
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. GST-MET was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_{mapp} / B_{in} : 27 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 67

IC₅₀ at 1 mM ATP (nM) : 730
