

---

## FGFR4[N535K]

Product code: 08-524

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and N535K 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	: 30 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 160
IC50 at 1 mM ATP (nM)	: 1200

