

---

# FGFR1

Product code 08-133

Human FGFR1, cytoplasmic domain [398-822(end) amino acids of accession number NP\_075598.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP ( $\mu$ M) Km app / Bin : 89 / 100

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.3

IC50 at 1 mM ATP (nM) : 12

---