

---

# FLT1

Product code: 08-189

Human FLT1, cytoplasmic domain [781-1338(end) amino acids of accession number NP\_002010.1] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CSKtide  
  
ATP ( $\mu$ M) Km app / Bin : 140 / 150  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 3.2  
IC50 at 1 mM ATP (nM) : 6.8

---