

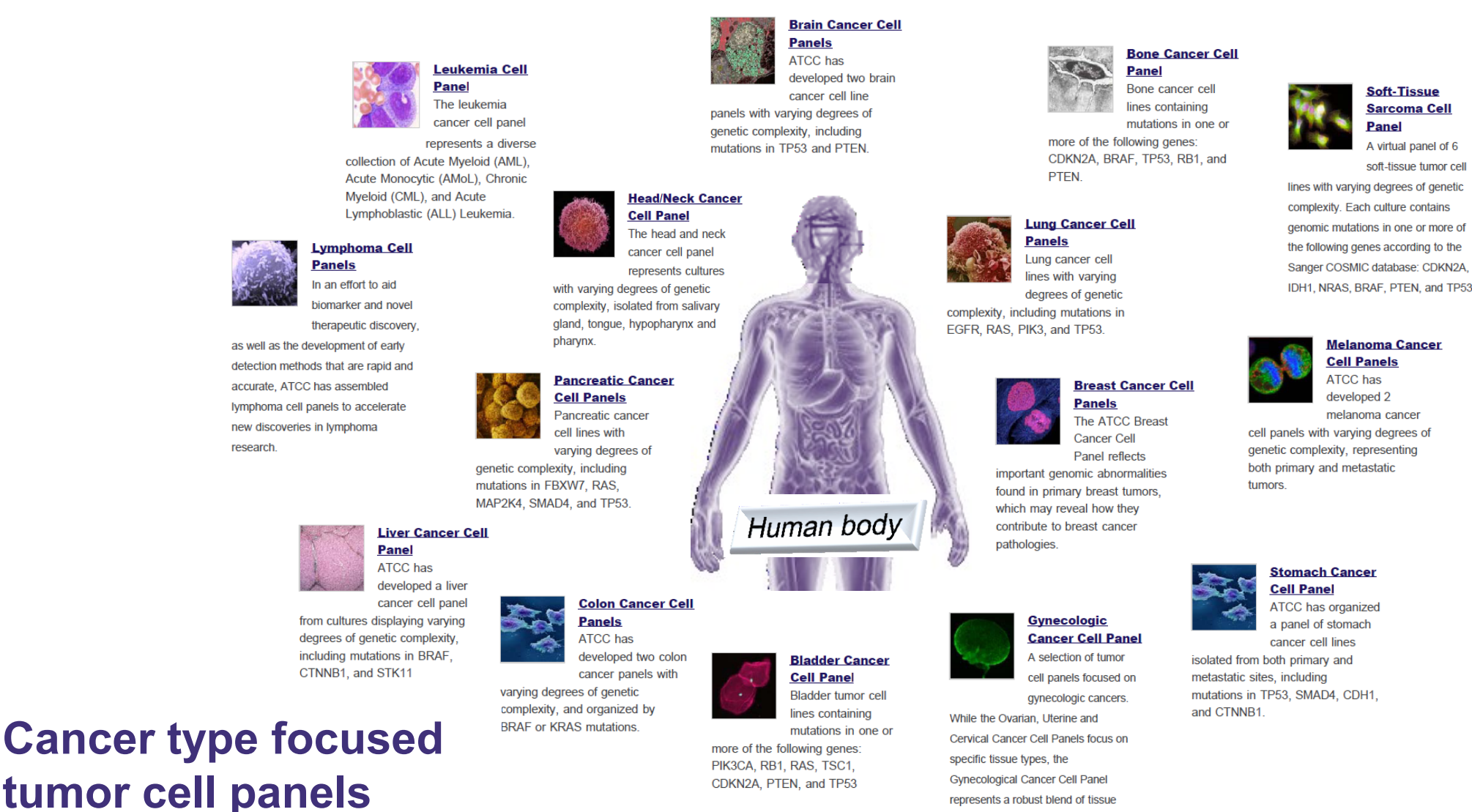
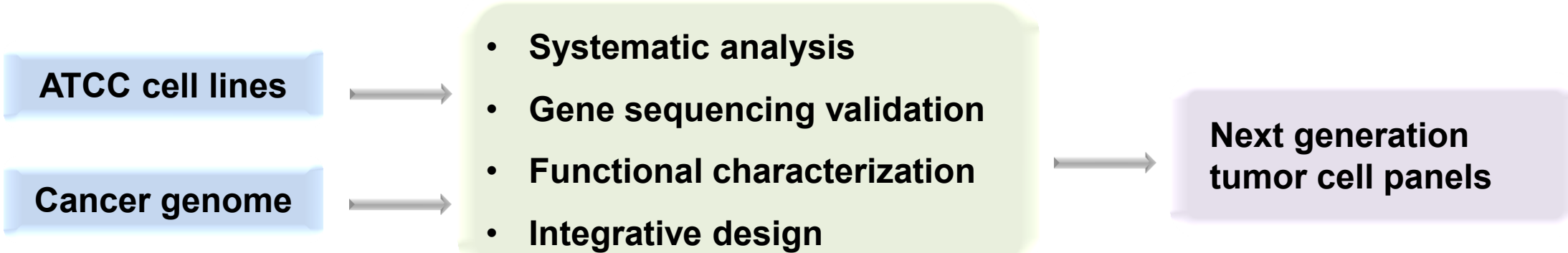
Abstract

