

## Product Information

### IKK $\alpha$ (CHUK)

 Product Number : **05-112**

#### Product description

Full-length human IKK $\alpha$  [1-745(end) amino acids of accession number NP\_001269.3] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-IKK $\alpha$  was purified by using glutathione sepharose chromatography.

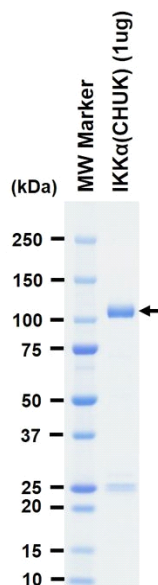
#### Storage buffer:

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.

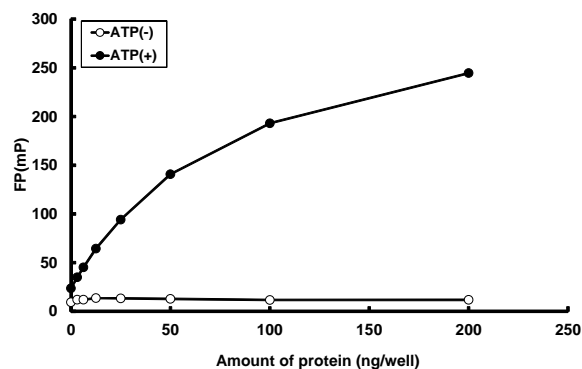
#### SDS-PAGE



Purity: 77 %

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 : I $\kappa$ B $\alpha$  peptide

 ATP : 100  $\mu$ M