

Product Information

BTN-MAPKAPK2

Product Number : **02-442-20N**

Product description

Full-length human MAPKAPK2 [1-400(end) amino acids of accession number NP_116584.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (49 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Notice : This product is especially sensitive to freeze/thaw cycles. After thawing the stock solution, we recommend freezing any left over kinase in single use aliquots to avoid activity loss caused by repeated freeze/thaw cycles.

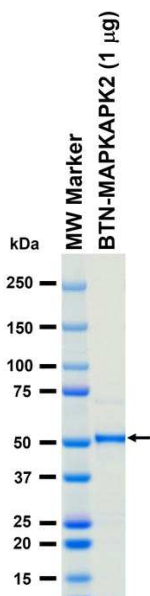
Storage buffer:

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.

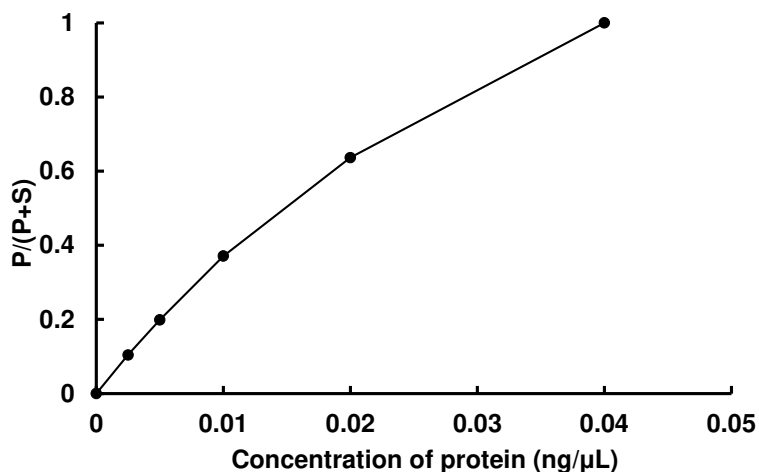
SDS-PAGE



Purity: 96 %

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

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : GS peptide

ATP : 100 µM