Abstract #2897



Discovery and characterization of driver MAPK and PI3K pathway mutations in tumors and association with drug response in cell lines

Mark Tomilo, Paul D. Williams, Emma T. Bowden, Supra R. Gajjala, Santhoshi Bandla, Sean F. Eddy, Seth E. Sadis, Peter J. Wyngaard, Nickolay A. Khazanov, Daniel R. Rhodes Compendia Bioscience, part of Life Technologies, Ann Arbor, MI, United States

Abstract:

The MAPK and PI3K pathways are frequently altered in human cancer and are targeted by dozens of agents in clinical trials. The efficacy of these therapies, alone or in combination, may depend on the activation status of both pathways. Next-generation sequencing of cancer exomes provides a unique opportunity to systematically survey pathway alterations in cancer. Using somatic mutation data obtained from The Cancer Genome Atlas, we sought to catalog the members of the MAPK and PI3K pathways with driver mutations, as well as the frequency of occurrence and co-occurrence of these mutations in common cancers. Furthermore, we sought to characterize the association of pathway mutation status with clinical response and with drug response in pre-clinical models.

While the MAPK and PI3K pathways were frequently altered, the frequency of single and dual pathway alteration and the altered genes varied substantially across cancer types. Predicted Gain of Function and Loss of Function mutations in MAPK pathway genes were most frequently observed in melanoma (77%), thyroid (69%), rectal (62%), colon (56%), lung adenocarcinoma (38%), and uterine cancers (24%). MAPK pathway genes were infrequently mutated in bladder (4%), breast (3%), prostate (1%), and renal cancer (1%). KRAS, BRAF and NRAS hotspot mutations were the most common pathway drivers. Predicted Gain of Function and Loss of Function mutations in PI3K pathway genes were most frequently observed in uterine (78%), breast cancer (36%), cervical squamous cell carcinoma (28%), glioblastoma (24%), head and neck squamous cell carcinoma (20%), squamous cell lung carcinoma (17%), and colon adenocarcinoma (17%). In contrast, PI3K pathway genes were rarely mutated in prostate cancer (1%), thyroid cancer (1%), ovarian serous adenocarcinoma (1%), and acute myeloid leukemia (0%). Hotspot mutations in PIK3CA as well as hotspot and deleterious mutations in *PTEN* were the most common pathway alterations. Predicted driver mutations were observed in *PIK3R1*, and *AKT1*. In uterine cancer, MAPK pathway mutations rarely occurred independently of PI3K pathway mutations while in colon and rectal adenocarcinoma, PI3K pathway mutations rarely occurred independently of MAPK pathway mutations. Co-occurring PI3K and MAPK pathway mutations in colorectal cancer led to worse overall survival compared to MAPK pathway mutations alone. Co-occurring MAPK and PI3K pathway mutations were also observed in melanoma (6%), and gastric (8%) cancers. In other cancer types one pathway was generally favored with the exception of prostate cancer, where pathway activating mutations in either the PI3K or MAPK pathway genes were uncommon.

To support the use of pre-clinical models as exhibiting responses representative of clinical populations, we assessed drug sensitivity to PI3K and MAPK pathway inhibitors in single and dual pathway activated cell lines. We integrated hybrid-capture sequencing data from the Cancer Cell Line Encyclopedia with pharmacological data from over 150 compounds. We found that MAPK and PI3K pathway mutations most significantly associated with sensitivity to MEK and PI3K/AKT/mTOR inhibitors, respectively. Notably, cell lines with co-occurring MAPK pathway and *PIK*3CA mutations were insensitive to MEK inhibitors and cell lines with co-occurring PI3K pathway and KRAS mutations were insensitive to PI3K inhibitors. Also, not all pathway mutations conferred equal sensitivity. For example, BRAF mutant cell lines were generally the most sensitive sensitive, followed by NRAS and KRAS mutant cell lines, while NF1 and RASA1 mutant cell lines were generally not responsive to MEK inhibitors. Taken together, our work highlights the need to consider pathways and co-occurrence in the development of standard of care and targeted therapies.

Methods:

Pathway mutation status of TCGA patients: Mutation data was integrated from TCGA. Broad mutation packager data as of the 20121024 stddata build were included as well as all DCC level 2 mutation annotation format (MAF) files as of November 15, 2012. A significant number of likely false positive mutations were identified in prostate adenocarcinoma MAF files; therefore, all calls for this disease were sourced from Compendia Bioscience's own mutation calling pipeline. The Compendia Bioscience mutation calls were made to conform to the MAF file format for integration.

Mutation classifications: In certain diseases, such as uterine corpus endometrioid cancer (UCEC) and stomach adenocarcinoma (STAD), several highly-mutated samples may dominate the overall mutation counts. Thirty-one patients with >5,000 mutations (ultra-mutators) were excluded from downstream annotation analysis. A mutation was classified as 'Recurrent' if it was observed in the same variant position in 3 or more tumor samples. A mutation was classified as 'Deleterious' if it is non-recurrent and was annotated as a frame shift insertion/deletion or nonsense mutation variant. All other mutations were classified as 'Other'. Genes were classified as 'Gain of Function' if >20% of the mutation calls within the gene were recurrent missense mutation, less than 10% of the mutations calls were deleterious, and was significantly enriched with mutations. Genes were classified as 'Loss of Function' if they contained more than 2 deleterious mutations across all samples analyzed if >20% of mutation calls were hotspot or deleterious, and mutations were significantly enriched within the gene.

Cell line mutation and drug response data: Hybrid-capture gene sequencing data from the Cancer Cell Line Encyclopedia (CCLE) was integrated with drug response data from multiple large cell line screening panels including CCLE, Sanger cell line panel, GlaxoSmithKline (GSK) cell line panel, and OncoPanel 240 (OP240, Eurofins).

Conclusions:

- Pre-clinical models representative of clinical populations can potentially predict outcome to individual targeted therapy.
- In clinical samples, co-occurring activating mutations in MAPK and PI3K pathway genes were observed in >5% of patients in colorectal cancers, gastric cancers, endometrial cancers, melanoma, and cervical squamous cell carcinoma.
- Colorectal cancer patients with activating mutations in both PI3K and MAPK pathways had worse outcome than patients with MAPK pathway mutations alone.

Results:

whole exome analysis.







Figure 4. MAPK and PI3K pathway gene classifications. MAPK (A) and PI3K (B) pathway genes were classified by Compendia Bioscience as Gain of Function (red) and Loss of Function (blue).



mutations in 17 cancer types.



Cell Line Model

Only somatic hotspot missense and Function were aggregated at the patients were also positive for PI3K events were reported within that

Cell Line Model

Cell Line Model