
EPHA5

Product code: 08-124

Human EPHA5, catalytic domain [662-948 amino acids of accession number NP_004430.4] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 34
IC50 at 1 mM ATP (nM)	: 220

