

---

---

## FGFR4[V550E]

Product code: 08-525

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and V550E 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	: 210 / 200
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 370
IC50 at 1 mM ATP (nM)	: 1300

