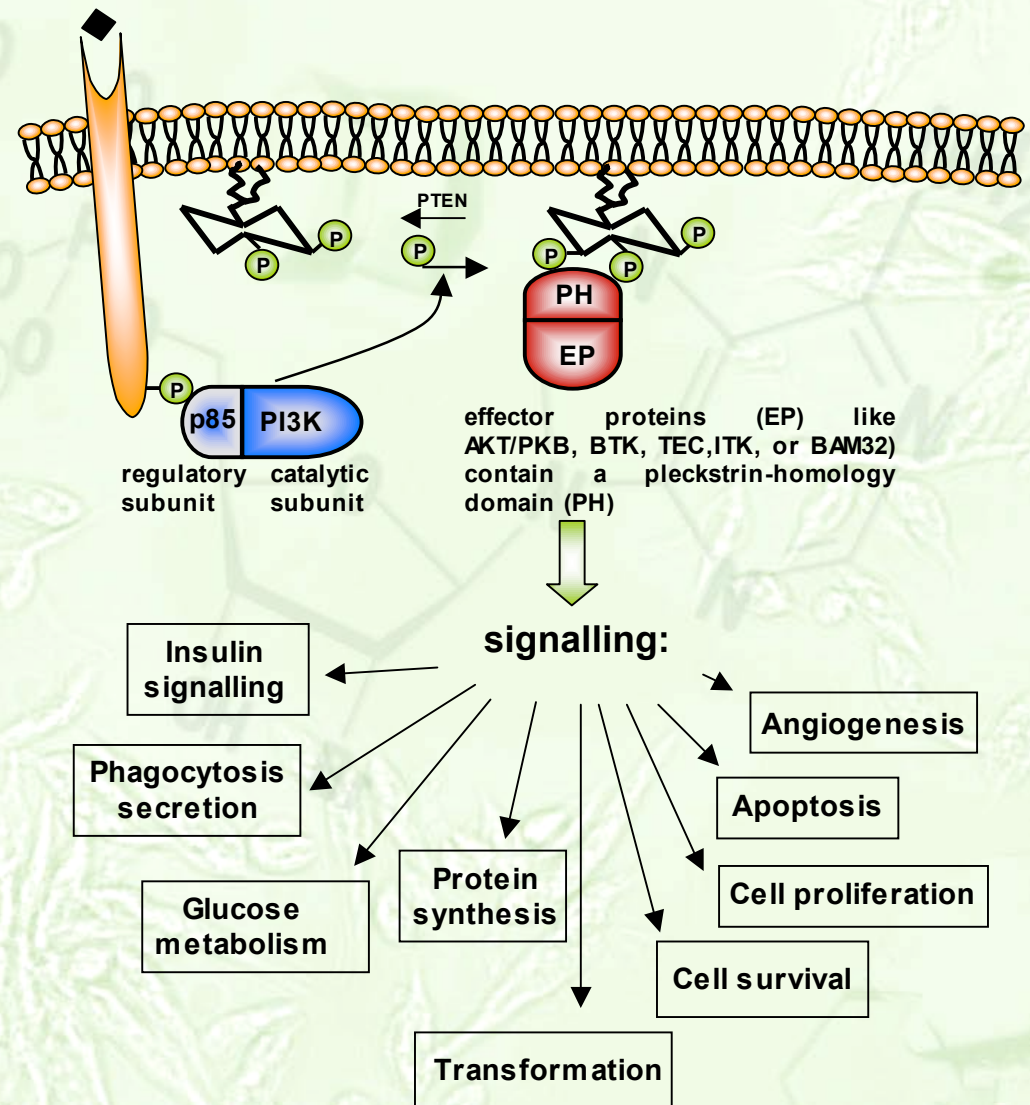


Phosphoinositide 3-kinases – a protein family of high medical relevance

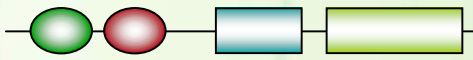



Phosphoinositide 3-kinases (PI3Ks):






