
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
