

---

# **MET[Y1235D]**

Product code: 08-198

Human MET, cytoplasmic domain [956-1390(end) amino acids and Y1235D 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	: 71 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 79
IC50 at 1 mM ATP (nM)	: 390

