
EPHA1

Product code 08-119

Human EPHA1, cytoplasmic domain [586-976(end) amino acids of accession number NP_005223.3] was expressed as N-terminal GST-fusion protein (71 kDa) using baculovirus expression system. GST-EPHA1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 22 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 20

IC50 at 1 mM ATP (nM) : 340
