
FGFR1

Product code: 08-133

Human FGFR1, cytoplasmic domain [398-822(end) amino acids of accession number NP_075598.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : CSKtide

ATP (μ M) Km app / Bin : 89 / 100

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 2.3

IC50 at 1 mM ATP (nM) : 12
