

---

# MET[D1228H]

Product code: 08-540

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.

Assay platform	: MSA
Substrate	: Srctide
ATP ( $\mu$ M) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 59
IC50 at 1 mM ATP (nM)	: 1200

