

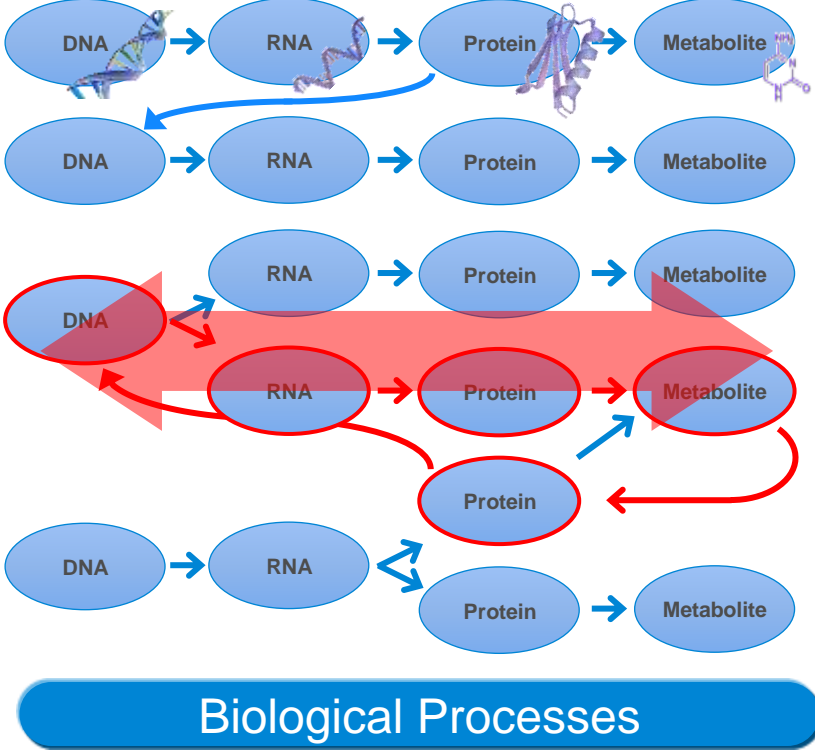
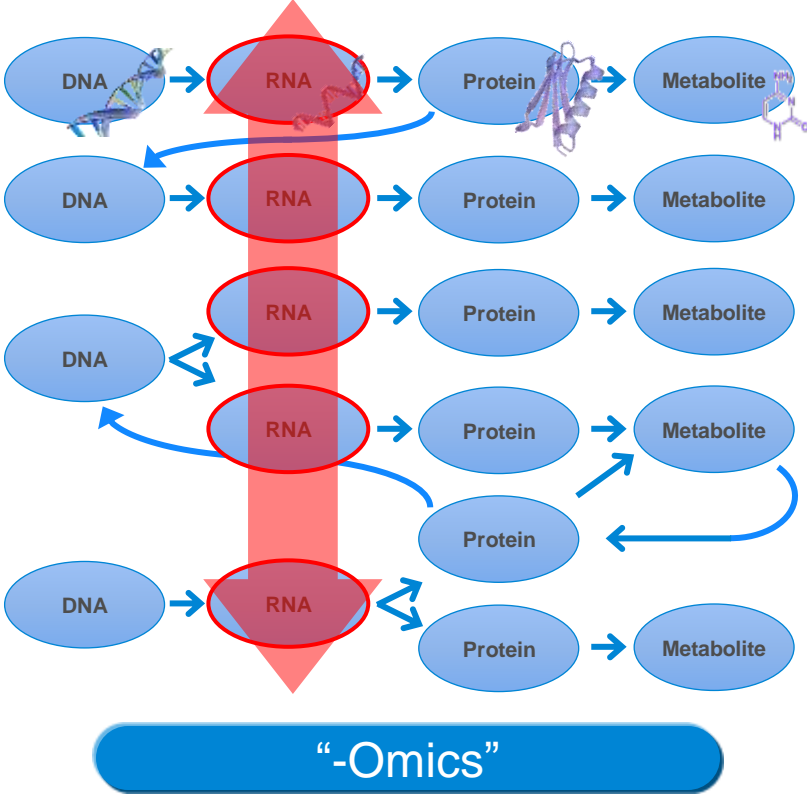


Extracting new insights from cancer genomics and proteomics data

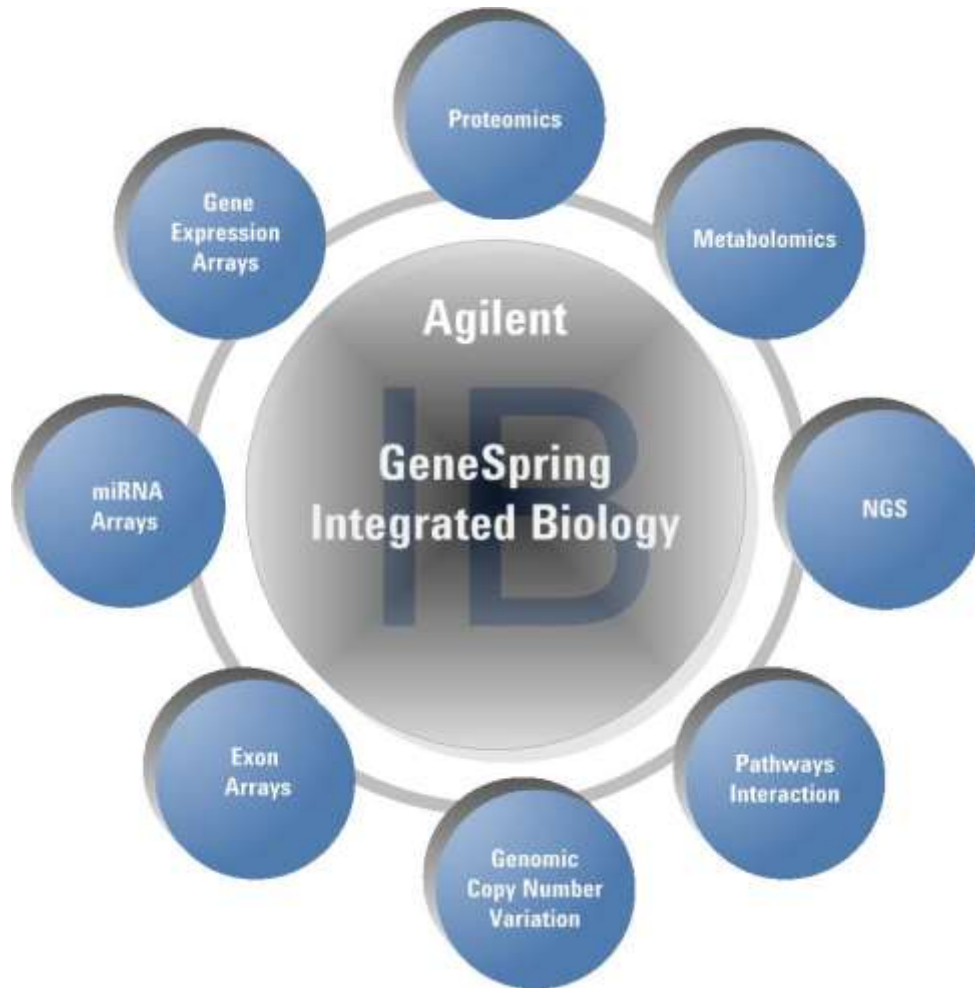
Shweta Shukradas
Product Manager

For Research Use Only. Not for use in diagnostic procedures

The Cross Technology Challenge



GeneSpring Suite of Products



GeneSpring GX

- mRNA Expression
- microRNA Expression
- Genome-wide Association
- Copy Number Variation
- Variant Analysis with vcf
- qPCR

Mass Profiler Professional

- Proteomics
- Metabolomics

Strand NGS

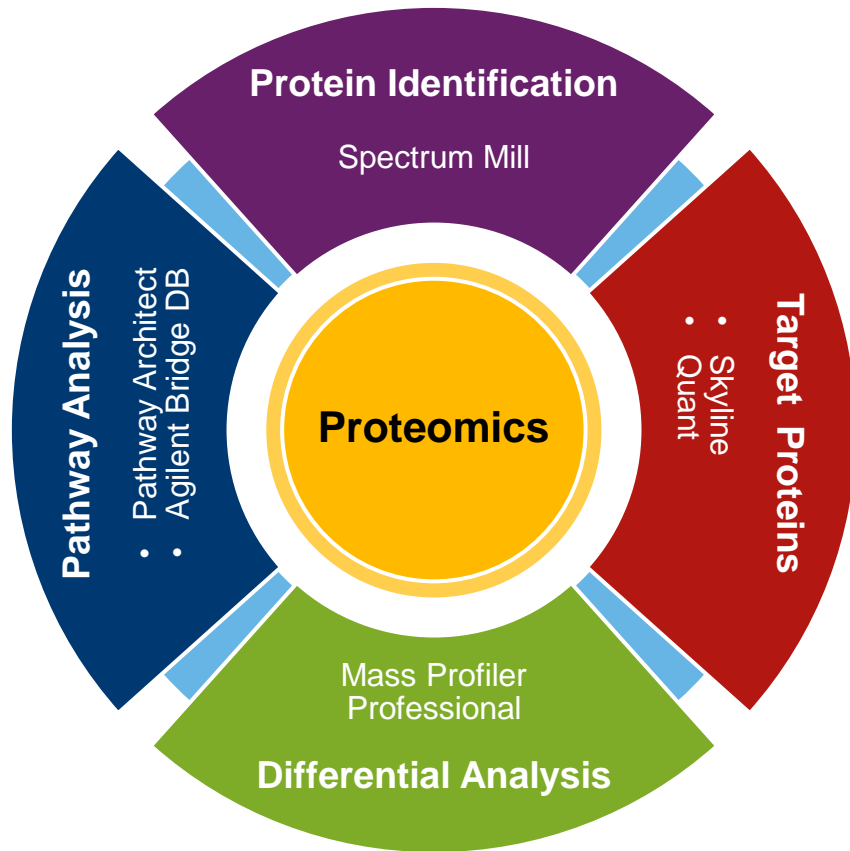
- DNA-Seq, RNA-Seq, ChIP-Seq,
- Methyl-Seq and small RNA-Seq

Pathway Architect

- Multi-omic Pathway Analysis
- NLP discovery

PROTEOMICS WORKFLOW IN MPP

Most Extensive Software Portfolio for Proteomics Data Analysis



Identify proteins from discovery data-dependent analysis

and/or

Target peptides for proteins of interest

Analyze data for statistically different features

Search and map annotated proteins to find interesting pathways

Visualize results on pathways

Agilent's Discovery Proteomics Workflow

Separate and Detect

6500 Series LC/Q-TOF



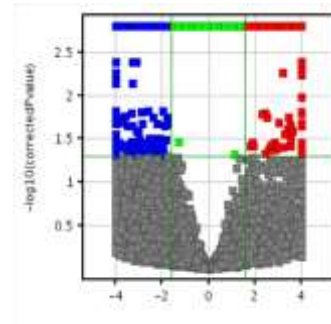
Identification and Quantitation

Spectrum Mill



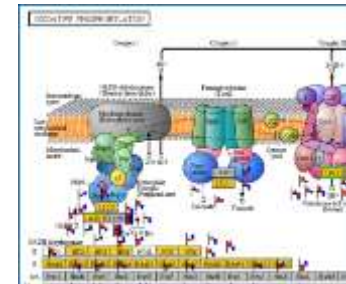
Statistical Analysis

Mass Profiler Professional



Pathway Analysis

Pathway Architect



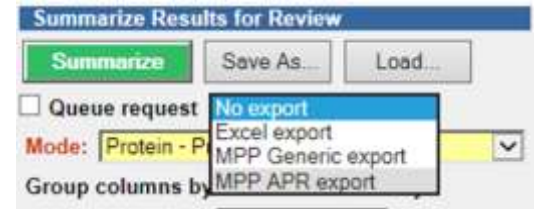
Pathway-directed targeted protein(s) experiment



Agilent Technologies

For Research Use Only. Not for use in diagnostic procedures

Two MPP Export Modes in Spectrum Mill



Generic

- Support all data
- Protein abundance information only



Group (#)	Subgroup (#)	Spectra (#)	Distinct Peptides (#)	Distinct Summed MS/MS Search Score	% AA Coverage	Total Protein Spectral Intensity	Database Accession #	Protein Name
2	2.1	1377	34	518.95	87	6.13e+009	P43490	Nicotinamide phosphoribosyl-transferase

#	Filename	#	Score	SPI (%)	Spectrum Intensity	Sequence	Precursor Average Charge Squared	m/z Measured (Da)
1	251T-L-000183428.83603.2	2	25.95	85.8	3.59e+007	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1079.5573
2	188T-L-000158456.88481.2	2	25.64	85.7	3.01e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1087.5513
3	195T-L-000158432.88630.2	2	25.52	85.4	3.02e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1087.5513
4	245T-L-000138829.39180.2	2	25.22	100.0	1.38e+007	(K)YLLETSMHLDGLEYK (L)	1.00	857.9312
5	202T-L-000160579.60758.2	2	25.01	85.4	4.45e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1087.5538
6	251T-L-000182702.83129.2	2	24.97	100.0	3.32e+007	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1079.5507
7	202T-L-000159780.59947.2	2	24.86	84.8	4.02e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1087.5535
8	245T-L-000138502.38928.2	2	24.83	81.2	1.43e+007	(K)YLLETSMHLDGLEYK (L)	0.97	857.9308
9	245T-L-000154982.55241.2	2	24.72	86.9	2.87e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1087.5543
10	251T-L-000183014.84438.2	2	24.02	87.0	3.88e+007	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1079.5574
11	245T-L-000153430.86602.2	2	24.00	82.8	4.70e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.99	1079.5574
12	202T-L-000170029.70092.2	2	23.78	84.7	4.56e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.97	1079.5564
13	234T-L-000167545.57911.2	2	23.75	87.0	4.48e+008	(R)VIQGGVDINTLQETVEGR (Q)	0.97	1087.5518
14	234T-L-000141804.42286.2	2	23.73	86.7	1.68e+007	(K)YLLETSMHLDGLEYK (L)	0.98	857.9308
15	234T-L-000142026.43130.2	2	23.58	87.0	1.89e+007	(K)YLLETSMHLDGLEYK (L)	0.98	857.9308
16	245T-L-000152087.52287.2	2	23.28	80.2	1.87e+008	(R)VIQGGVDINTLQETVEGR (Q)	1.00	1087.5547

