
FGFR3[V555M]

Product code: 08-543

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and V555M 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	: 37 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.21
IC50 at 1 mM ATP (nM)	: 1.8

