

## Product Information

### DYRK4

Product Number : **04-434**

#### Product description

Full-length human DYRK4 [1-520(end) amino acids of accession number NP\_003836.1] was expressed as N-terminal DYKDDDDK tagged protein (61 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

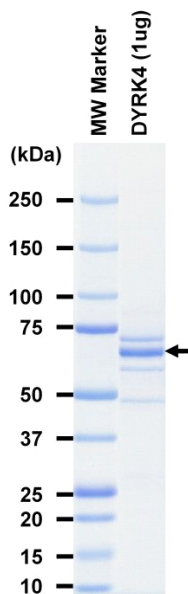
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

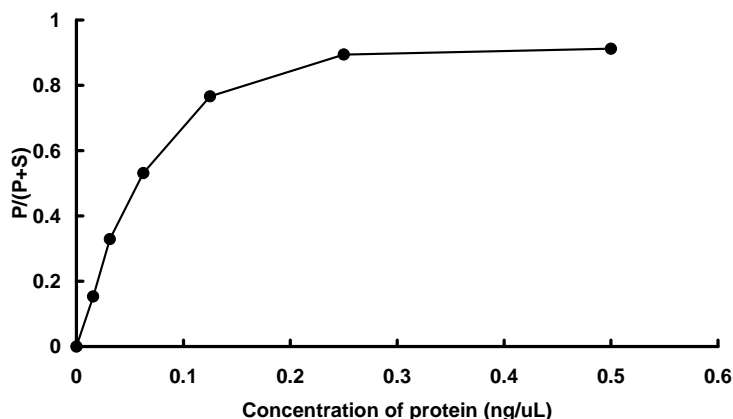
#### SDS-PAGE



Purity: 60 %

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

#### Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : DYRKtide

ATP : 100  $\mu$ M