

---

---

# FGFR2

Product code: 08-134

Human FGFR2, cytoplasmic domain [399-821(end) amino acids of accession number NP\_000132.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 66 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 5.4

