

Product Information

MET[D1228H]

Product Number : **08-540**

Product description

Human MET, cytoplasmic domain [956-1390(end) amino acids and D1228H of accession number NP_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-MET[D1228H] was purified by using glutathione sepharose chromatography.

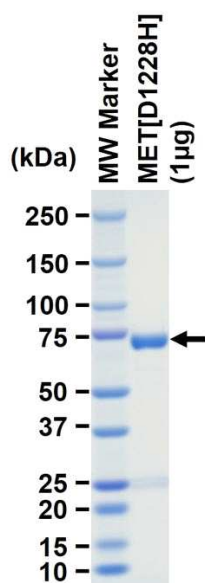
Storage buffer:

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.

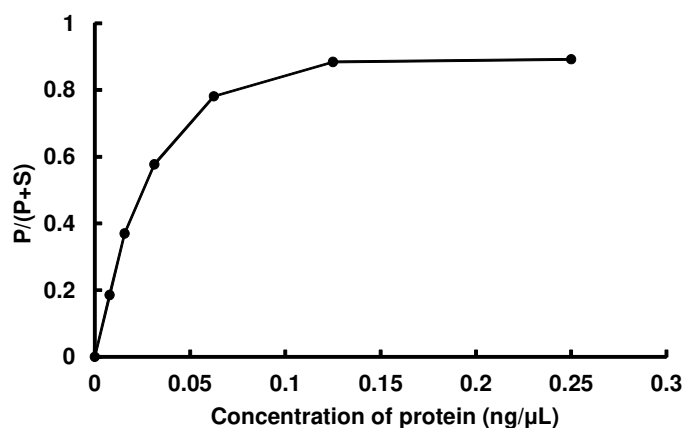
SDS-PAGE



Purity: 87 %

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 : Src tide

ATP : 100 µM