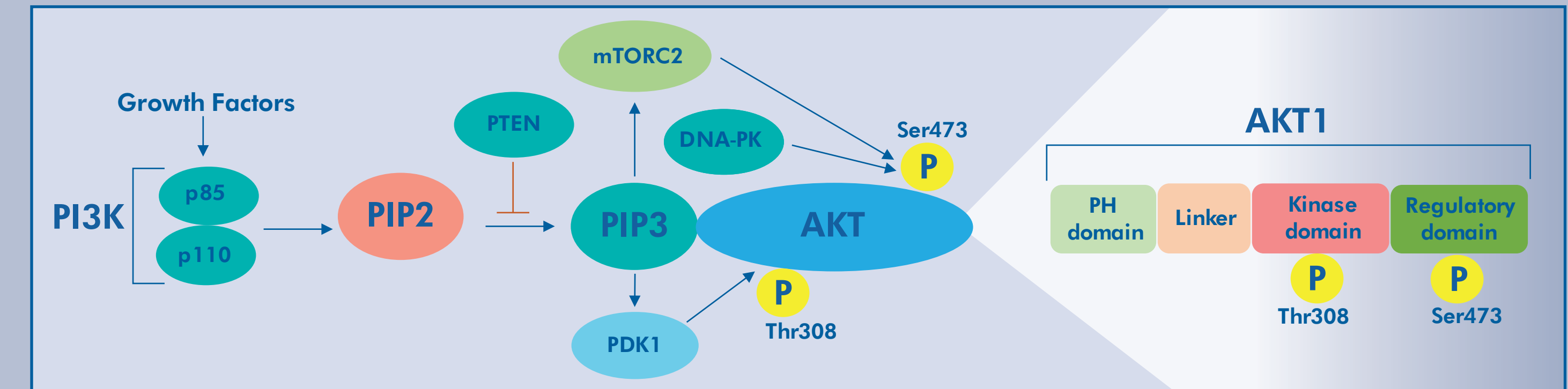
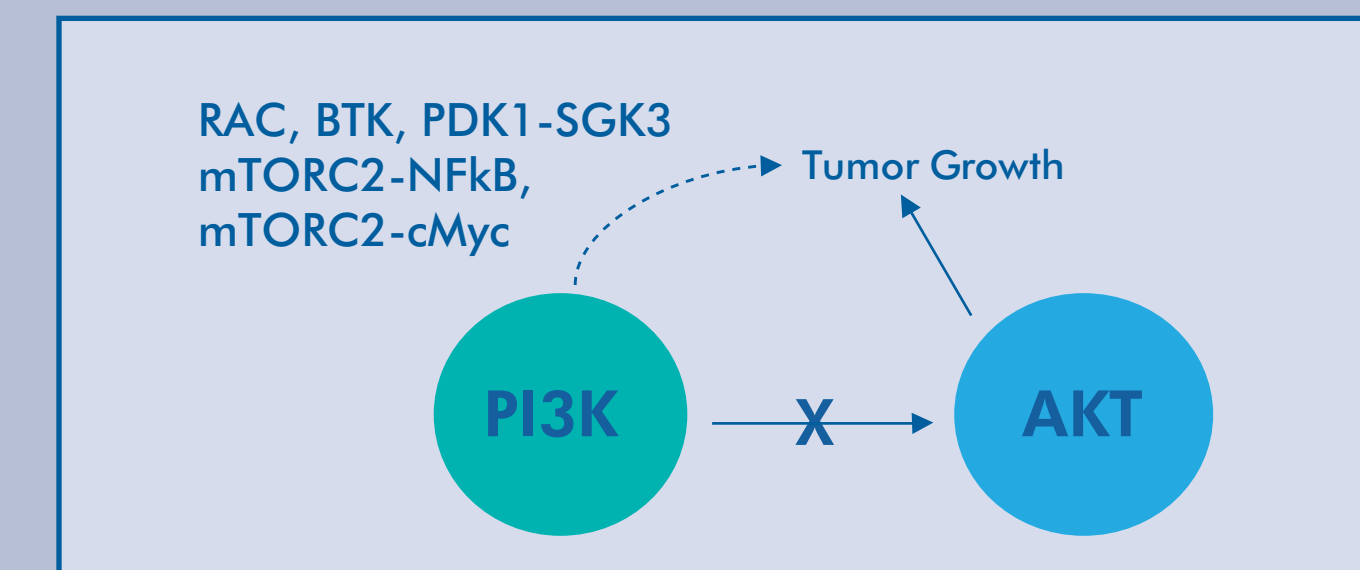


