

---

## MST2(STK3)

Product code: 07-117

Full-length human MST2 [1-491(end) amino acids of accession number NP\_006272.2] 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 : 69 / 75

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 6.7

IC50 at 1 mM ATP (nM) : 3.1

---