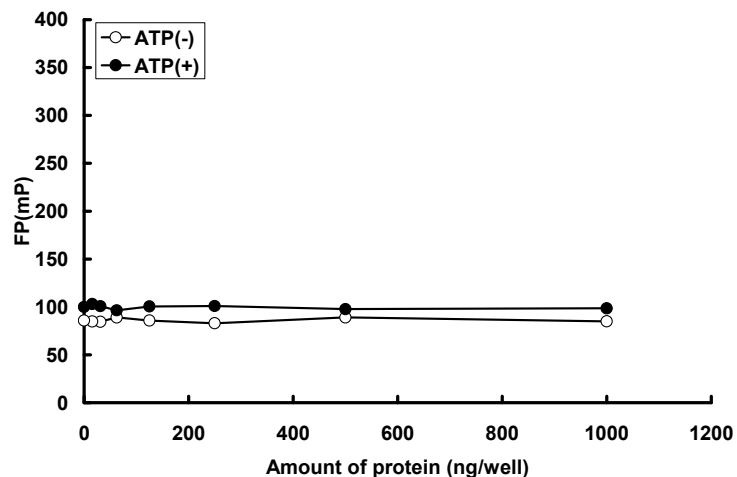
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 80 %  
 The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : IkBa peptide

ATP : 100  $\mu$ M