

---

# 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.

Assay platform : MSA

Substrate : IRS1

ATP ( $\mu$ M) Km app / Bin : 50 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 1.0

IC50 at 1 mM ATP (nM) : 0.55

---