
FGFR3[K650M]

Product code: 08-199

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and K650M 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	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.68
IC50 at 1 mM ATP (nM)	: 17
