

Product Information

COT(MAP3K8)

 Product Number : **07-301**

Product description

Human COT, catalytic domain [30-397 amino acids of accession number NP_005195.2] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-COT was purified by using glutathione sepharose chromatography. GST-COT was cleaved by PreScission protease and GST-free COT (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

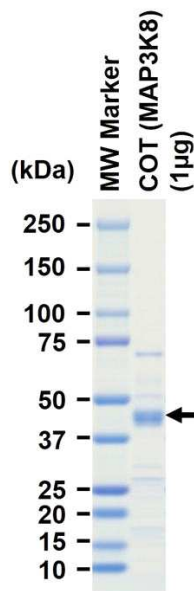
Storage buffer:

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.

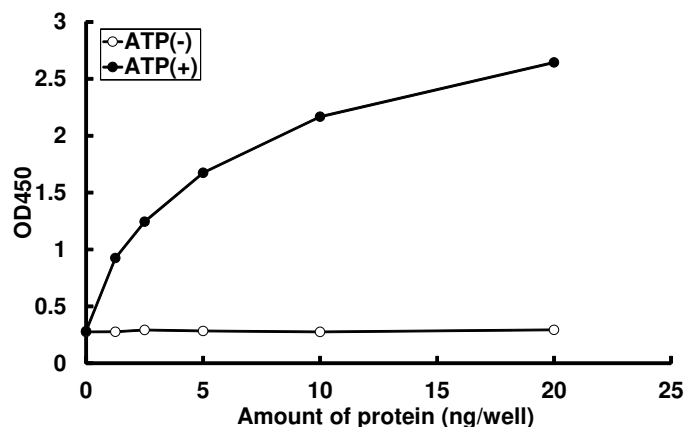
SDS-PAGE



Purity: 68 %

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 : MAP2K1 peptide

 ATP : 100 μ M