

---

## **FGFR1[V561M]**

Product code: 08-536

Human FGFR1, cytoplasmic domain [398-822(end) amino acids and V561M 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 : 33 / 50  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.14  
IC50 at 1 mM ATP (nM) : 1.3

---