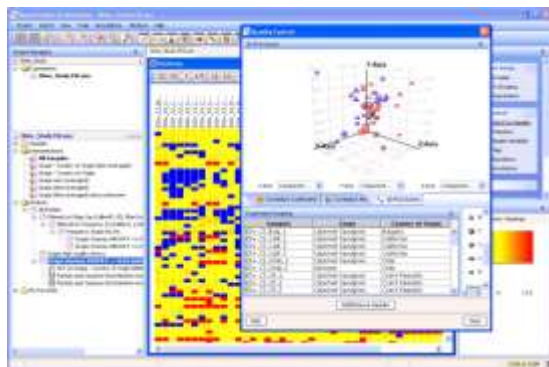
Agilent Proteomics Results (.apr)

- Support only Agilent data
- Protein abundance and peptide abundance included in export
- Supports label-free, SILAC, iTRAQ, TMT
- Requires MPP 14.0 release

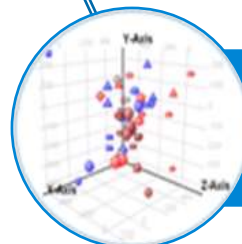


Mass Profiler Professional

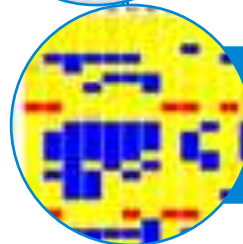
Find differences in mass spec data sets and reach statistically valid conclusions



Statistical Analysis



Get from data to answers using statistical tests such as PCA



Dive deeper into results using visualizations e.g. heat maps, volcano plots



Generate target lists for Q-TOF and QQQ to do targeted proteomics



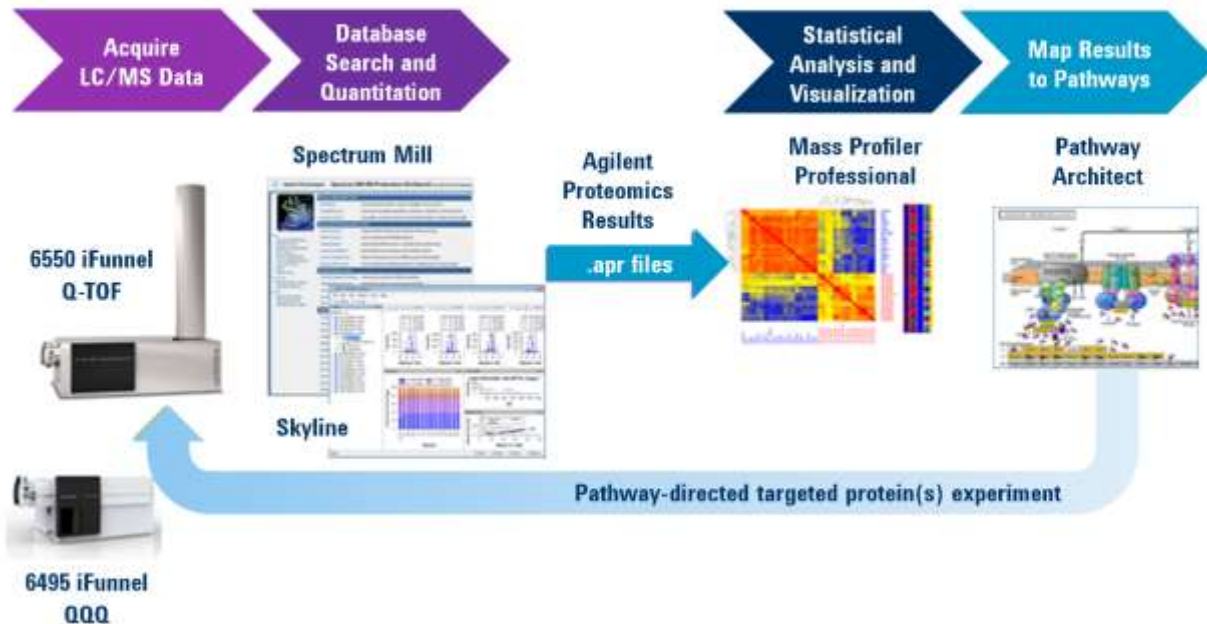
Bring genomics, proteomics, and metabolomics together



Agilent Technologies

For Research Use Only. Not for use in diagnostic procedures

Proteomics Workflow in MPP



- Designed for new .apr format export with protein and peptide-level information
- Filters available for peptide-level and protein-level (abundance, frequency, PTMs)
- Protein/Peptide Entity Inspector

Protein List Inspector

Protein List Inspector

Name: All Entries
 Notes: All Entries

Creation date: Tue Feb 07 11:51:06 PST 2017
 Last modified date: Tue Feb 07 11:51:07 PST 2017
 Owner: grouer
 Technology: SpectrumML IDENTIFIED PROTEINS toxome_2016_Jan_20_17_34_18
 Number of proteins: 3999
 Experiments: toxome

Entities Attributes

Protein	Database	NCBI ID	Protein Name	Species	Swiss-Prot ID
Q95JF1	SwissProt		Dynamin family chain B, axonal	HUMAN	Q95JF1
Q92820	SwissProt		Gamma-glutamyl hydrolase	HUMAN	Q92820
P55145	SwissProt		Mesencephalic astrocyte-deri...	HUMAN	P55145
Q14318	SwissProt		Peptidyl-prolyl cis-trans isom...	HUMAN	Q14318
Q96008	SwissProt		Mitochondrial import receptor	HUMAN	Q96008
P09012	SwissProt		U1 small nuclear ribonucleop...	HUMAN	P09012
P05161	SwissProt		Ubiquitin-like protein E3C15	HUMAN	P05161
Q9V445	SwissProt		Transformation/transcription	HUMAN	Q9V445
Q13425	SwissProt		Beta-2-sitrophen	HUMAN	Q13425
Q9H1E3	SwissProt		Nuclear ubiquituous casein an...	HUMAN	Q9H1E3
Q5JFE7	SwissProt		Nodal modulator 2	HUMAN	Q5JFE7
Q13492	SwissProt		Phosphatidylinositol-binding	HUMAN	Q13492
P78344	SwissProt		Eukaryotic translation initiato...	HUMAN	P78344
Q03135	SwissProt		Caveolin-1	HUMAN	Q03135
P68400	SwissProt		Casein kinase II subunit alpha	HUMAN	P68400
Q15427	SwissProt		Splicing factor 38 subunit 4	HUMAN	Q15427
P28340	SwissProt		DNA polymerase delta catalytic subunit	HUMAN	P28340
Q94826	SwissProt		Mitochondrial import receptor subunit TOM70	HUMAN	Q94826
Q9NV59	SwissProt		Pyridoxine-5'-phosphate oxidase	HUMAN	Q9NV59
Q9HCY8	SwissProt		Protein S100-A14	HUMAN	Q9HCY8

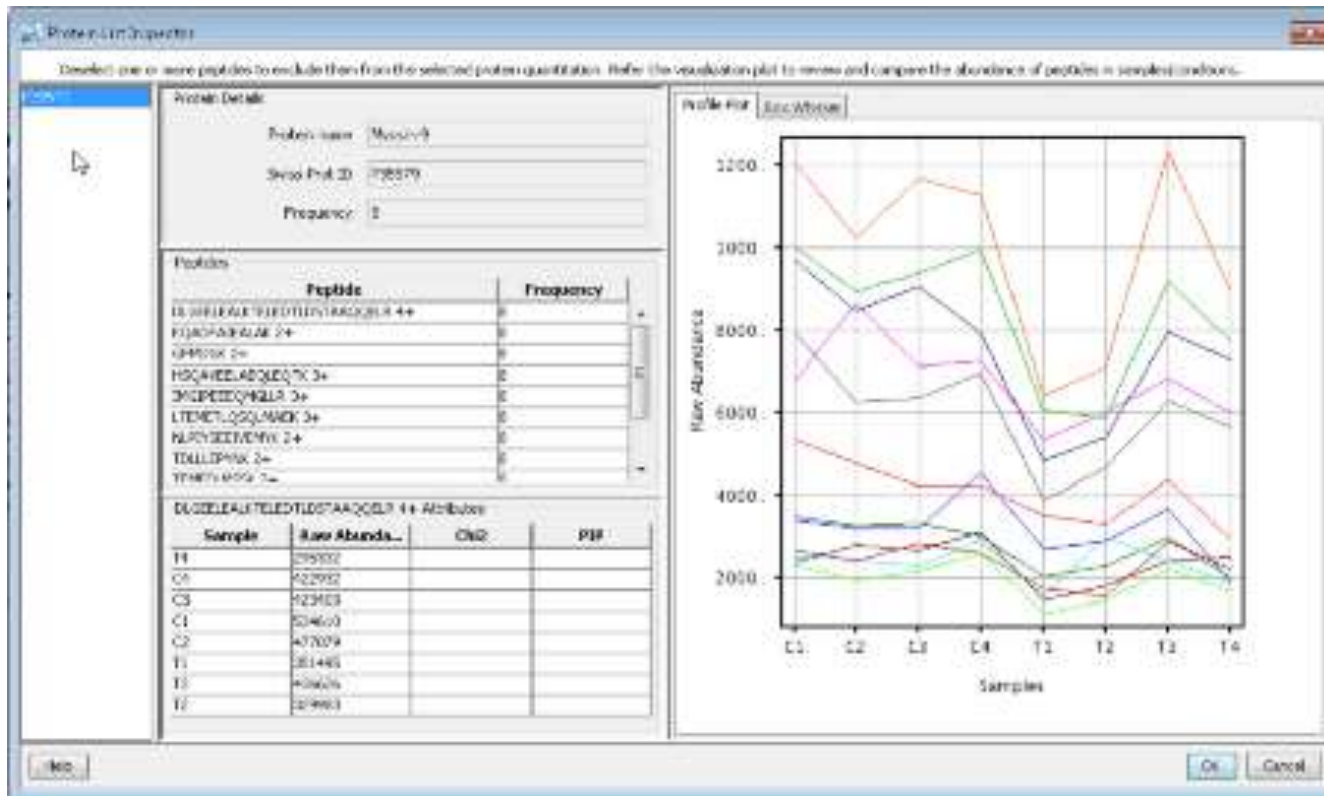
Find: Q03135 Find Next Find Previous Match Case

Help

Protein	Protein Name	Swiss-Prot ID	Species
Q13492	Phosphatidylinositol-binding clathrin assembly protein	Q13492 Q60641	HUMAN
P78344	Eukaryotic translation initiation factor 4 gamma 2	P78344	HUMAN
Q03135	Caveolin-1	P66539 Q03135	HUMAN
P68400	Casein kinase II subunit alpha	P68400	HUMAN
Q15427	Splicing factor 38 subunit 4	Q15427	HUMAN
P28340	DNA polymerase delta catalytic subunit	P28340	HUMAN
Q94826	Mitochondrial import receptor subunit TOM70	Q94826	HUMAN
Q9NV59	Pyridoxine-5'-phosphate oxidase	Q9NV59	HUMAN
Q9HCY8	Protein S100-A14	Q9HCY8	HUMAN

Find: Q03135 Find Next Find Previous Match Case

Protein Inspector Visualization



- Works at **both the protein and the peptide level simultaneously**
- This workflow design truly answers the biologist's questions of **"what's changing in my proteins?"** as well as **"what's changing at the peptide level?"**

Protein Inspector

Review peptide level information

Protein Inspector

Select a protein to review the corresponding peptides and their attributes for the chosen entity list and interpretation. Refer the visualization plot to review and compare the abundance of peptides in samples/conditions. Protein not included in chosen entity list or missing data in the chosen interpretation will be grayed.

Inspect selected protein in list: All Entities Select interpretation: Group (Non-averaged)

Protein	Protein Name	Frequency
Q03135	Caveolin-1	4

Find: Find Next Find Previous Match Case

Peptides

Peptide	Charge	Frequency	Alt. Accession	Vmod	Mod	Mass
AMADELSEK 2+	2	4				993.456
EIDLVRDPK 2+	2	3	P56539			1,198.643
HLNDDVVK 2+	2	4				939.489
IDFEDVIAEPEGTHSF...	3	4				2,405.13
IFSNWR 2+	2	3				735.415
YVDSEGHLYTVPIR 3+	3	4				1,648.833
EIDLVRDPK 3+	3	2	P56539			1,198.643
qYDAHTK 2+	2	2		Q40q	q:Pyroglutamic acid	961.474

EIDLVRDPK 2+ Attributes

Sample	Flag	Raw Abundance	Chi2	PIP
C1	P	178,715	1	55.9
C2	A			
T1	P	102,668	0.99	62.3
T2	P	135,962	1	71.8

Peptides: Profile plot Peptides: BoxWhisker plot Protein: Profile plot Protein: BoxWhisker plot

Group (Non-averaged)

Help OK Cancel

Protein Inspector Visualization

Protein Inspector

Select a protein to review the corresponding peptides and their attributes for the chosen entity list and interpretation. Refer the visualization plot to review and compare the abundance of peptides in samples/conditions. Protein not included in chosen entity list or missing data in the chosen interpretation will be grayed.

