

---

# FLT4

Product code: 08-190

Human FLT4, cytoplasmic domain [798-1298(end) amino acids and H890Q of accession number NP\_002011.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CSKtide  
  
ATP ( $\mu$ M) Km app / Bin : 72 / 75  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.66  
IC50 at 1 mM ATP (nM) : 2.4

---