
PIK3CA[E542K]/PIK3R1

Product code: 11-413-20N

Full-length human PIK3CA[1-1068(end) amino acids and E542K 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
Substrate	: PI(4,5)P2
ATP (μ M) Km app / Bin	: 42 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.
