

Product Information

NLK

Product Number : **04-151**

Product description

Truncated human NLK [133-527(end) amino acids of accession number NP_057315.3] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-NLK was purified by using glutathione sepharose chromatography.

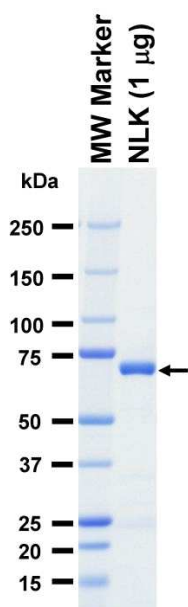
Storage buffer:

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.

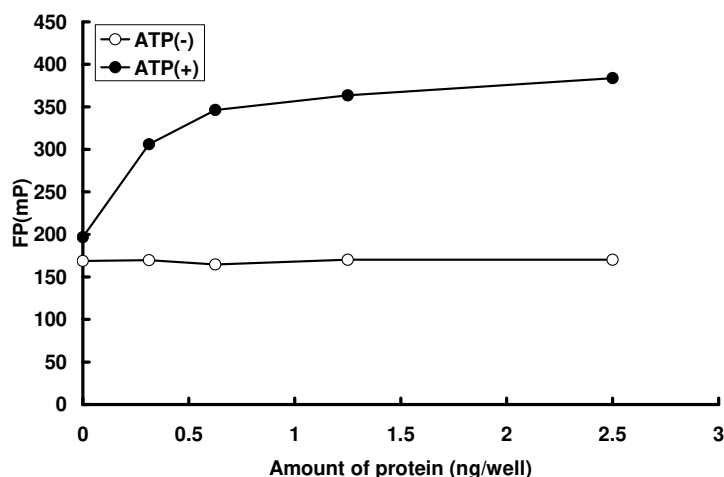
SDS-PAGE



Purity: 92 %

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