
FGFR2[V564I]

Product code 08-546

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and V564I of accession number NP_000132] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR2[V564I] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CSKtide

ATP (μ M) Km app / Bin : 21 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.0

IC50 at 1 mM ATP (nM) : 47
