
FGFR3[V555L]

Product code: 08-548

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and V555L 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	: 29 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.49
IC50 at 1 mM ATP (nM)	: 9.4