Inspect selected protein in list: **Moderated T-Test p ([Control] vs [Tumor]) out-off = 0.05** Select interpretation: **Group (Non-averaged)**

Protein	Protein Name	Frequency	p (Contr...	Log FC	FC (abs)	Regulation	FC	Abund Diff ...	Abund Dif...
Q8IY67	Ribonucleoprotein PTB-binding 1	4	4.95E-02	1.13	2.19	up	2.19	111668.32	16.77
Q86UQ4	ATP-binding cassette sub-family A member 13	4	3.74E-02	1.46	2.75	up	2.75	361821.97	18.46
P63151	Serine/threonine-protein phosphatase 2A 55 kDa regulatory s...	4	4.27E-02	1.48	2.80	up	2.80	122410.20	16.90
Q98UH6	Uncharacterized protein C9orf142	4	1.78E-02	1.62	3.07	up	3.07	71737.02	16.13
Q9H187	Interferon regulatory factor 2-binding protein-like	4	4.77E-02	1.43	2.70	up	2.70	180707.22	17.46
Q86U86	Protein polybromo-1	4	1.89E-02	-2.30	4.94	down	-4.94	-478220.22	-18.87
Q6UXG2	UPF0577 protein KIAA1324	4	3.04E-02	2.56	5.89	up	5.89	280704.00	18.00
P98179	Putative RNA-binding protein 3	4	4.32E-02	1.03	2.04	up	2.04	251621.00	14.00
Q02166	DNA damage-binding protein 2	4	2.57E-02	1.72	2.22	up	2.22	60603.26	11.26

Find: Find Next Find Previous Match Case

Peptides

Peptide	Charge	Freq...	Alt. Acc...	Vmod	Mod	Mass
LLGDLLEAVR 2+	2	4				1,195.704
NAEALAESESRLR 2+	2	2				1,402.717
TVLSAK 2+	2	2				689.419
VFAPDAVSKPGMDEYELK 3+	3	4				1,916.946
YGLSAAAAAAAAAAVEQKR 3+	3	4				1,803.935
EKQRNAEALAESESRLR 3+	3	1				1,944.015

LLGDLLEAVR 2+ Attributes

Sample	Flag	Raw Abundance	Chi2	PIP
T1	P	93,890	0.96	52.8
C1	P	243,138	0.96	54.4
T2	P	149,262	0.95	60.4
C2	P	356,079	0.97	50.4

Peptides: Profile plot Peptides: BoxWhisker plot Protein: Profile plot Protein: BoxWhisker plot

Raw Abundance

Group (Non-averaged)

Control: C1, C2 Tumor: T1, T2

Help OK Cancel

Filtering on Peptides

Filter on peptides
Select an entity list and interpretation to filter peptides based on Frequency, Sample Variability, Abundance, Modifications, Charge, and Properties.
To apply the filter, click on Preview button.

Entity List: All Entities Choose...
Interpretation: Group (Non-averaged) Choose...

Frequency Sample Variability Abundance **Modifications** Charge Properties

Filter by Modifications

Retain peptides with atleast one of the selected modifications

Allowed Disallowed

Oxidized Methionine Allowed Disallowed

Pyroglutamic acid(only at N-terminus of peptide) Allowed Disallowed

Phosphorylated Threonine Allowed Disallowed

Deamidated Asparagine Allowed Disallowed

Phosphorylated Serine Allowed Disallowed

Phosphorylated Tyrosine Allowed Disallowed

Select/Deselect all modifications

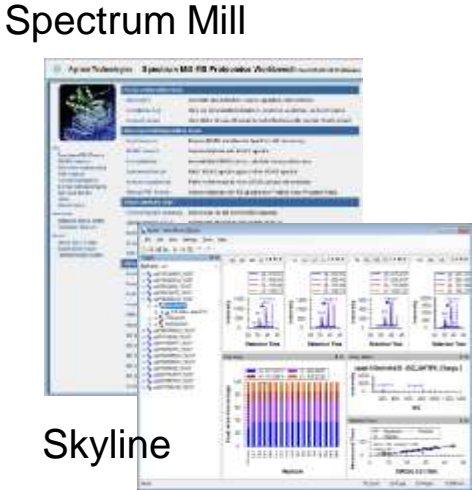
Include corresponding unmodified peptide

Preview

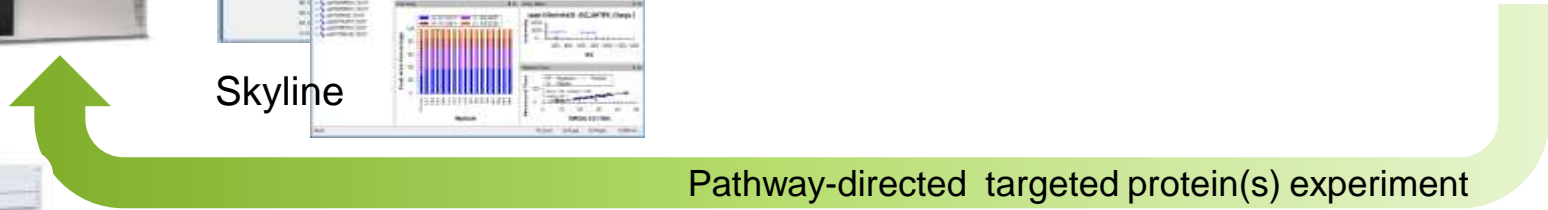
Help OK Cancel

- Modifications filter at the level of peptides
- The filter dialog lists all the available peptide modifications that are present in the selected entity list
- Filtering results in a list of proteins that only consist of specific peptides selected as allowed

Accelerate Your Research With a Complete Pathway Centric Workflow



6495 iFunnel QQQ



GENOMICS WORKFLOWS IN GENESPRING GX

Genomics Research Solutions



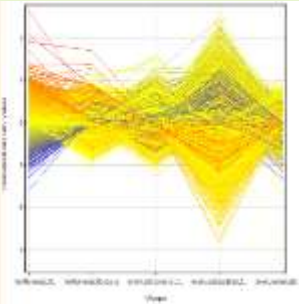
Gene Expression Analysis Workflow

Filtering

Statistical Analysis

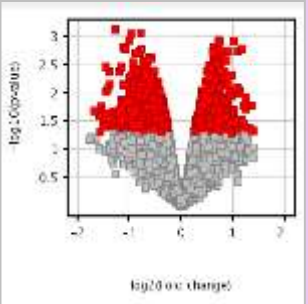
Pathway Analysis

Filtering Probe sets on expression



Probe sets filtered based on the percentile criteria

Statistical analysis



Performed to find differentially expressed probes

Pathway Analysis



Enrichment test performed with pathways of interest

Clustering - Hierarchical with preview window

Clustering Analysis

Select an entity list, interpretation, and a clustering algorithm. The available parameters change according to the selected clustering algorithm and whether you choose Normalized intensity values or Associated values for the hierarchical clustering algorithm. When you have selected all the parameters, click Preview. The clustering results are displayed in the results pane on the right. To save your results in the experiment, click OK.

Entity List: 13 entities E2 and PPT col
Interpretation: PPT dose - Time Point - In
Clustering Algorithm: Hierarchical

Parameters

- Normalized intensity values
- Associated values
- Cluster on: Both entities and conditions
- Distance metric: Eukclidean
- Linkage rule: Ward's
- Cluster within conditions

Preview

Heatmap visualization showing hierarchical clustering results. The top panel shows a dendrogram and a heatmap with a color scale from 0 to 30. The bottom panel shows a more detailed heatmap with a dendrogram on the left and labels for conditions: PPR, EGRC, RABD, GREE, FHL2, FHL2, MYB, SGK1, GRHL, HSPE, MYC, CXCL, and NC41. The rows are labeled with entity IDs such as NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_J1, NC_MCF7_J1, NC_MCF7_J1, NC_MCF7_J1, NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_B1, NC_MCF7_J1, NC_MCF7_J1, NC_MCF7_J1, NC_MCF7_J1.

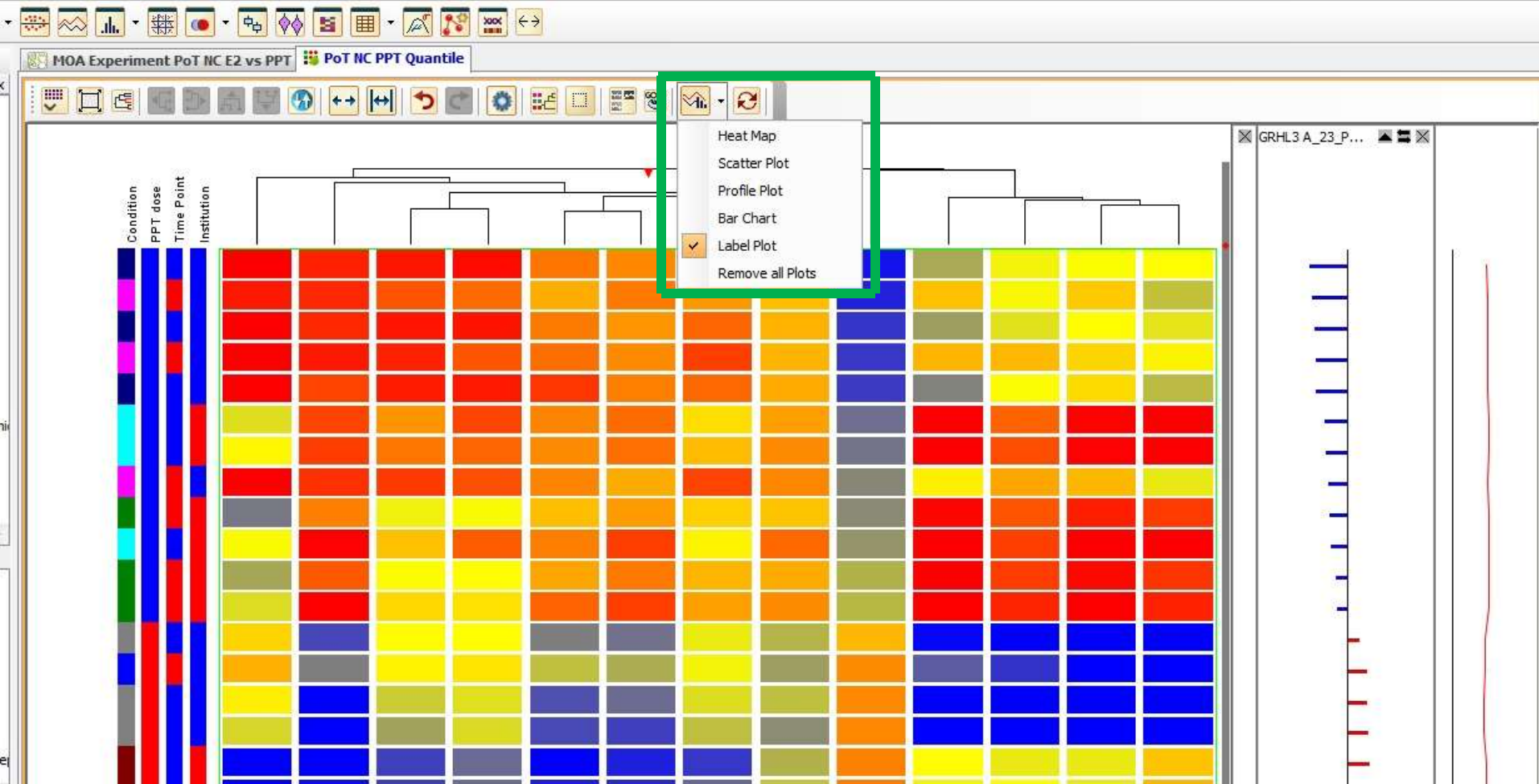
Workflow

- Experiment Setup**
 - Quick Start Guide
 - Experiment Grouping
 - Create Interpretation
 - Create New Gene-level Experiment
- Quality Control**
 - Quality Control on Samples
 - Filter Probesets by Expression
 - Filter Probesets by Flags
 - Filter Probesets on Data Files
 - Filter Probesets by Error
- Analysis**
 - Statistical Analysis
 - Filter on Volcano Plot
 - Find Similar Entities
 - Filter on Parameters
 - Principal Component Analysis
- Class Prediction**
- Results Interpretations**

