

## Product Information

### WEE1

Product Number : **05-177**

#### Product description

Human WEE1, catalytic domain [215-646(end) amino acids of accession number NP\_003381.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-WEE1 was purified by using glutathione sepharose chromatography.

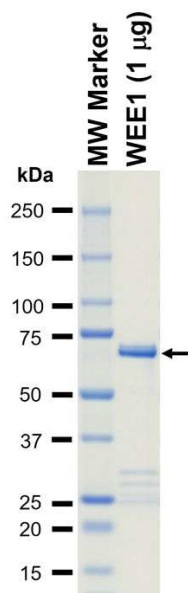
#### Storage buffer:

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.

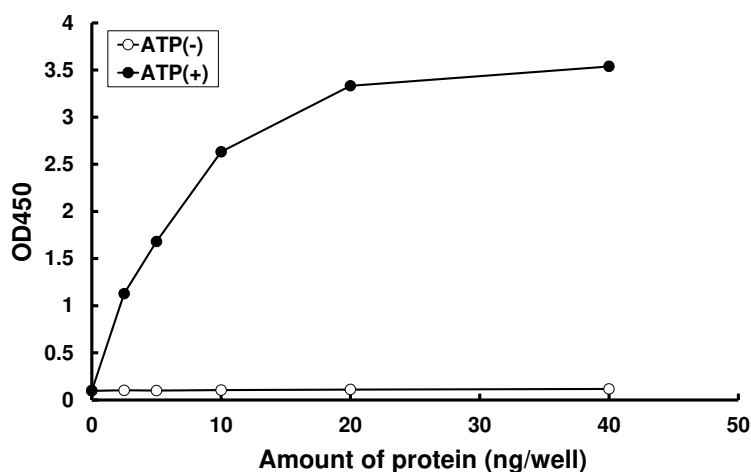
#### SDS-PAGE



Purity: 79 %

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 : CDC2 peptide

ATP : 100 µM