Increased understanding of cancer genome is affecting every corner of cancer research. Although human tumor cell lines have been used as essential tools for decades, there are only a few cell line panels have been developed for the drug screening. There is a gap between the new knowledge of cancer genome and the cell line based platforms for both basic and translational research.

Here, we show that new generation tumor cell panels are filling the gap. The panels were generated by selecting authenticated cell lines derived from variant cancer types, and annotated with genetic alteration information generated by large scale sequencing projects such as the Catalog of Somatic Mutations in Cancer (COSMIC) and the Cancer Cell Line Encyclopedia (CCLE). To capture the genetic diversity of cancer, each panel includes cell lines with varying gene mutation complexity. To further facilitate targeted drug discovery, the molecular signature tumor cell line panels focus on individual driver genes, critical protein kinases, transcription factors and cell signaling pathways. Those panels have been analyzed to verify gene mutation, gene expression, protein expression and bio-functions.

Introduction

ATCC tumor cell panels program initiative



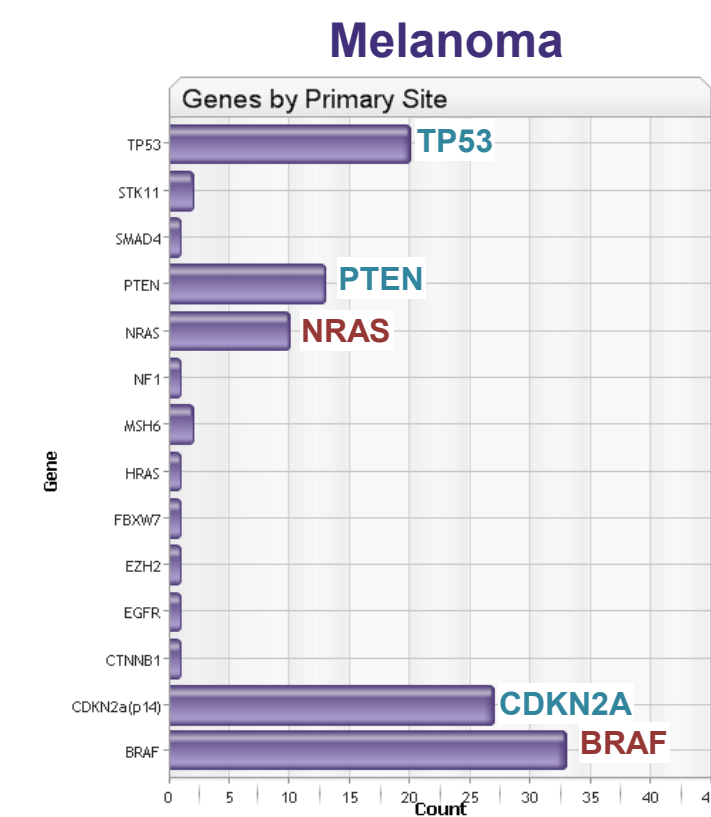
Cancer type focused tumor cell panels

The mutation data was obtained from the Sanger Institute Catalogue Of Somatic Mutations In Cancer database

Results

Bioinformatics analysis and clustering of ATCC cell lines

Diseased focused clustering



Mutation frequencies in melanoma cell lines

Metastatic melanoma cancer cell panel

Item No.	Name	Mutant	Zygoty	DNA Sequence	AA Sequence
HTB-69	SK-MEL-3	BRAF	heterozygous	c.1799T>A	p.V600E
		TP53	homozygous	c.799C>T	p.R267W
CRL-7724	SH-4	BRAF	homozygous	c.1799T>A	p.V600E
		CDKN2A	homozygous	c.1_471del471	p.0?
HTB-71	SK-MEL-24	BRAF	heterozygous	c.1799T>A	p.V600E
		CDKN2A	homozygous	c.1_471del471	p.0?
		PTEN	homozygous	c.80_164del85	p.?
HTB-66	RPMI-7951	BRAF	heterozygous	c.1799T>A	p.V600E
		CDKN2A	homozygous	c.47T>G	p.L16R
		PTEN	homozygous	c.1_79del79	p.?
		TP53	homozygous	c.497C>A	p.S166*

