
PIK3CA[P539R]/PIK3R1

Product code: 11-412-20N

Full-length human PIK3CA[1-1068(end) amino acids and P539R of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo_Lipid

Substrate : PI(4,5)P2

ATP (μ M) Km app / Bin : 37 / 50

Metal : Mg

Reference compound : PI-103

IC50 at ATP Bin (nM) : 8.7

IC50 at 1 mM ATP (nM) : n.a.