
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