Gene focused clustering

PIK3CA mutation

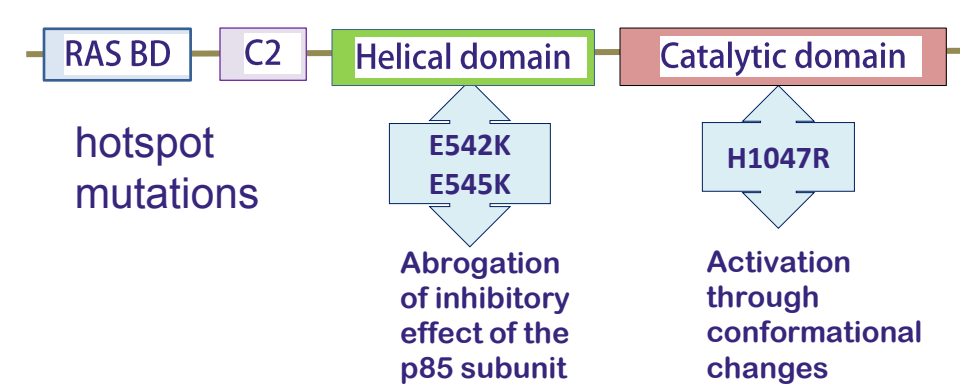
Mutant	frequency
p.E545K	28%
p.E545D	3%
p.H1047R	28%
p.H1047L	3%
p.E542K	5%
p.R88Q	5%
p.K111E	3%
p.K111N	3%
p.K111R	3%
p.P539R	3%
p.Q546R	3%
p.D549N	3%
p.E453K	3%
p.G118D	3%
p.P124L	3%
p.P449T	3%
p.G106_R108del	3%

Tissue source

Breast
Caecum
Cervix
Colon
Lung
Lymphoid
Ovary
Pharynx
Prostate
Stomach
Urinary bladder
Uterus

