

Product Information

MTOR(FRAP)/MLST8

Product Number : 11-431

Product description

Catalytic domain of human MTOR [1362-2549(end) amino acids of accession number NP_004949.1] was co-expressed as N-terminal DYKDDDDK tagged protein (138 kDa) with MLST8[1-326(end) amino acids of accession number NP_071767.3] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

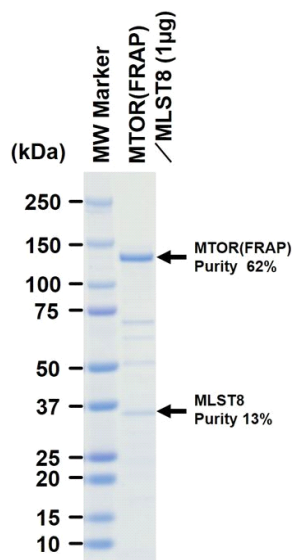
Storage buffer:

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.

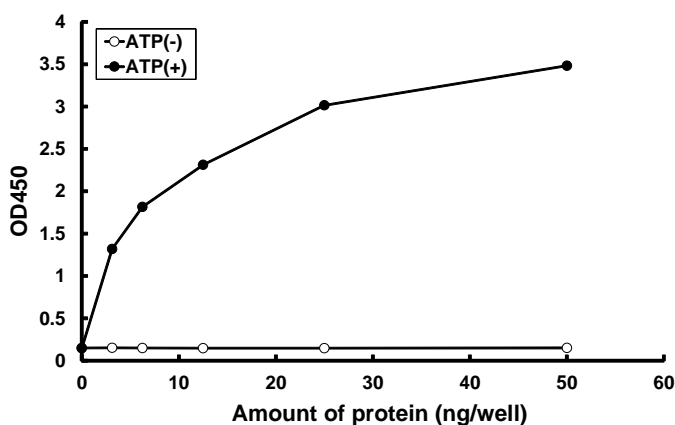
SDS-PAGE



Purity: 75 %

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

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated substrate protein/peptide, and after stopping kinase reaction by EDTA, the reaction solution was transferred into streptavidin-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG).

Substrate : EIF4EBP1

ATP : 100 μ M