

---

---

# 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.

Assay platform	: MSA
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

