
FGFR3[K650E]

Product code: 08-501

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and K650E of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CSKtide

ATP (μ M) Km app / Bin : 41 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.2
IC50 at 1 mM ATP (nM) : 14
