
EPHA2

Product code 08-121

Human EPHA2, cytoplasmic domain [572-976(end) amino acids of accession number NP_004422.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHA2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 67 / 75

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 160

IC50 at 1 mM ATP (nM) : 530
