
TIE2(TEK)

Product code: 08-185

Human TIE2, cytoplasmic domain [771-1124(end) amino acids of accession number NP_000450.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 94 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 100
IC50 at 1 mM ATP (nM)	: 190