Cell lines: 36

PIK3CA somatic mutation



Cell death pathway BCL-2 family cell panels

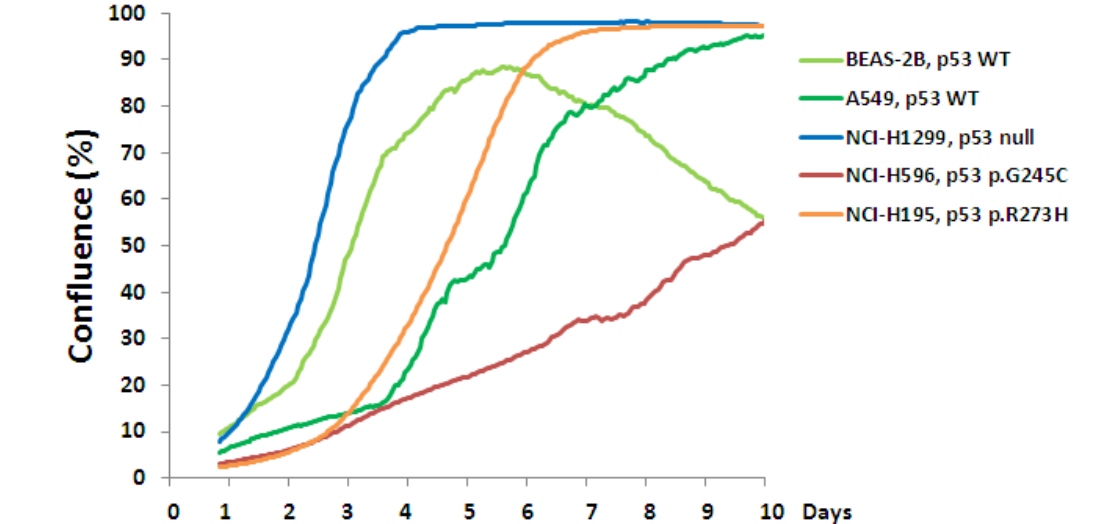
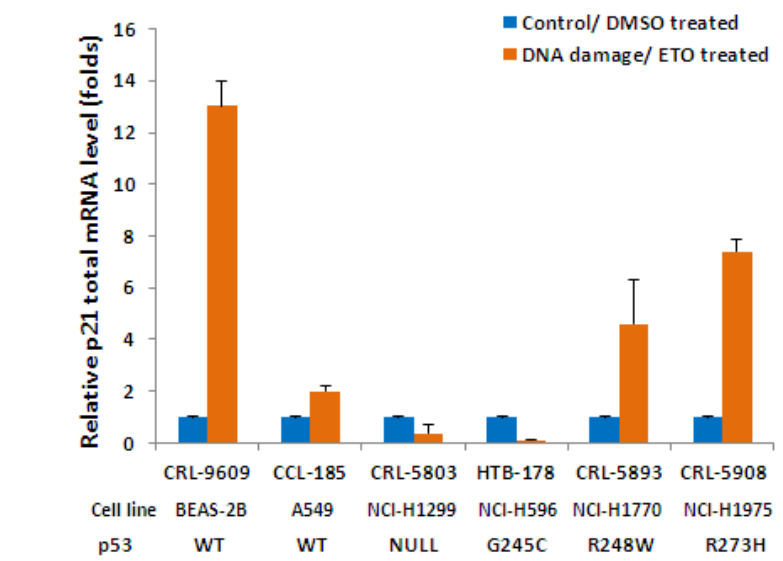
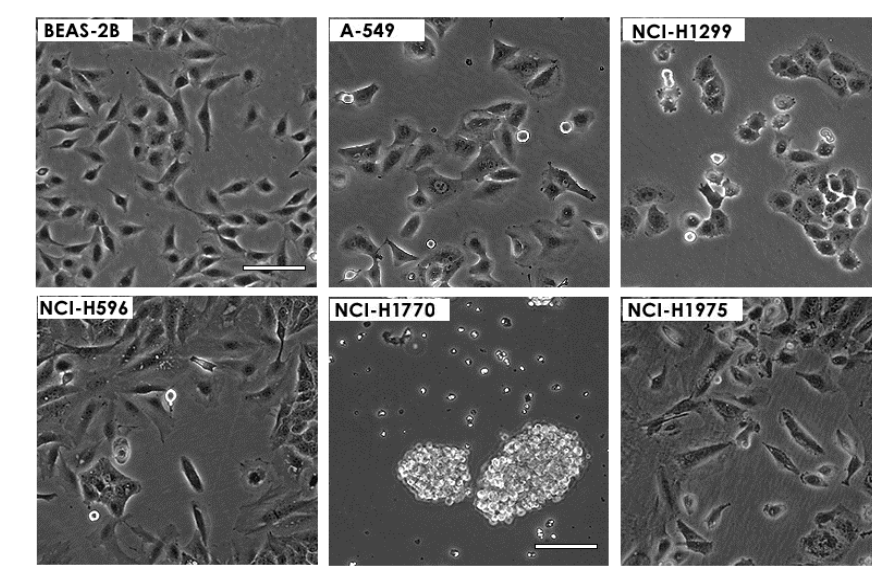
Item No.	Cell line name	Expressing /Signature
CRL-2898	Neo Jurkat	empty vector
CRL-2899	BCL2 Jurkat	Bcl-2
CRL-2900	BCL2 (S70A) Jurkat	phosphorylation deficient mutants of Bcl-2 (S70A)
CRL-2901	BCL2 (S87A) Jurkat	phosphorylation deficient mutants of Bcl-2 (S87A)
CRL-2902	BCL2 (AAA) Jurkat	multi-site T69A/S70A/S87A (AAA) mutants of Bcl-2

Item No.	Cell line name	Derived from
CRL-2907	WT SV40 MEF	WT mouse
CRL-2908	Bcl2 KO SV40 MEF	Bcl2 KO mouse
CRL-2909	Bad KO SV40 MEF	Bad KO mouse
CRL-2910	Bax KO SV40 MEF	Bax KO mouse
CRL-2911	Bid KO SV40 MEF	Bid KO mouse
CRL-2912	Bak KO SV40 MEF	Bak KO mouse
CRL-2913	Bax Bak DKO SV40 MEF	Bax and Bak double KO mouse

Molecular signature tumor cell panels

Non-small cell lung cancer p53 hotspot mutation tumor cell lines

ATCC® No.	Name	Tissue	Histology	Tumor Source	TP53 status	Zygoty	CDS mutation	AA mutation
CRL-9609™	BEAS-2B	lung	normal tissue,SV-40 immortalized	NA	WT	-	-	-
CCL-185™	A549	lung	non-small cell lung carcinoma	primary	WT	-	-	-
CRL-5803™	NCI-H1299	lung	non-small cell lung carcinoma	metastasis (lymph node)	NULL	homozygous	c.(del)	-
HTB-178™	NCI-H596	lung	adenosquamous carcinoma	primary	MUT	homozygous	c.733G>T	p.G245C
CRL-5893™	NCI-H1770	lung	non-small cell lung carcinoma	metastasis (lymph node)	MUT	homozygous	c.741 742CC>TT	p.R248W
CRL-5908™	NCI-H1975	lung	adenocarcinoma	primary	MUT	homozygous	c.818G>A	p.R273H

