

---

## 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