Legend

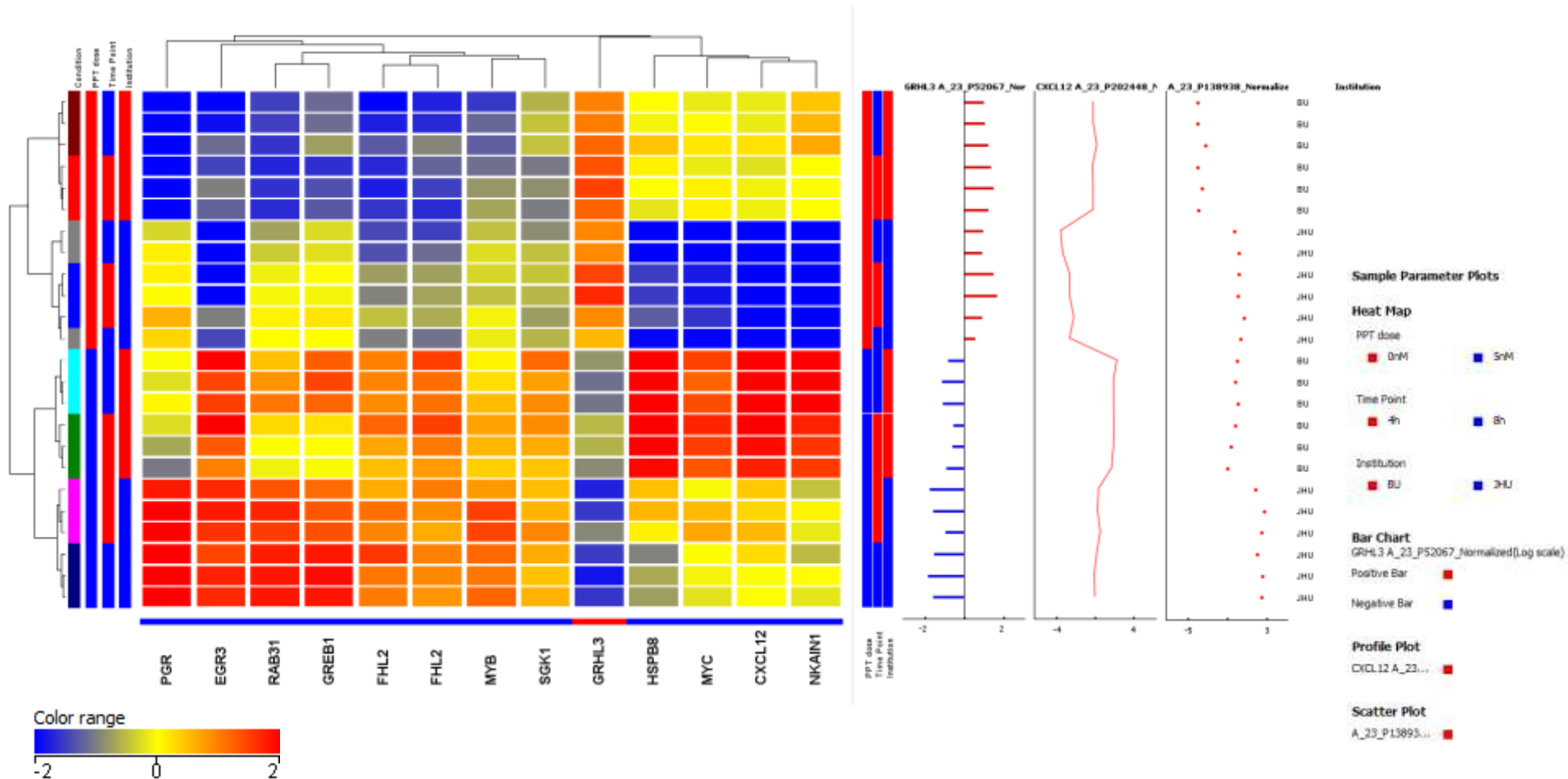
Legend - BoxWhisker Plot

Metadata Framework: adding plots to clustered heatmap

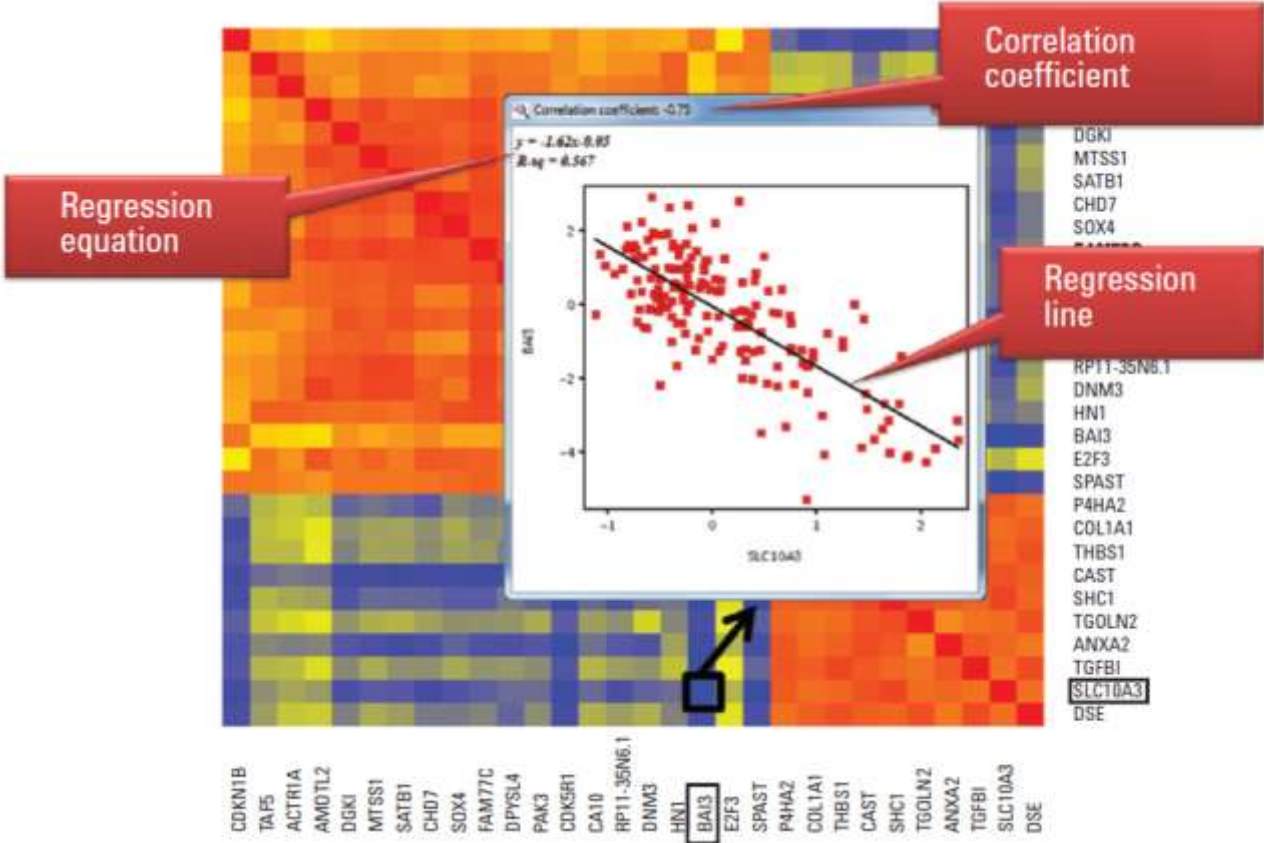


Metadata Framework

From Experimental Grouping Table information



Correlation Analysis



Researchers can view relationships between entities (compounds or proteins) or samples. Clicking on a cell of a heat map, they can quickly view the specific parameters of the correlation

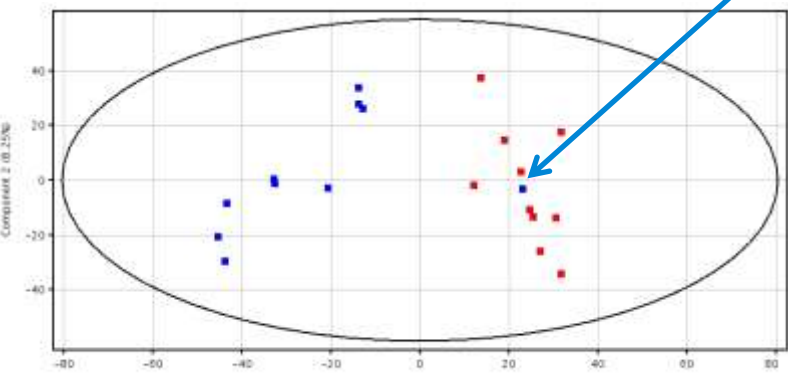
INTEGRATED ANALYSIS

Genomics and Proteomics

Brain Tumor Study - Quality Control of Discovery Proteomics Data

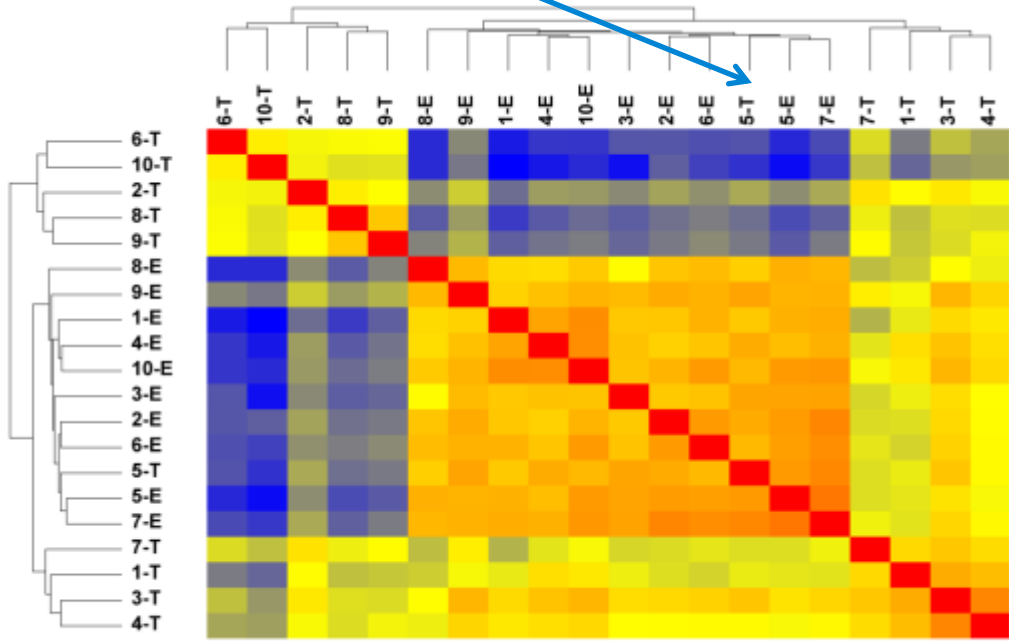
Sample 5-T

Principle Component Analysis Plot



■ control
■ tumor

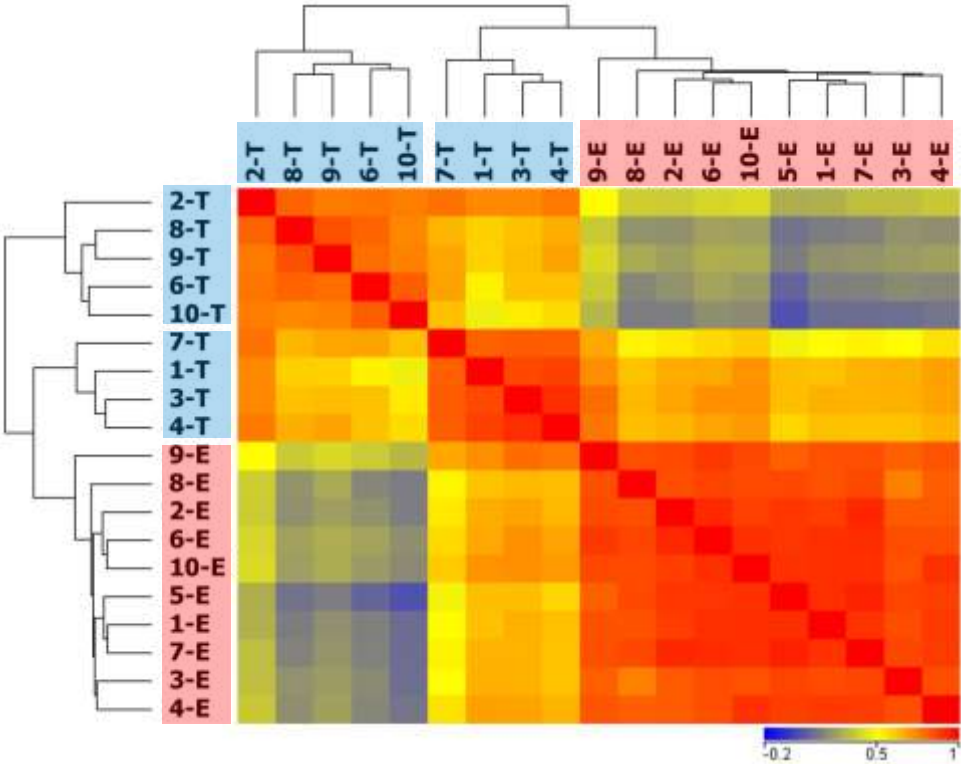
Sample Clustered Correlation Plot



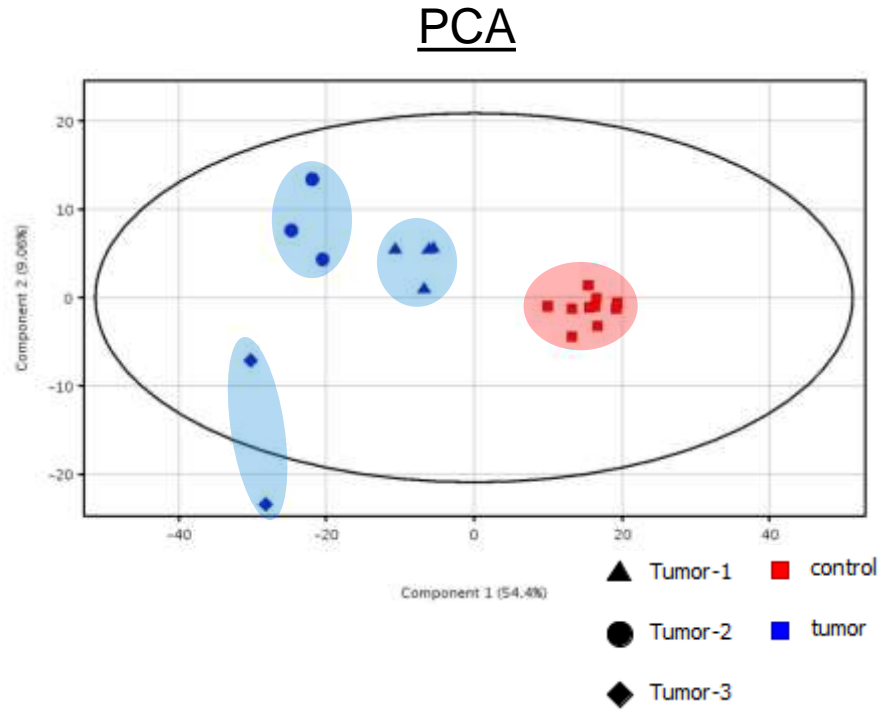
T – Tumor samples E – Epilepsy samples

Brain Tumor Study - Sample-sample Correlation and PCA of Discovery Proteomics Data

Tumor subgroups demonstrated, using **587 differential proteins**, in both PCA and sample correlation showing the sample heterogeneity within tumors

