

---

---

## FGFR2[V564I]

Product code: 08-546

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and V564I 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	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 47

---

---