

Original Papers:

- (1) Ho LK et al. (1997) Identification of four novel human phosphoinositide 3-kinases defines a multi-isoform subfamily. *Biochem Biophys Res Commun.* **235**:130
- (2) Vanhaesebroeck B et al. (1997) P110delta, a novel phosphoinositide 3-kinase in leukocytes. *Proc Natl Acad Sci U S A*, **94**:433
- (3) Stoyanov B et al. (1995) Cloning and characterization of a G protein-activated human phosphoinositide-3 kinase. *Science.* **269**:690
- (4) Volinia S et al. (1994) Molecular cloning, cDNA sequence, and chromosomal localization of the human phosphatidylinositol 3-kinase p110 alpha (PIK3CA) gene. *Genomics.* **24**:472
- (5) Hu P et al. (1993) Cloning of a novel, ubiquitously expressed human phosphatidylinositol 3-kinase and identification of its binding site on p85. *Mol Cell Biol.* **13**:7677
- (6) Escobedo JA et al. (1991) cDNA cloning of a novel 85 kd protein that has SH2 domains and regulates binding of PI3-kinase to the PDGF beta-receptor. *Cell.* **65**:75