
FLT1

Product code: 08-189

Human FLT1, cytoplasmic domain [781-1338(end) amino acids of accession number NP_002010.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 140 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: 6.8
