

## Product Information

### GSK3 $\beta$ (GSK3B)

Product Number : **04-041**

#### Product description

Full-length human GSK3 $\beta$  [1-420(end) amino acids of accession number P49841] was expressed as N-terminal His-tagged protein (50 kDa) using baculovirus expression system. His-tagged GSK3 $\beta$  was purified by using Ni-NTA affinity chromatography.

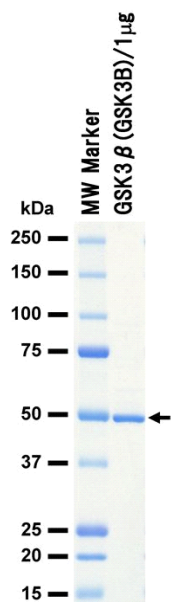
#### Storage buffer:

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.

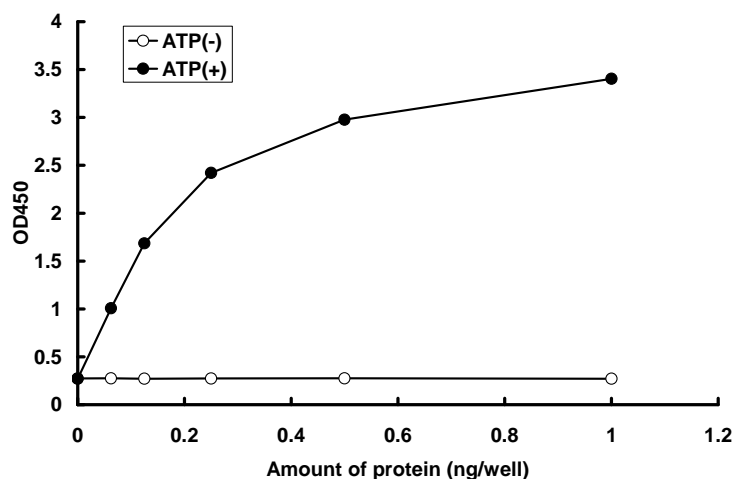
#### SDS-PAGE



Purity: 98 %

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

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated peptide in streptavidin-coated ELISA plate. Phosphorylation was detected by HRP-labeled anti-phospho antibody.

Substrate : GSK3 substrate peptide2

ATP : 5  $\mu$ M