
EPHA7

Product code 08-126

Human EPHA7, cytoplasmic domain [595-998(end) amino acids of accession number NP_004431.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHA7 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 58 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 48

IC50 at 1 mM ATP (nM) : 480
