
MET

Product code: 08-151

Human MET, cytoplasmic domain [956-1390(end) amino acids 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 (μ M) Km app / Bin : 27 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 67
IC50 at 1 mM ATP (nM) : 730
