

---

# 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

---