
FLT4

Product code: 08-190

Human FLT4, cytoplasmic domain [798-1298(end) amino acids and H890Q of accession number NP_002011.2] 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

