

---

---

# 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\_Lipid

Substrate : PI(4,5)P2

ATP ( $\mu$ 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.