

---

# 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] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo  
Substrate : PI(4,5)P2  
  
ATP ( $\mu$ M) Km app / Bin : 88 / 100  
Metal : Mg  
Reference compound : PI-103  
IC50 at ATP Bin (nM) : 22  
IC50 at 1 mM ATP (nM) : n.a.

---