- ▶ play a pivotal role in important cellular regulatory mechanisms such as cell growth, proliferation and survival/apoptosis
- ▶ play key roles in cardiac growth (PI3K α), cardiac contractability and the function of β -adrenergic receptors (PI3K γ)
- ▶ are thus of outstanding medical interest
 - diabetes
 - various types of cancer
 - chronic inflammation
 - cardiovascular disease
- ▶ are targets for novel cancer therapeutics and/or immunosuppressants



Phosphoinositide 3-kinases are divided into 3 classes

- ▶ Phosphoinositide 3-kinases are lipid kinases that catalyze the transfer of the γ - phosphate from ATP to the 3'-hydroxyl group of Phosphatidylinositol (PtdIns) and its derivatives (collectively called phosphoinositides)
- ▶ 8 isoforms of the PI3K family have been isolated from mammalian cells and grouped into *three classes* depending on their primary structure and substrate specificity

class	primary structure	subunits		regulated by	substrate	
		catalytic	regulatory		<i>in vitro</i>	<i>in vivo</i>
I		p110 α,β,δ	p85 α,β p55 γ	tyrosine kinases, Ras	PI PI4P	PI4,5P ₂
						
II		PI3K-C2 α,β,γ	not known	clathrin, chemokines, integrins	PI PI4P	not known
III		Vps34p-Analogs	p150	constitutively active	PI	PI

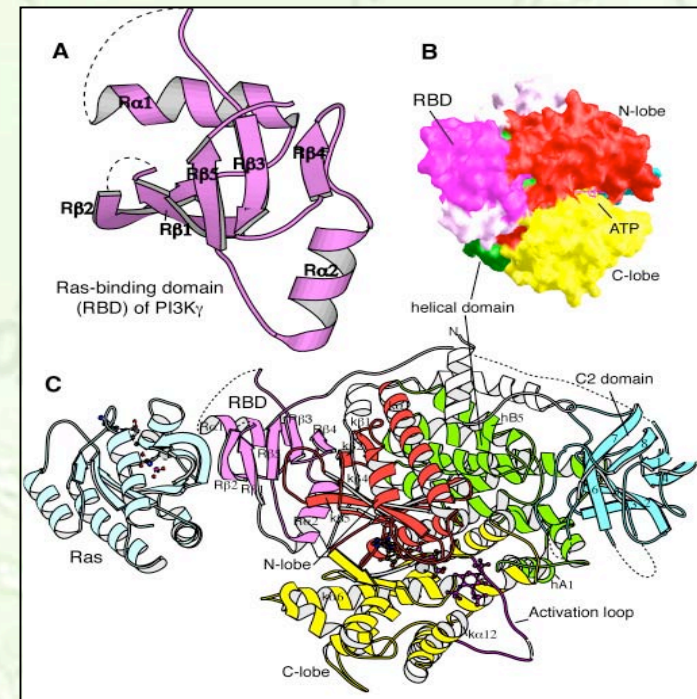
 Adaptor-binding site
  Ras-binding domain
  PI3K accessory domain (PIK)
  catalytic domain
  C2-domain

Phosphoinositide 3-kinases now available from Jena Bioscience

- ▶ produced in insect cells
- ▶ native proteins, affinity-tagged variants and important mutants
- ▶ purity > 95 %
- ▶ quality monitored by SDS-Page and *in vitro* activity assays

Recombinant PI3-kinases are used for:

- ▶ screening strategies to identify PI3K inhibitors and activators
- ▶ target identification for cancer prevention and therapy
- ▶ interaction studies with PI3K signaling modulators
- ▶ phosphorylation assays (lipid and protein kinase assays)



taken from: Walker et al. (1999) Structural insights into phosphoinositide 3-kinase catalysis and signalling. *Nature*. **402**:313.