

WEE2 Protein

Full-length recombinant mouse protein expressed in sf9 cells

Catalog # **W06-30BG**

Lot # U1739-8

Product Description

Full-length recombinant mouse WEE2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC052883](#).

Gene Aliases

WEE1B

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

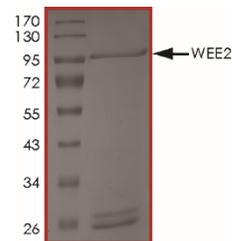
Scientific Background

WEE2, also known as WEE1B, is a dual-specificity kinases that uniquely function to phosphorylate specific Tyr/Thr residues on CDKs. It belongs to WEE1 family that also includes WEE1 homolog 1 (WEE1, also designated as somatic WEE1A) and PKMYT1 (previously known as MYT1). It is conserved from yeast to humans and has been shown to suppress CDK1 activity by phosphorylating its inhibitory sites. A key M phase-promoting factor (MPF) inhibitory kinase in oocytes and is required for maintaining meiotic arrest.

References

1. Leise W 3rd, et al: Multiple Cdk1 inhibitory kinases regulate the cell cycle during development. Dev Biol. 2002 Sep 1;249(1):156-73.
2. Liu C, et al: Ser 15 of WEE1B is a potential PKA phosphorylation target in G2/M transition in one-cell stage mouse embryos. Mol Med Rep. 7(6):1929-37, 2013.

Purity



The purity of WEE2 protein was determined to be **>70%** by densitometry, WEE2 approx. MW **96kDa**.

WEE2 Protein

Full-length recombinant mouse protein expressed in sf9 cells

Catalog # W06-30BG

Lot # U1739-8

Purity >70%

Concentration 0.05 µg/µl

Stability 1yr at -70°C from date of shipment

Storage & Shipping Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.