
EPHA2

Product code: 08-121

Human EPHA2, cytoplasmic domain [572-976(end) amino acids of accession number NP_004422.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 67 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 160
IC50 at 1 mM ATP (nM)	: 530