
FGFR4[V550L]

Product code: 08-526

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and V550L of accession number NP_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 160 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 44
