
MST1(STK4)

Product code 07-116

Full-length human MST1 [1-487(end) amino acids of accession number NP_006273.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-MST1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : IRS1

ATP (μ M) K_{mapp} / B_{in} : 50 / 50

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 1.0

IC₅₀ at 1 mM ATP (nM) : 0.55