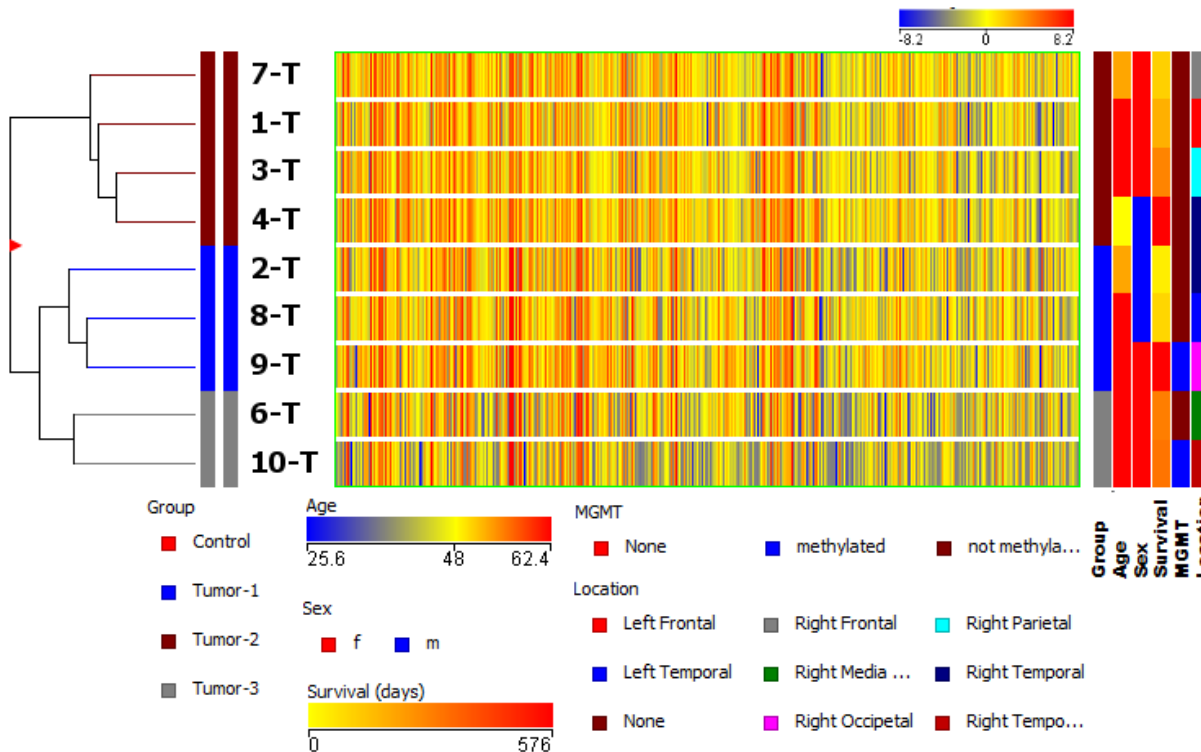


Sample-Sample Correlation



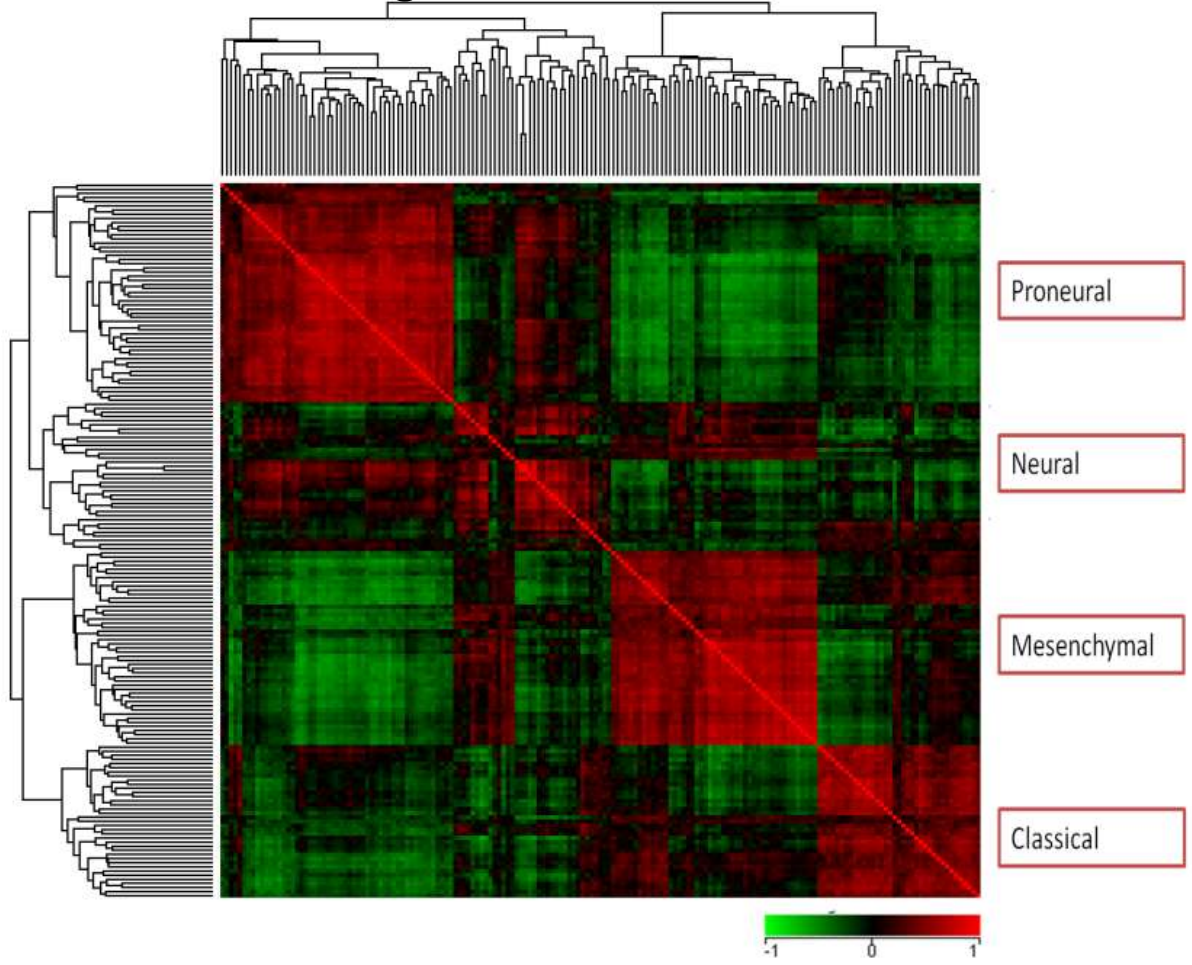
PCA

Brain Tumor Study - Metadata Analysis of Discovery Proteomics Data



- Hierarchical Clustering of samples provided similar classification as PCA
- Tumor subgroups aligned to sample metadata such as age, sex, tumor location and methylation status of MGMT
- The metadata did not explain the sub-grouping observed within the tumors based on the differentially expressed proteins

Brain Tumor Study - Hierarchical Clustering of Genomic Signatures

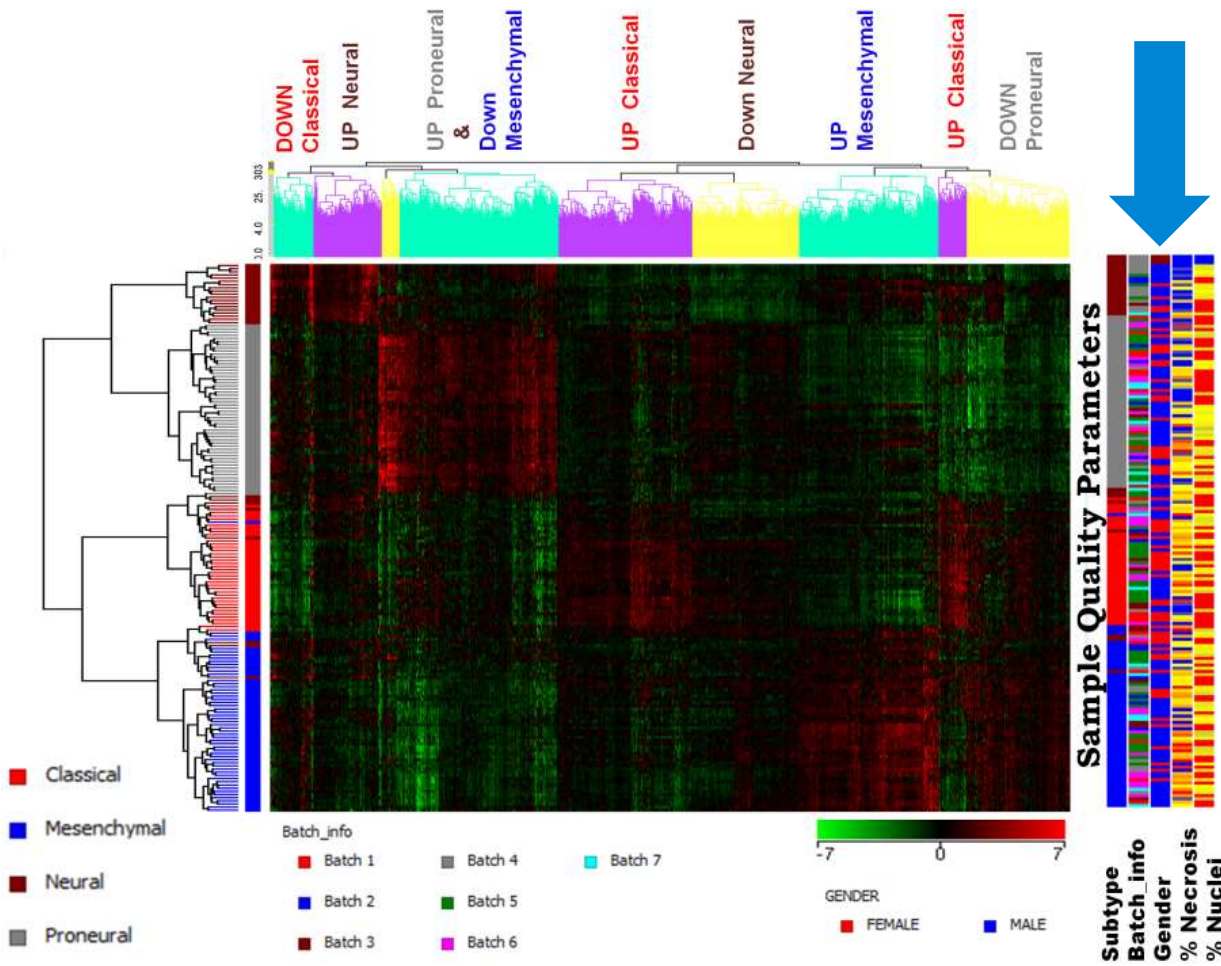


Hierarchical clustering of 840 mRNA profiles from 173 GBM tumors

The four GBM subtypes were re-created

Cancer Cell. 2010 January 19; 17(1): 98

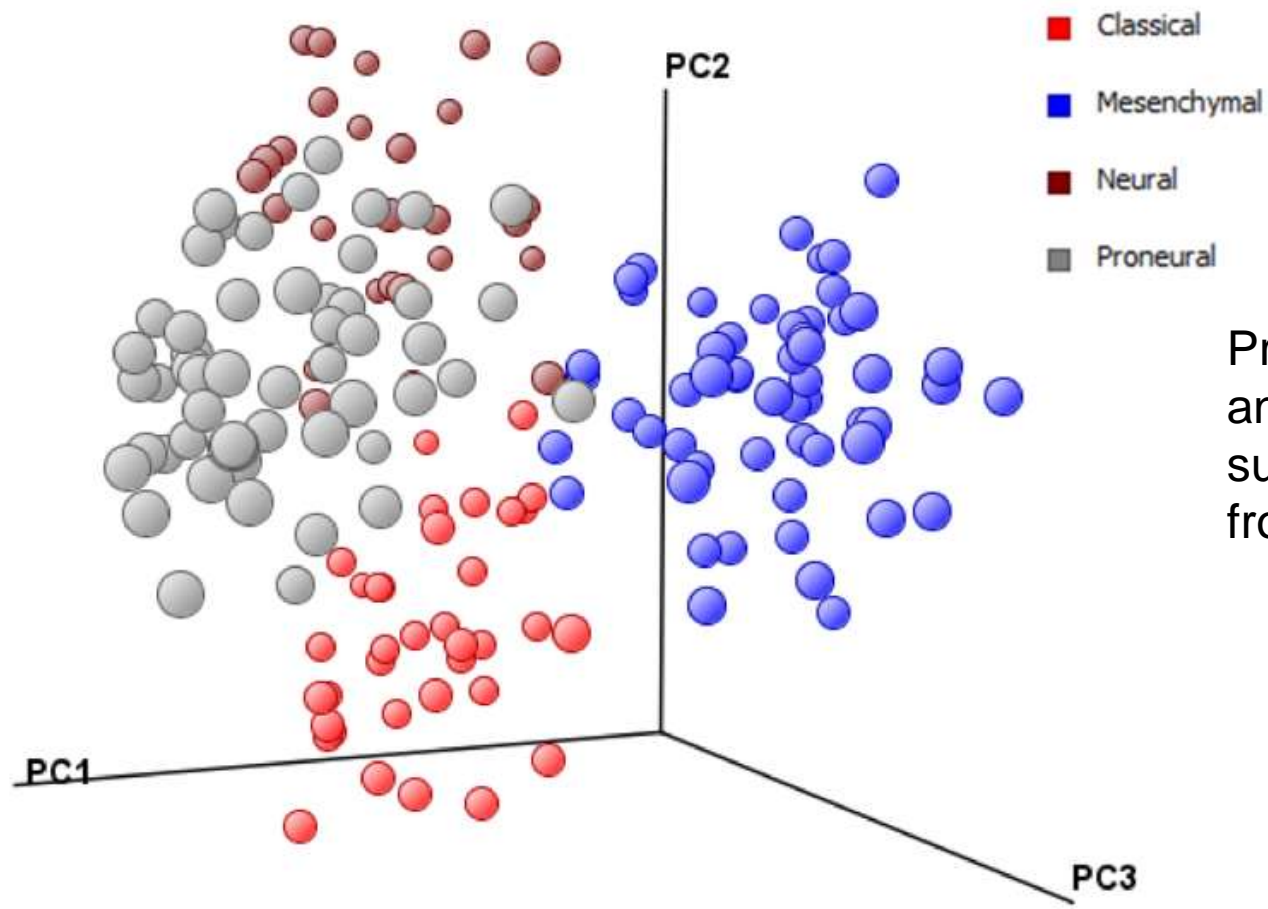
Brain Tumor Study - Metadata Analysis of Genomic Signatures



