
EPHA4

Product code: 08-123

Human EPHA4, cytoplasmic domain [586-986(end) amino acids of accession number NP_004429.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 52 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 50
IC50 at 1 mM ATP (nM)	: 330