

Product Information

NLK

 Product Number : **04-151**

Product description

Truncated human NLK [121-515(end) amino acids of accession number NP_057315.2] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-NLK was purified by using glutathione sepharose chromatography .

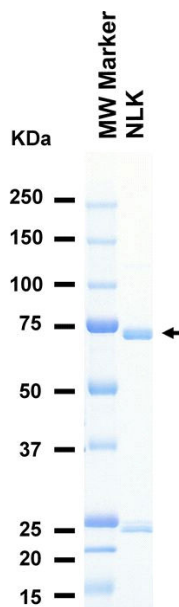
Storage buffer:

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

Storage and Handling:

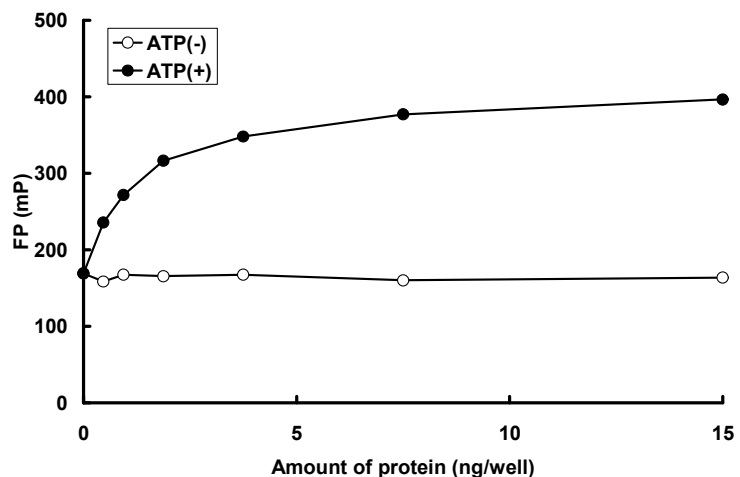
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 62 %
 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 : Multide

ATP : 100 μM