

Product Information

BTN-IKK α (CHUK)

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

Product description

Full-length human IKK α [1-745(end) amino acids 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:

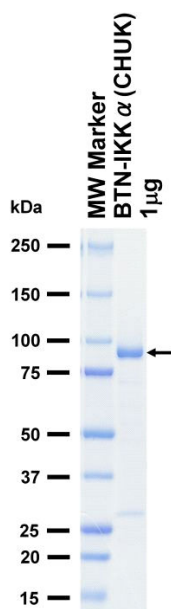
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

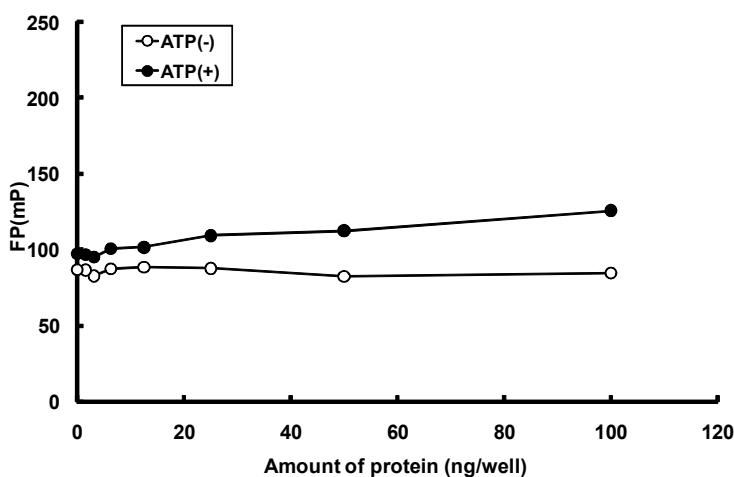
SDS-PAGE



Purity: 88 %

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 μ M