Metadata obtained from TCGA was imported into the project

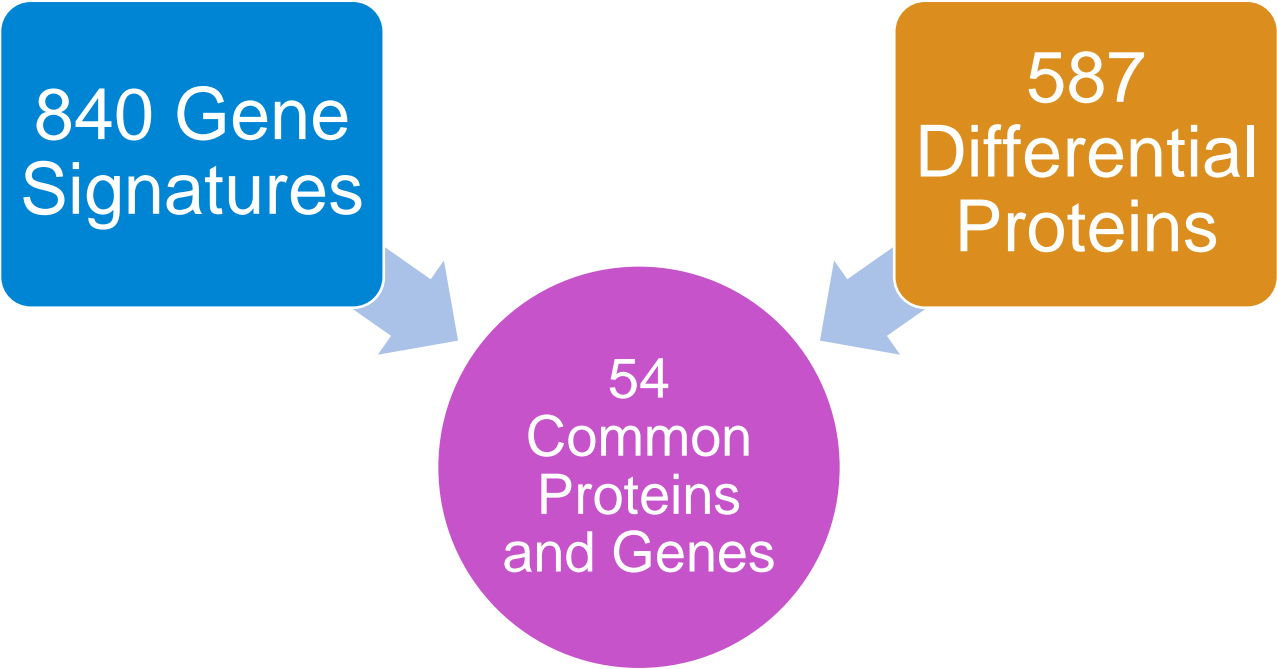
Metadata framework analysis of batch and sex was used to assess batch affects or bias

Brain Tumor Study - Principal Component Analysis of Genomic Signatures



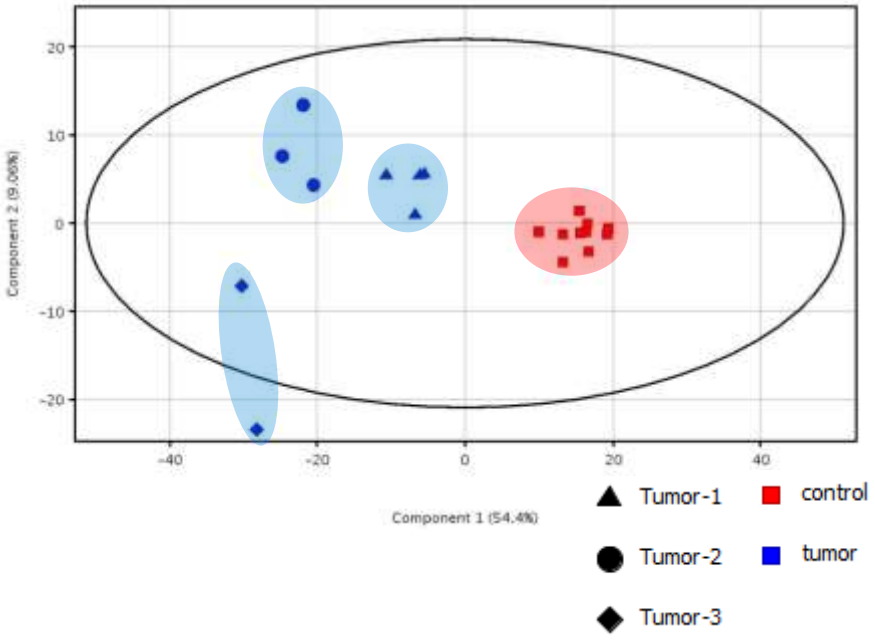
Principal component analysis of confirms subtypes in GBM samples from TCGA

Brain Tumor Study - Venn Analysis of Genomic and Proteomic Data Set

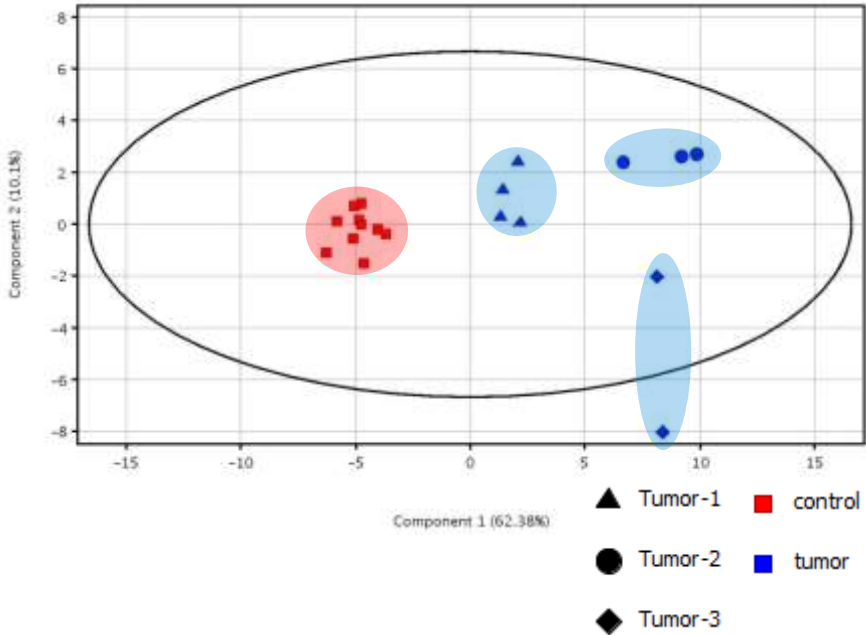


Brain Tumor Study - PCA Comparison of Original 587 Proteins to 54 Proteins Identified Using Integration

587 Differential Proteins



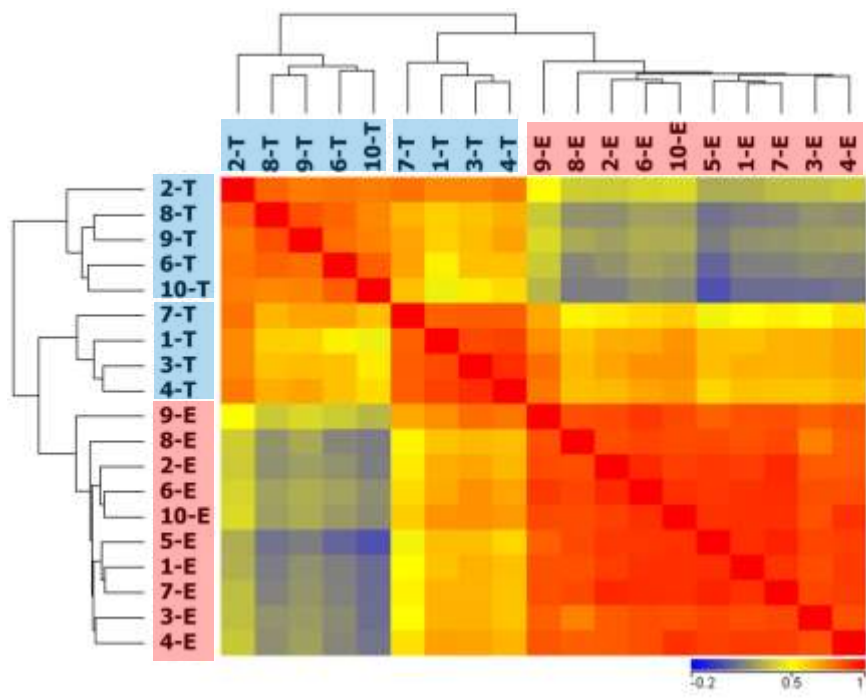
54 Proteins from Integration



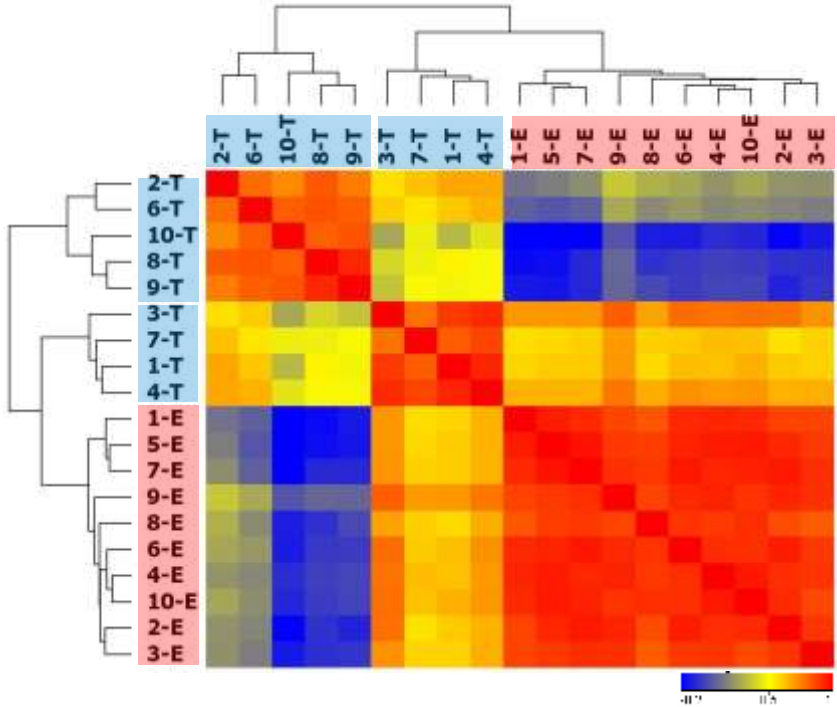
54 of 587 proteins that overlap the genomics signatures were good enough to capture the sub grouping in tumor samples, indicating the relevance between the molecular subtypes from genomics data and the subgroup pattern in proteomics data

Brain Tumor Study - Comparison of 587 Differential Proteins to 54 Proteins Identified Using Integration

587 Differential Proteins

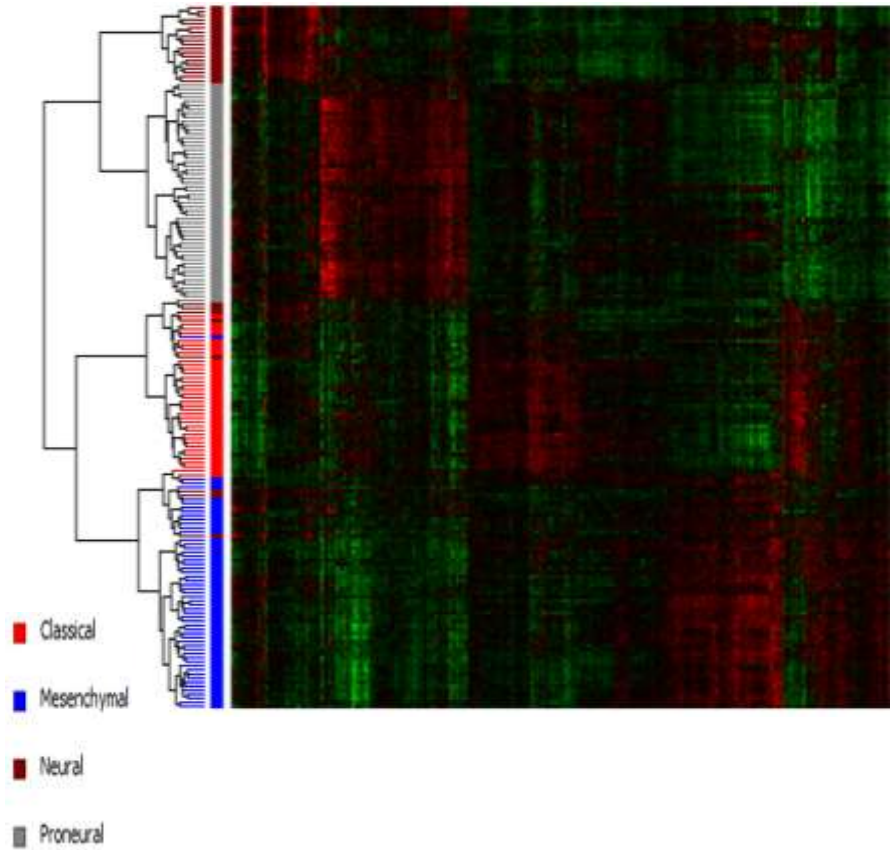


54 Proteins from Integration

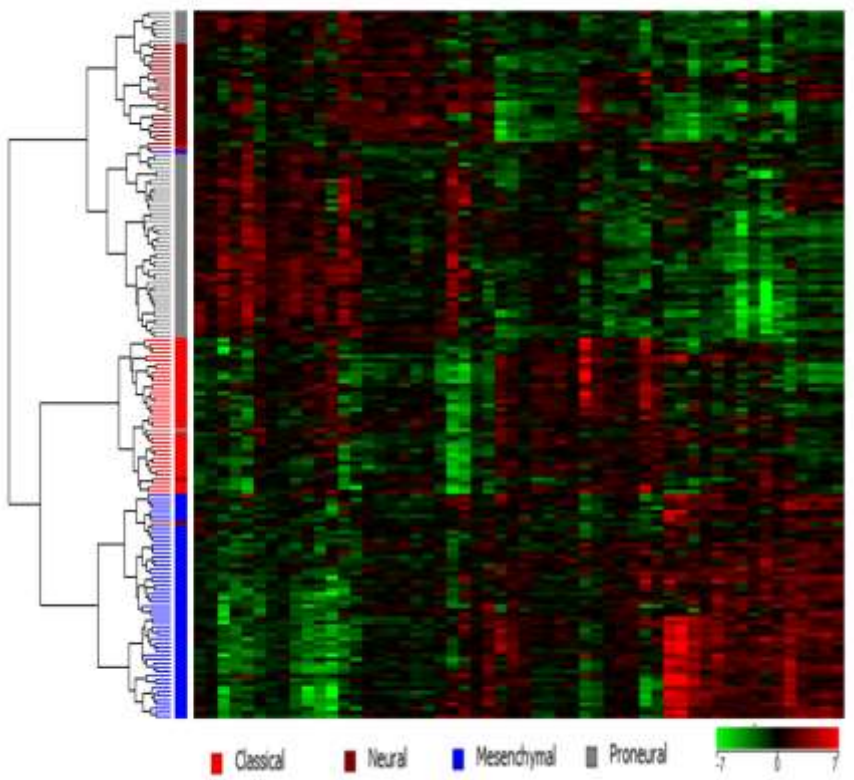


Brain Tumor Study - Performance of 54 Genes Identified Using Proteomics Integration to Original 840 Signature

