

Product Information

WNK4

Product Number : **05-182**

Product description

Human WNK4, catalytic domain [1-444 amino acids of accession number NP_115763.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-WNK4 was purified by using glutathione sepharose chromatography.

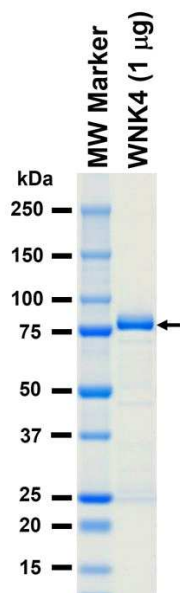
Storage buffer:

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

Storage and Handling:

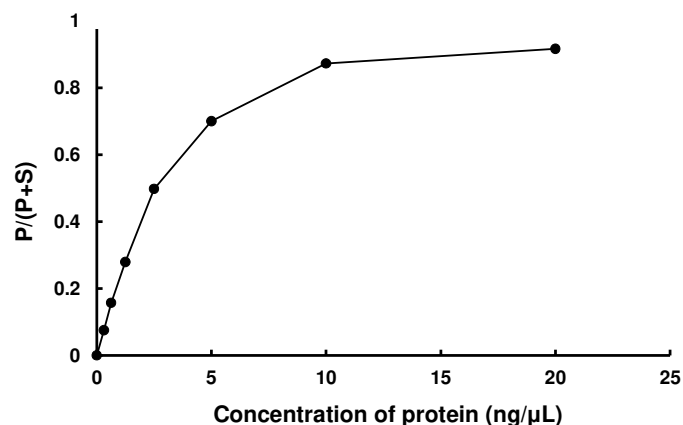
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 86 %
 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 : 200 µM