

Product Information

BTN-WEE1

Product Number : **05-477-20N**

Product description

Human WEE1, catalytic domain [215-646(end) amino acids of accession number NP_003381.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (53 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

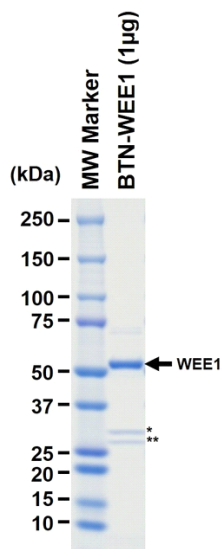
Storage buffer:

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

Storage and Handling:

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

SDS-PAGE

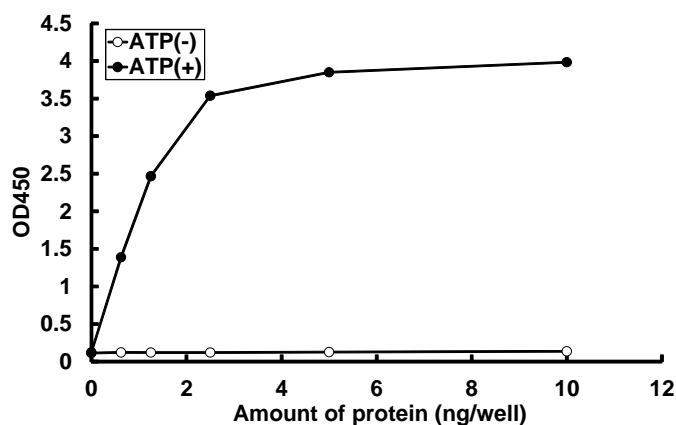


* 14-3-3 epsilon protein from Sf21 cells
** 14-3-3 zeta protein from Sf21 cells

Purity: 74 %

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 μ M