
EPHA6

Product code: 08-125

Human EPHA6, cytoplasmic domain [683-1130(end) amino acids of accession number NP_001073917.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 17
IC50 at 1 mM ATP (nM)	: 60
