

---

---

## 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

