

Product Information

MYT1(PKMYT1)

 Product Number : **05-176**

Product description

Full-length human MYT1 [1-499(end) amino acids of accession number NP_004194.3] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system. GST-MYT1 was purified by using glutathione sepharose chromatography.

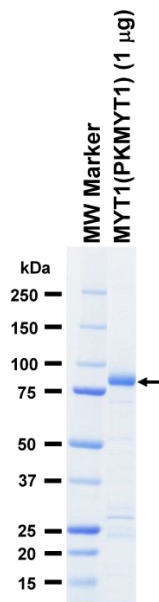
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.2% Brij35, 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

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

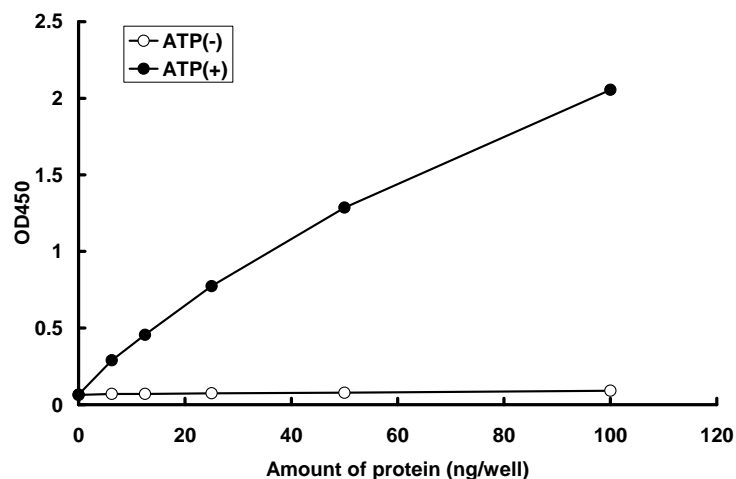
SDS-PAGE



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 μ M