
EPHA1

Product code: 08-119

Human EPHA1, cytoplasmic domain [586-976(end) amino acids and M900V of accession number NP_005223.4] was expressed as N-terminal GST-fusion protein (71 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 22 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 20
IC50 at 1 mM ATP (nM) : 340
