

---

## ***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] (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 ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 37 / 50

Metal : Mg

Reference compound : PI-103

IC50 at ATP  $B_{in}$  (nM) : 24

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