

## Recommended Reading:

- (1) Larribere L et al. (2004) PI3K mediates protection against TRAIL-induced apoptosis in primary human melanocytes. *Cell Death Differ.* **11**:1084
- (2) Viard P et al. (2004) PI3K promotes voltage-dependent calcium channel trafficking to the plasma membrane. *Nature Neuroscience.* **7**:939
- (3) Yip SC et al. (2004) Over-expression of the p110beta but not p110alpha isoform of PI 3-kinase inhibits motility in breast cancer cells. *Cell Motil Cytoskeleton.* **59**:180
- (4) Siddiqui IA et al. (2004) Modulation of phosphatidylinositol-3-kinase/protein kinase B-and mitogen-activated protein kinase-pathways by tea polyphenols in human prostate cancer cells. *J Cell Biochem.* **91**:232
- (5) Huang W et al. (2004) Impaired activation of phosphatidylinositol 3-kinase by leptin is a novel mechanism of hepatic leptin resistance in diet-induced obesity. *J Biol Chem.* **279**:21695
- (6) Hess KL et al. (2004) Frontline: The p85alpha isoform of phosphoinositide 3-kinase is essential for a subset of B cell receptor-initiated signaling responses. *Eur J Immunol.* (Epub ahead of print)
- (7) Weiss-Haljiti C et al. (2004) Involvement of Phosphoinositide 3-Kinase {gamma}, Rac, and PAK Signaling in Chemokine-induced Macrophage Migration. *J Biol Chem.* **279**:43273
- (8) Ali K et al. (2004) Essential role for the p110delta phosphoinositide 3-kinase in the allergic response. *Nature.* **431**:1007