

---

# FGFR4

Product code: 08-136

Human FGFR4, cytoplasmic domain [460-802(end) amino acids of accession number NP\_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CSKtide  
  
ATP ( $\mu$ M) Km app / Bin : 230 / 250  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 43  
IC50 at 1 mM ATP (nM) : 120

---