Phenotypic Characterization and Quality Control of Cells from Bacteria to Human Cells

Barry R. Bochner, Ph.D. CEO & CSO, Biolog, Inc. October 29, 2015





About ATCC

- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with and supports the scientific community with industry-standard biological products and innovative solutions
- Strong team of 400+ employees; over onethird with advanced degrees





Certification and Accreditation

ISO 9001:2008 Certification for quality management system

 Demonstrates commitment to quality products, customer service, and continued improvement



ISO 13485:2003 Certification for the design, development, production, testing, and distribution of medical devices

 Applies to synthetic molecular standards, the HIV surveillance kit, and other diagnostic and research kits

ISO Guide 34:2009 accreditation for production

 Applies to Certified Reference Materials (CRMs)



ISO/IEC 17025:2005 accreditation for testing

 Applies to all ATCC cultures, derivatives, and bioproducts tested in our laboratories





FM 610678



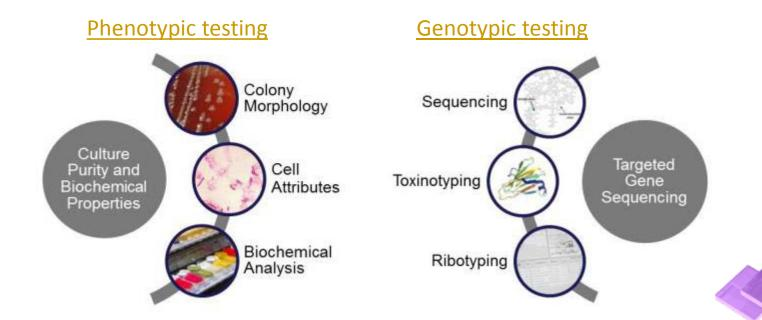
Highly Characterized Microbial Reference Strains



ATCC utilizes both classical and modern techniques

- Phenotypic analysis
- Genotypic analysis
- Proteotypic analysis
- Functional analysis

No single method of identification is sufficient



Phenotypic Characterization and QC of Cells

From Bacteria to Human Cells

Barry Bochner, Biolog, Inc., <u>bbochner@biolog.com</u>

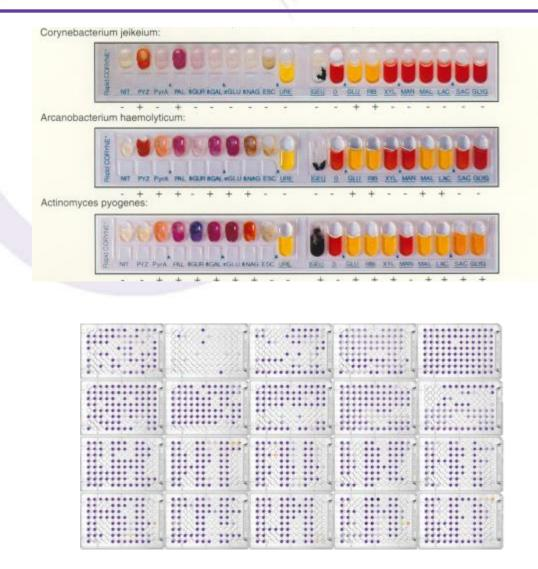
Biolog

Part One:

Phenotypic Characterization of Cells



Tremendous Advances in Phenotyping Technology



1970s Technology 20 tests

21st Century Technology

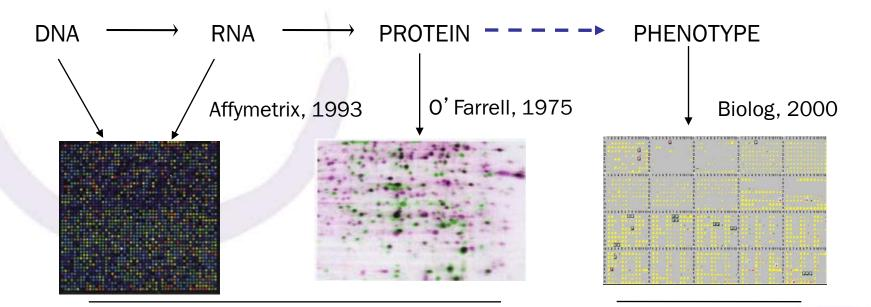
2,000 tests

Biolog

Tremendous Advances in Phenotyping Technology

- Hundred fold expansion in tests
- Much broader range of tests
- Single dye, single color chemistry
- Bacteria
- Yeast
- Filamentous fungi
- Algae
- Human and other Animal Cells
- Kinetic phenotypes

