
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
