
MST3(STK24)

Product code: 07-118

Full-length human MST3 [1-431(end) amino acids of accession number NP_001027467.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Moesin-derived peptide
ATP (μ M) Km app / Bin	: 66 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.9
IC50 at 1 mM ATP (nM)	: n.a.