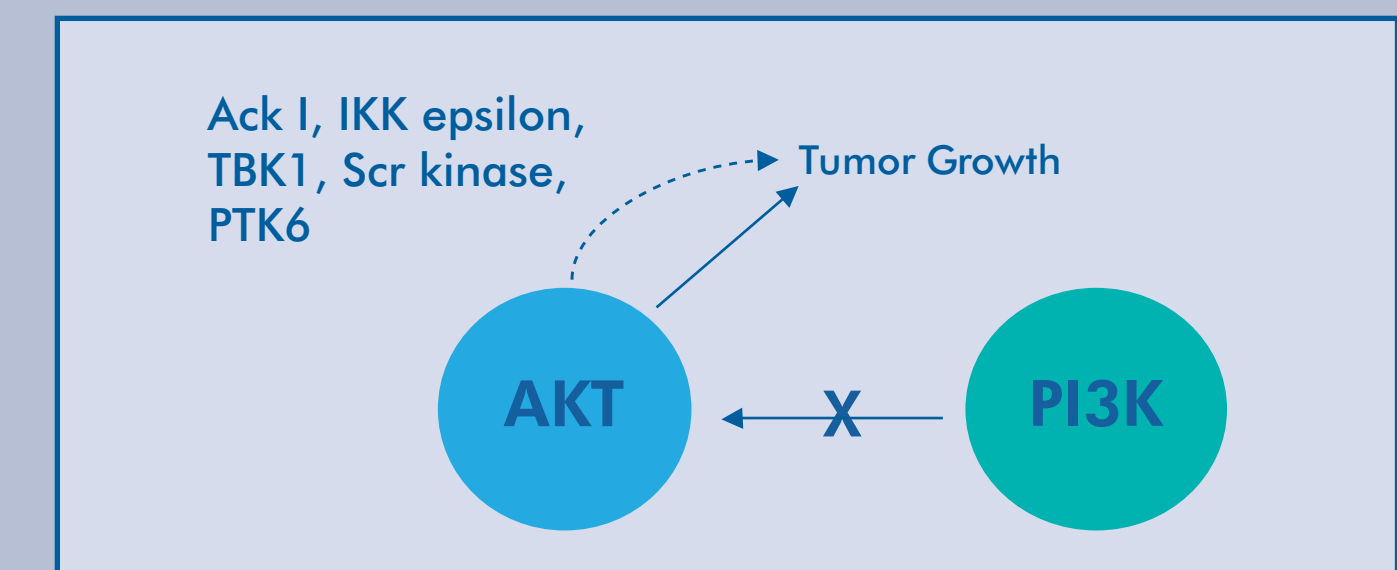
Mechanism of Activation and Major Phosphorylation Sites of AKT1



AKT Independent PI3K Activity



PI3K Independent AKT Activity



PI3K-AKT Signaling: Processes & Targets*

| Cell proliferation | Survival signaling |
|--|--|
| Cyclin D1 p21/Cip p27/Kip Myt1 Wee1 | Bad MDM2/p53 FOXO1 Bcl-2 Bax |
| Metabolism | Translational Control |
| GSK-3 PFKFB2 PIP5K AS160 TBC1D4 PDE3B | S6 kinase RPS6 TSC1 mTOR PDCD4 |
| Neuronal Function | Immunity |
| Huntingtin Ataxin-1 GABAA receptor NfκB | TSC-mTORC1 FOXO1 GSK3β ACLY |
| Invasion/Migration | Others |
| Palladin Girdin Rac/GTP | WNK1 eNOS LMNA |

Small Molecule Inhibitors of AKT*

| | |
|--------------|---|
| Akt (PKB) | Akti-1/2, API-1, API-2, 10-DEBC hydrochloride, FPA 124, GSK 690693, Perifosine, SC 66 |
| Akt and PDK1 | PHT 427 |
| mTOR | Rapamycin, Torin 1, AZD 3147, eCF 309 |
| PDK1 | GSK 2334470, BX 795, OSU 03012 |
| PIP3 | PIT 1 |
| PI3K | LY 294002, Wortmannin, PI 828 |
| PI3K p110α | A66 |
| PI3Kβ | AZD 6482 |
| PI3K p110δ | PI 3065 |
| PI3K γ | AS 605240 |

Other molecules for AKT Signaling*

| Small molecule | Description |
|---------------------------|----------------------------------|
| SC 79 | Activator of Akt |
| 740 YP | Cell permeable activator of PI3K |
| Akt/SKG Substrate Peptide | Synthetic substrate of Akt/PKB |
| AKTide-2T | Peptide substrate for Akt/PKB |

PI3K-AKT Signaling in Diseases

Overgrowth syndromes, cancer (growth, angiogenesis and metastasis), cardiovascular diseases, glucose metabolism related disorders (insulin resistance and diabetes), autoimmune diseases and inflammation, and neurodegeneration related disorders.

*Visit www.novusbio.com, www.rndsystems.com and www.tocris.com to learn more.