

## Product Information

### GSK3 $\beta$ (GSK3B)

 Product Number : **04-141**

#### Product description

Full-length human GSK3 $\beta$  [1-420(end) amino acids of accession number NP\_001139628.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-GSK3 $\beta$  was purified by using glutathione sepharose chromatography.

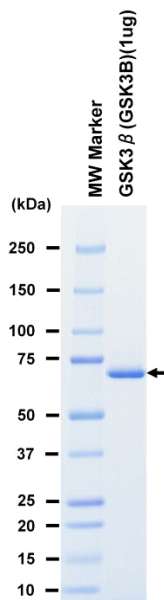
#### Storage buffer:

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.

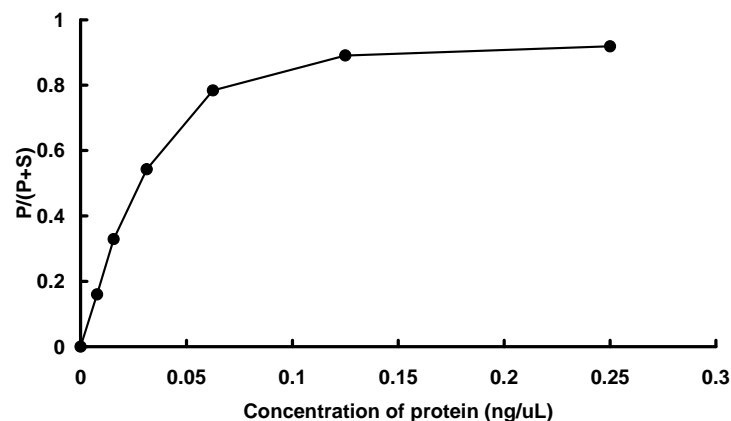
#### SDS-PAGE



Purity: 99 %

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 Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : CREBtide-p

 ATP : 100  $\mu$ M