Original Genomic 840 Signature



Genomic-Proteomic 54 Signature



Brain Tumor Study : Summary

Proteomics

- Use sample correlation & PCA on discovery proteomics data to identify outlier sample
- Identify a set of differential proteins for the tumor samples
- Use correlation and PCA to identify subgroups within tumor samples
- Clinical metadata visualized to reason the sub grouping within tumor samples

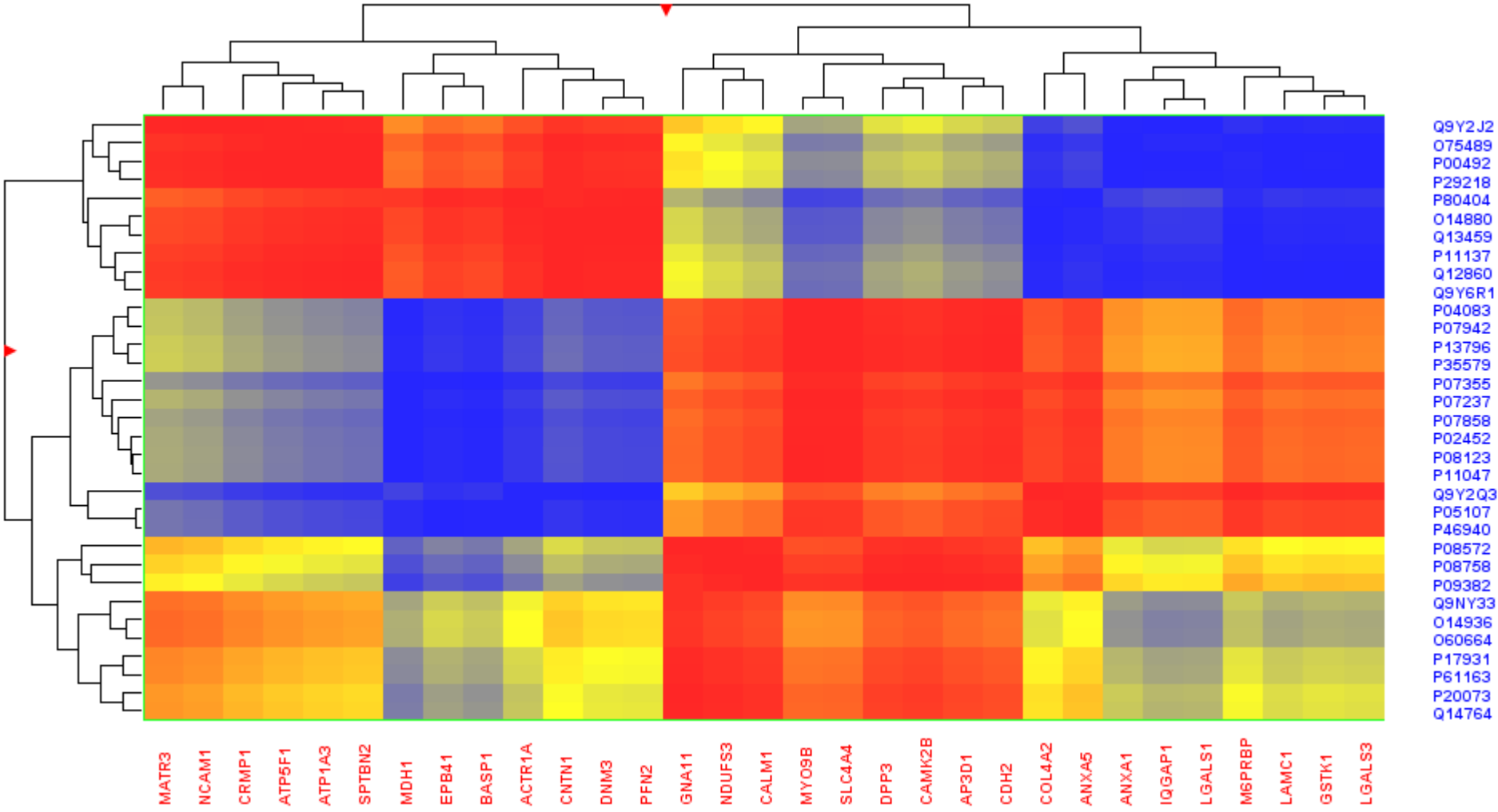
Genomics

- Use clustering and PCA on mRNA profiles to confirm subtype specific genomic signatures
- Assess data quality and bias using meta data analysis framework

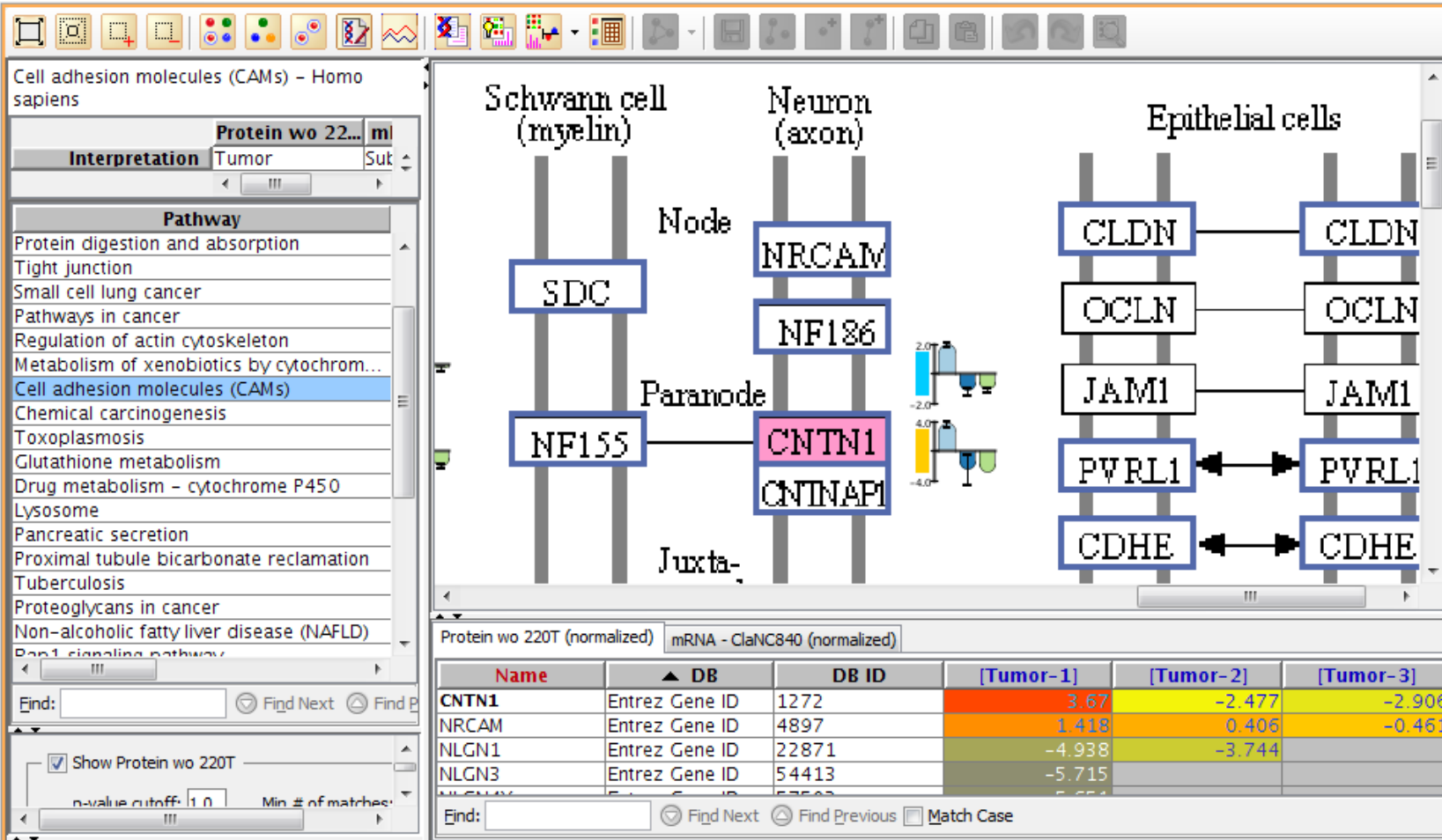
Integration

- Map gene id to protein accession number
- Perform Venn analysis of differential proteins and gene signatures
- Reduced set of gene signatures based on proteomics classified tumors in larger TCGA cohort

MOA Correlation Analysis - Concordant Genes and Proteins



Common Pathways - Concordant Genes and Proteins



Proteomics

- [Tumor-1]
- [Tumor-2]
- [Tumor-3]

Genomics

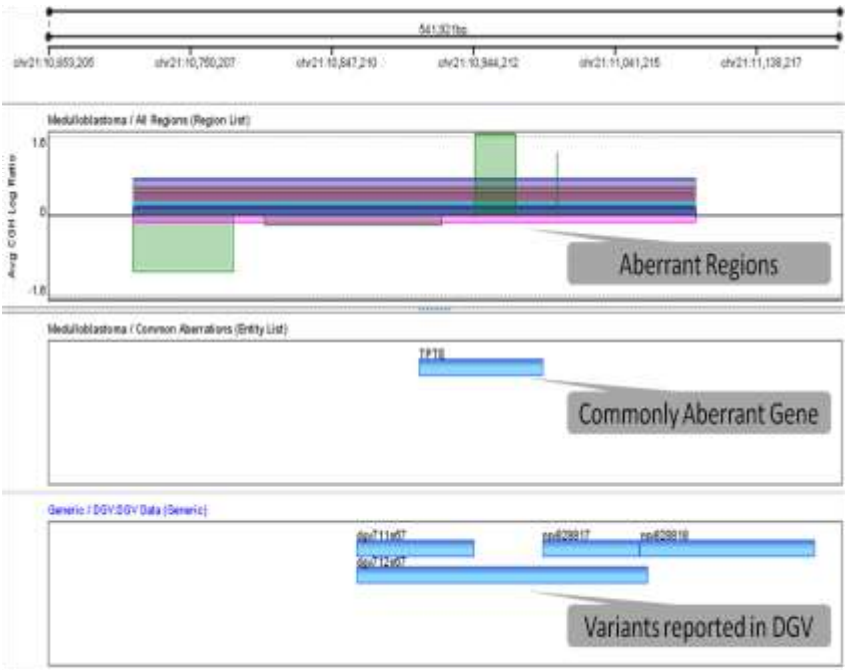
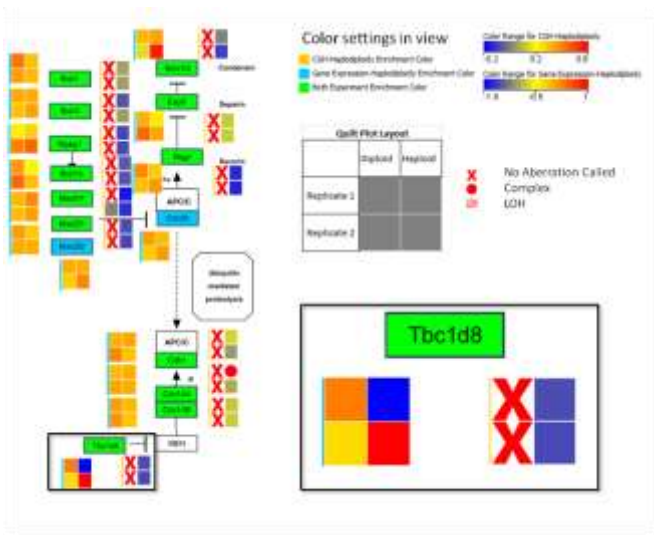
- [Proneural]
- [Classical]
- [Mesenchymal]

DOWNSTREAM PROCESSING AND BIOLOGICAL CONTEXTUALIZATION

CGH Analysis Workflow

- Reports from Cytogenomics and AGW are imported in GeneSpring GX for further downstream analysis and data integration
- Overlay on pathways will help cancer researchers visualize aberrations with other enriched entities

