
EPHA8

Product code: 08-127

Human EPHA8, catalytic domain [571-924 amino acids of accession number NP_065387.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 69 / 75
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 61
IC50 at 1 mM ATP (nM) : 240