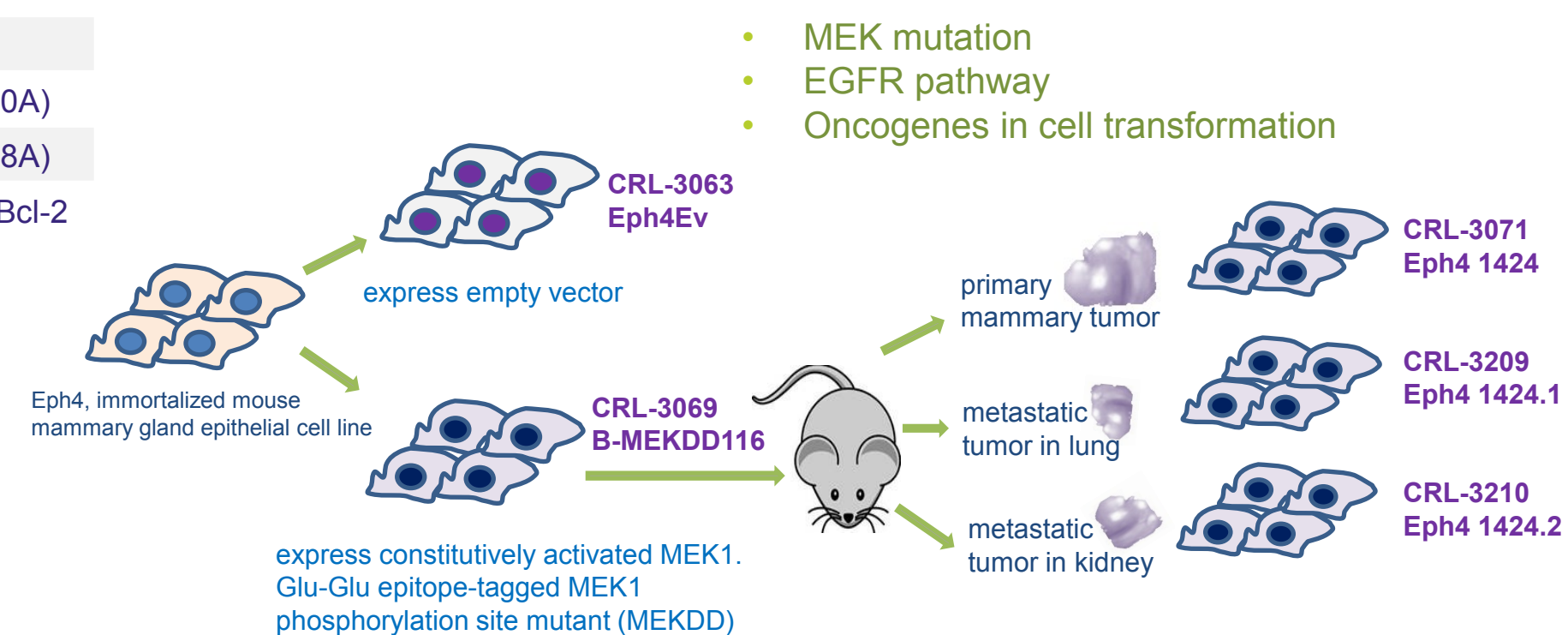


Seven p53 hotspot mutation panels include cell lines

- p53 wild type
- p53 null
- p53 hotspot codon 175 mutation
- p53 hotspot codon 248 mutation
- p53 hotspot codon 273 mutation
- p53 hotspot codon 282 mutation
- Other p53 hotspot: codon 220, 249

- Breast cancer cells
- Colon cancer cells
- NSCLC cells
- SCLC cells
- Pancreatic cancer cells
- Leukemia cells
- Lymphoma cells

Breast cancer mouse model cell panel



Summary

- ATCC is harnessing the combined forces of genomic data and our highly reliable, authenticated cell lines to generate useful tools for cancer research
- Focused on disease, driver genes and signaling pathways, various cancer cell panels and normal cell controls facilitate both basic and translational research

Reference

- Cancer Cell (2002) 2, 29-42
- Oncogene (1998) 16, 737 - 746
- Science (2001) 292, 727-730
- Nature (2010) 463, 899-905
- Nature reviews Cancer (2010) 10, 241-253
- Nature reviews drug discovery (2009) 8, 627-644