
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
