

---

## ***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. GST-FLT1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 140 / 150

Metal : Mg

Reference compound : Staurosporine

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 3.2

IC<sub>50</sub> at 1 mM ATP (nM) : 6.8

---