

Product Information

WNK2

Product Number : **05-180**

Product description

Human WNK2, catalytic domain [166-489 amino acids of accession number NP_006639.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-WNK2 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

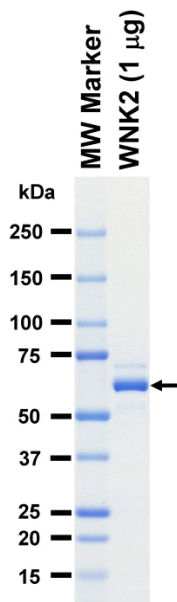
Storage buffer:

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.

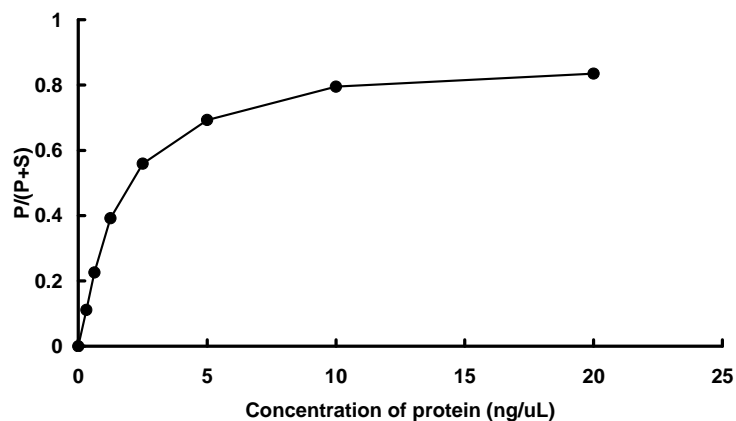
SDS-PAGE



Purity: 93 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mn/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : SPAKtide

ATP : 100 μ M