
EPHA6

Product code 08-125

Human EPHA6, cytoplasmic domain [683-1130(end) amino acids of accession number NP_001073917.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-EPHA6 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 27 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 17

IC50 at 1 mM ATP (nM) : 60
