
EPHA3

Product code: 08-122

Human EPHA3, cytoplasmic domain [579-983(end) amino acids of accession number NP_005224.2] 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	: 170 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 37
IC50 at 1 mM ATP (nM)	: 76

