
PIK3CB/PIK3R1

Product code 11-402-20N

Full-length human PIK3CB[1-1070(end) amino acids of accession number NP_006210.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (126 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] (84kDa) 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) K_{mapp} / B_{in} : 88 / 100

Metal : Mg

Reference compound : PI-103

IC50 at ATP B_{in} (nM) : 22

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