
EPHB4

Product code 08-131

Human EPHB4, cytoplasmic domain [577-987(end) amino acids of accession number NP_004435.3] was expressed as N-terminal GST-protein (73 kDa) using baculovirus expression system. GST-EPHB4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Blk/Lyntide

ATP (μ M) Kmapp / Bin : 56 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 230

IC50 at 1 mM ATP (nM) : 1500
