

---

## ***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. GST-TIE2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : Blk/Lyntide

ATP ( $\mu$ M) Kmapp / Bin : 94 / 100  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 100  
IC50 at 1 mM ATP (nM) : 190

---