
FLT4

Product code: 08-190

Human FLT4, cytoplasmic domain [798-1298(end) amino acids of accession number NP_002011.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CSKtide

ATP (μ M) Km app / Bin : 72 / 75
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 0.66
IC50 at 1 mM ATP (nM) : 2.4
