
FLT3

Product code 08-154

Human FLT3, cytoplasmic domain [564-993(end) amino acids of accession number NP_004110.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-FLT3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Src tide

ATP (μ M) K_m app / B_{in} : 94 / 100

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 0.20

IC₅₀ at 1 mM ATP (nM) : 0.34
