
FGFR2

Product code 08-134

Human FGFR2, cytoplasmic domain [399-821(end) amino acids of accession number NP_000132.1] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP (μ M) K_{mapp} / B_{in} : 66 / 75

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 1.4

IC₅₀ at 1 mM ATP (nM) : 5.4
