

---

## FGFR3[K650E]

Product code: 08-501

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and K650E of accession number NP\_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP ( $\mu$ M) Km app / Bin	: 41 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 14

---