- Genome browser enhancements to support CGH data



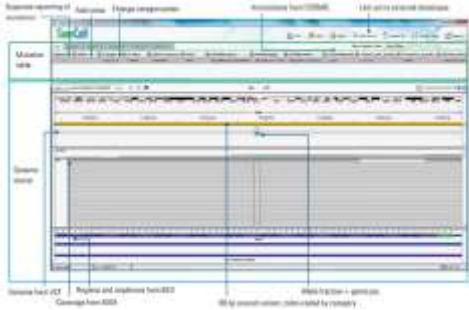
Variant Analysis with VCF Import

process

discover

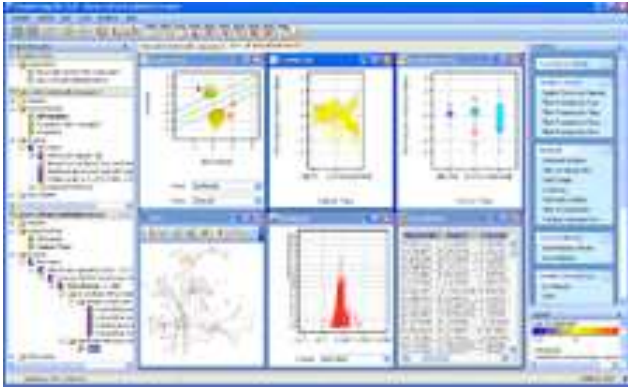


Strand NGS



SureCall

Any other NGS analysis tool



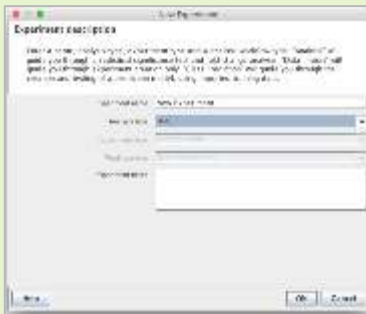
GeneSpring GX

Variant Analysis in GeneSpring GX

Data Import and Filtering

Visualization
and Results Interpretation

Pathway Analysis



- Import files in VCF format
- Annotate region list
- Region list operations

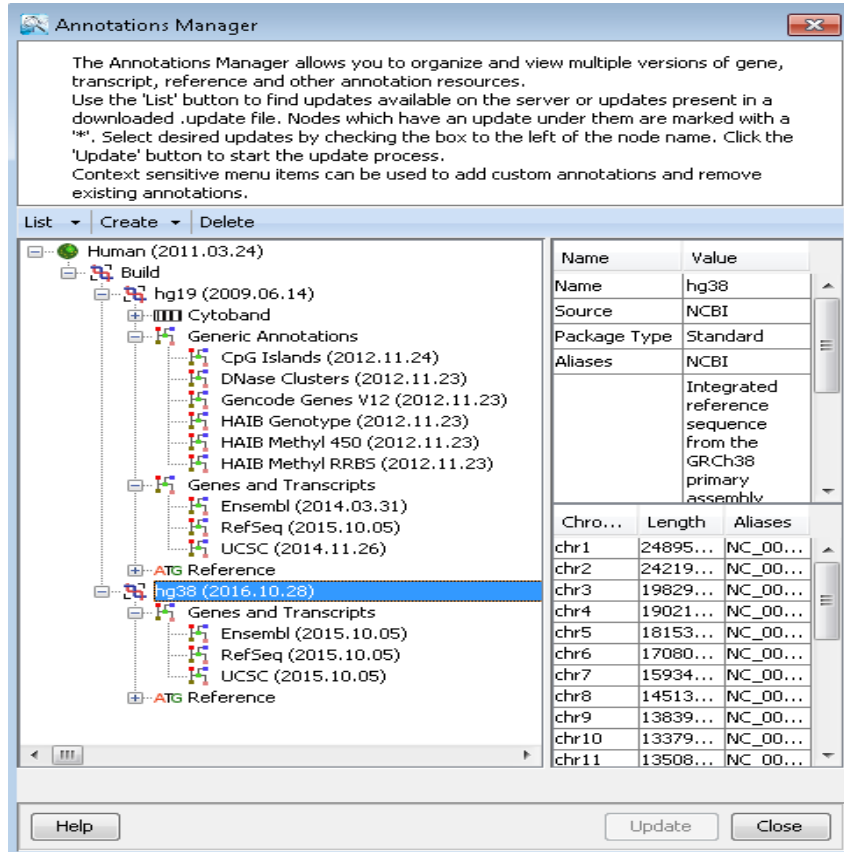


- Genome Browser
- Find Genic Parts
- Gene Ontology analysis



- Single experiment analysis
- Multi-omic analysis

Annotations Manager – Use available annotations or import from external sources



- Import annotations from ClinVar, COSMIC, dbNSFP, 1000 Genomes and more
- Annotate region list using annotations from Annotations Manager
- View annotations in Genome Browser

Filter genetic variants

The screenshot shows the 'Edit Region List' window with a table of genetic variants and a filter panel on the right. The table has columns for Chromosome, Start, End, Reference, Variant A, Variant T, VCF ID, Qual, and Fi. The filter panel on the right includes sections for Chromosome, Variant Allele, Variant Type, Alleles, Mapping Quality, and BPF.

Chromos...	Start	End	Reference	Variant A...	Variant T...	VCF ID	Qual	Fi
chr1	20915589	20915590	TC	T	Deletion		255	0/Pass
chr1	98349390	98349390	C	CA	Insertion		255	0/Pass
chr1	162750...	162750...	CTT	C	Deletion		255	0/Pass
chr1	165378...	165378...	TAC	T	Deletion		255	0/Pass
chr1	205640...	205640...	ACT	A	Deletion		255	0/Pass
chr2	25497684	25497684	A	AC	Insertion		255	0/Pass
chr2	29416025	29416025	C	CATTG	Insertion		255	0/Pass
chr2	29430942	29430944	GTC	G	Deletion		255	0/Pass
chr2	39224615	39224615	G	GT	Insertion		255	0/Pass
chr2	39224747	39224747	A	ATT	Insertion		255	0/Pass
chr2	39251375	39251375	T	TCAC	Insertion		255	0/Pass
chr2	209101...	209101...	GA	G	Deletion		255	0/Pass
chr2	212285...	212285...	AAAAG	A	Deletion		255	0/Pass
chr2	212544...	212544...	CT	C	Deletion		255	0/Pass
chr2	234668...	234668...	C	CAT	Insertion		255	0/Pass
chr3	37089840	37089841	CT	C	Deletion	rs63751	255	0/Pass
chr4	25678531	25678531	T	TG	Insertion		255	0/Pass
chr4	55131001	55131003	TAA	T	Deletion		255	0/Pass
chr4	55151958	55151958	T	TA	Insertion		255	0/Pass
chr4	55962545	55962545	T	TG	Insertion		255	0/Pass
chr4	55976947	55976947	T	TA	Insertion		255	0/Pass
chr4	143003...	143003...	G	GA	Insertion		255	0/Pass
chr4	162380...	162380...	T	TA	Insertion		255	0/Pass
chr4	162841...	162841...	T	TA	Insertion		255	0/Pass
chr5	56177848	56177851	TCAA	T	Deletion		255	0/Pass
chr5	67575323	67575325	ATT	A	Deletion		255	0/Pass
chr5	149441...	149441...	AC	A	Deletion		255	0/Pass
chr5	180038...	180038...	T	TC	Insertion		255	0/Pass
chr5	180046...	180046...	CAG	C	Deletion		255	0/Pass
chr5	180049...	180049...	A	AC	Insertion		255	0/Pass
chr6	117681...	117681...	G	GAC	Insertion		255	0/Pass
chr6	152265...	152265...	A	AT	Insertion		255	0/Pass
chr6	160645...	160645...	C	CT	Insertion		255	0/Pass
chr6	160666...	160666...	A	AC	Insertion		255	0/Pass
chr7	87257900	87257902	CGT	C	Deletion		255	0/Pass

- Create a new region list that only contains selected columns
- Filter the input region list based on specified conditions on the retained columns
- Add columns that are functions of selected existing columns

Find Genic Parts

Find Genic Parts

Select transcript model: **RefSeq Transcripts (2014.11.26)**

Select desired region types

- Exons
- Introns
- Upstream
- Downstream

Entity List: **All Entities** Choose...

Padding for Upstream: **5000**

Padding for Downstream: **5000**

Help OK Cancel



Region List Inspector

Name: **Genic Regions of RefSeq Transcripts (2014.11.26)**

Organism: **Homo sapiens**

Notes: **Source Entity List: All Entities
transcript model = RefSeq Transcripts (2014.11.26)
selected region types = [Exonic]**

Creation date: **Mon Feb 13 14:45:41 PST 2017**

Last modified date: **Mon Feb 13 14:45:41 PST 2017**

Owner: **gxuser**

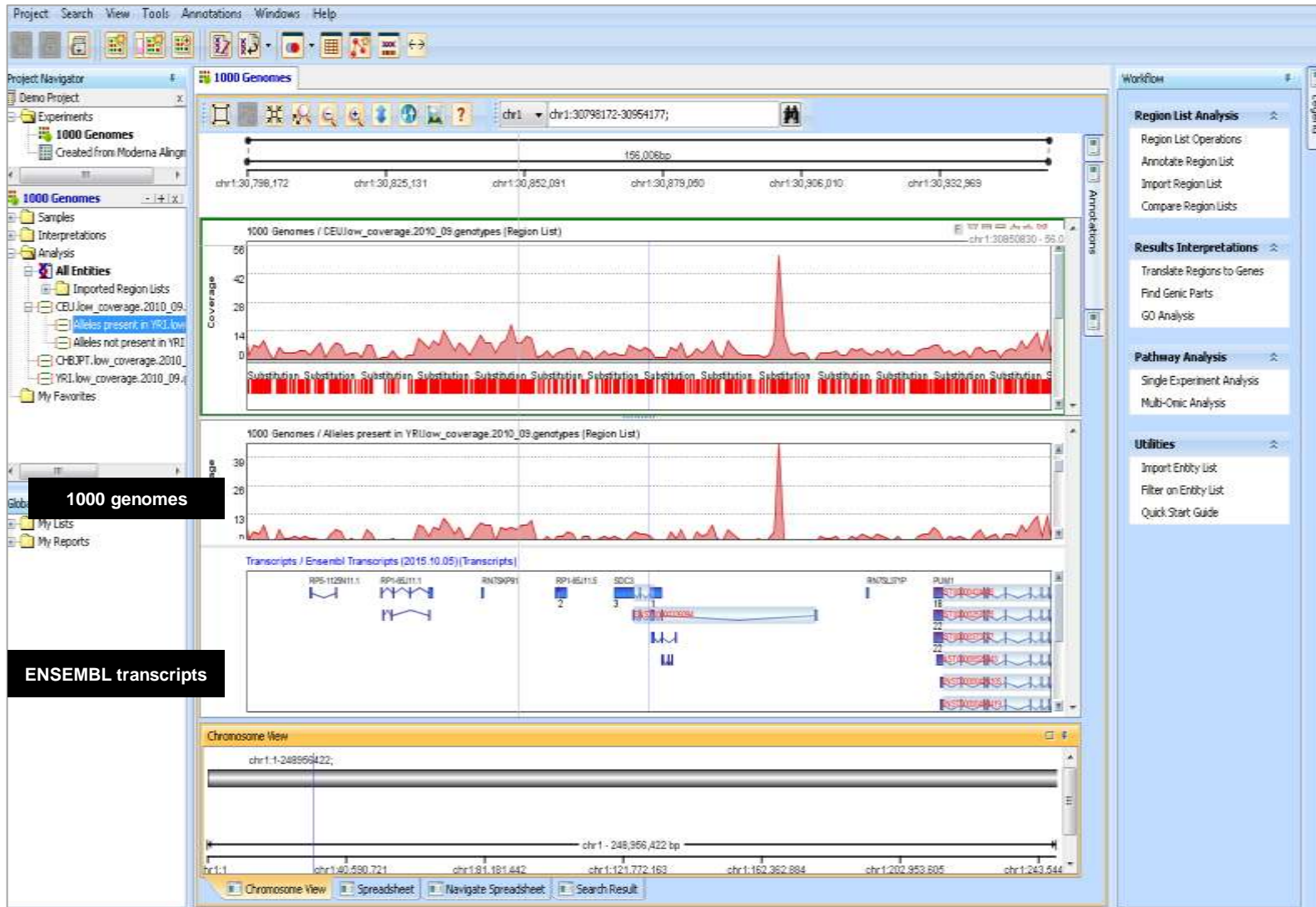
Number of Regions: **222,116**

Regions | Histogram | Summary Statistics

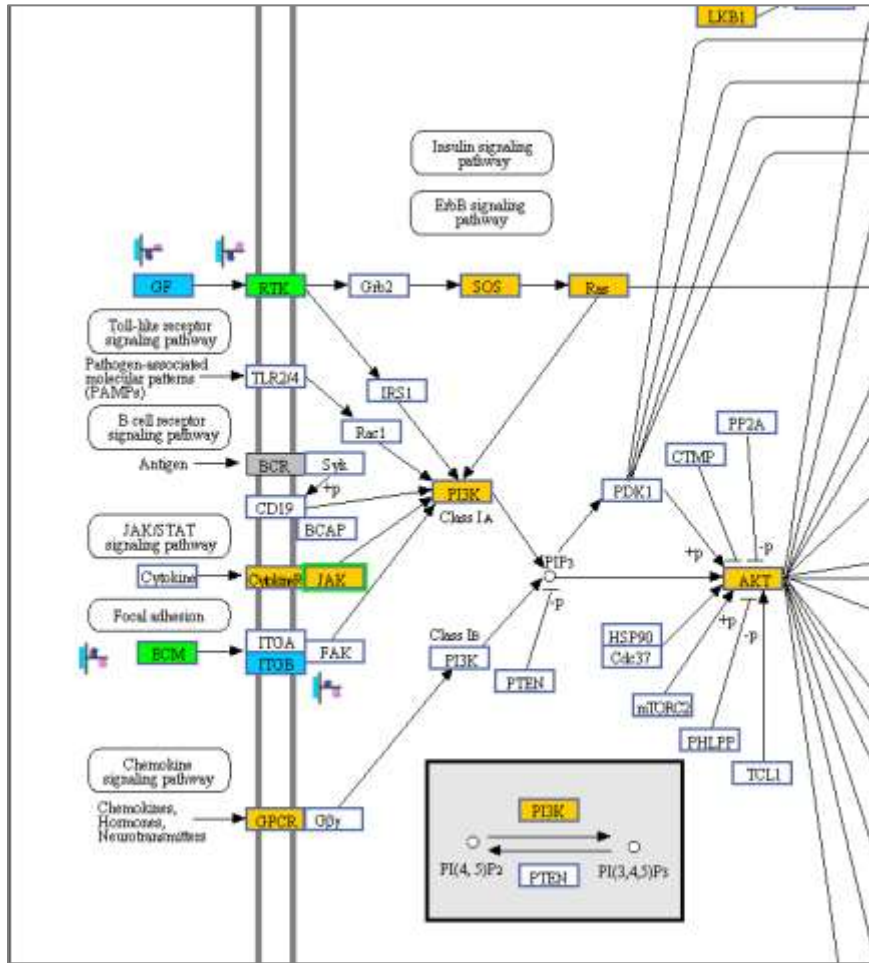
Chromos...	Start	End	Strand	Region T...	Gene ID	Gene Na...	Regi...
chr1	11874	12227	+	Exonic	100287...	DDX11L1	
chr1	12613	12721	+	Exonic	100287...	DDX11L1	
chr1	13221	14409	+	Exonic	100287...	DDX11L1	
chr1	14362	14829	-	Exonic	653635	WASH7P	
chr1	14970	15038	-	Exonic	653635	WASH7P	
chr1	15796	15947	-	Exonic	653635	WASH7P	
chr1	16607	16765	-	Exonic	653635	WASH7P	
chr1	16858	17055	-	Exonic	653635	WASH7P	
chr1	17233	17368	-	Exonic	653635	WASH7P	
chr1	17606	17742	-	Exonic	653635	WASH7P	
chr1	17915	18061	-	Exonic	653635	WASH7P	
chr1	18268	18366	-	Exonic	653635	WASH7P	
chr1	24738	24891	-	Exonic	653635	WASH7P	
chr1	29321	29370	-	Exonic	653635	WASH7P	

Help OK Cancel

Data Visualization



Multi-Omic Pathway Analysis



- Multi-omic experiment for data integration
- Overlay gene level information for variants
- Overlay other genomics, transcriptomics, proteomics or metabolomics data

