
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.

Assay platform : MSA
Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 56 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 230
IC50 at 1 mM ATP (nM) : 1500
