
PIK3CD/PIK3R1

Product code: 11-403-20N

Full-length human PIK3CD[1-1044(end) amino acids of accession number NP_005017.3] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (123 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) : 24

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