

Reviews:

- (1) Wetzker R et al. (2004) Phosphoinositide 3-kinases as targets for therapeutic intervention. *Curr Pharm Des.* **10**:1915
- (2) Fresno Vara JA et al. (2004) PI3K/Akt signalling pathway and cancer. *Cancer Treat Rev.* **30**:193
- (3) Cooray S (2004) The pivotal role of phosphatidylinositol 3-kinase-Akt signal transduction in virus survival. *J Gen Virol.* **85**:1065
- (4) Hannigan MO et al. (2004) Roles of PI3K in neutrophil function. *Curr Top Microbiol Immunol.* **282**:165
- (5) Wymann MP et al. (2003) Phosphoinositide 3-kinase gamma: a key modulator in inflammation and allergy. *Biochem Soc Trans.* **31**:275
- (6) Anderson KE et al. (2003) Class I phosphoinositide 3-kinases. *Int J Biochem Cell Biol.* **35**:1028
- (7) Koyasu S (2003) The role of PI3K in immune cells. *Nat Immunol.* **4**:313
- (8) Franke TF (2003) PI3K/Akt and apoptosis: size matters. *Oncogene* **22**:8983
- (9) Rodgers EE et al. (2002) Functions of PI3-kinase in development of the nervous system. *Int J Dev Neurosci.* **20**:187
- (10) Vivanco I et al. (2002) The phosphatidylinositol 3-Kinase AKT pathway in human cancer. *Nat Rev cancer.* **2**:489
- (11) Cantley LC (2002) The phosphoinositide 3-kinase pathway. *Science.* **296**:1655
- (12) Vanhaesebroeck B et al. (1997) Phosphoinositide 3-kinases: a conserved family of signal transducers. *Trends Biochem Sci.* **22**:267
- (13) Parker PJ et al. (1992) Phosphatidylinositol 3-kinase: a novel effector. *Cell Growth Differ.* **3**:747