

---

# MET[M1250T]

Product code: 08-545

Human MET, cytoplasmic domain [956-1390(end) amino acids and M1250T of accession number NP\_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : Srctide  
  
ATP ( $\mu$ M) Km app / Bin : 17 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 86  
IC50 at 1 mM ATP (nM) : 1900

---