Phenomics = High Content Cell Phenotyping



Molecular Analyses

Transcriptomics

Proteomics

Phenomics

Biolog

Cellular Analysis

Complex Metabolic Circuitry of Cells

Biochemical Pathways	
	BIOLO

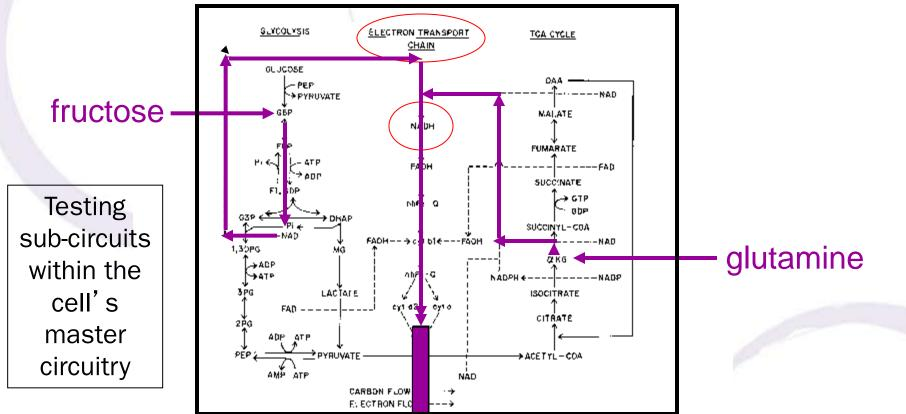
Biolog

Assay Principal:

Colorimetric Analysis of Energy Production



Metabolism of C-sources Produces an Electron Flow

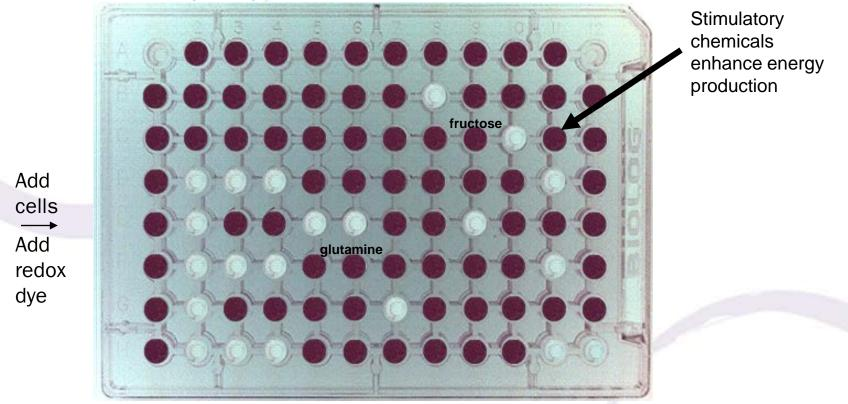


Fuc. 3. The pathways of central metabolism in are used: glucose-6-P (G6P), fructose-6-P (F6P), f diP-glycerate (1,3DPG), 3-P-glycerate (3PG), 2-P-f P (DHAP), methyl glycard (MG), non-heme iron late (OAA), and a-heloglutarate (aKG). coli and S. typhimurium. The following abbreviations tose-1.6-diP (FI 6DP), glyceraldehyde-3-P (G3P), 1,3craw (2PG), P-enolpyravate (PEP), dihydraxyacetoneyme Q complex (uhFe-Q), cytochrome (cyt), oxuloace-



Redox Chemistry Measures Cell Energetics

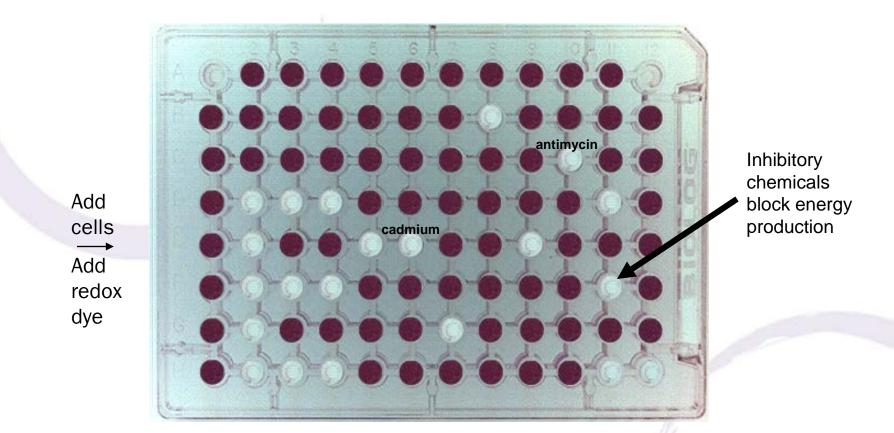
Microplate containing a negative control well and 95 different carbon substrates



Biolog

Wells contain different tests and measure different pathway activities and phenotypes of cells

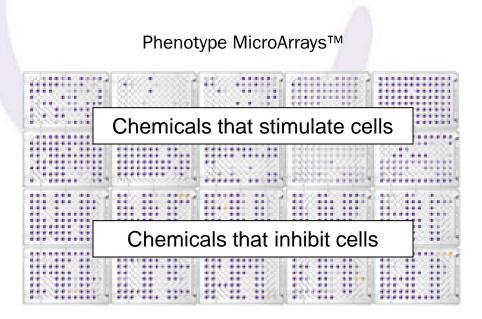
Redox Chemistry Measures Cell Energetics



Wells contain different tests and measure different pathway activities and phenotypes of cells

BIOLOG

2 Components of the Phenotyping Assay Platform



colorimetric cell assays in 96-well microplates

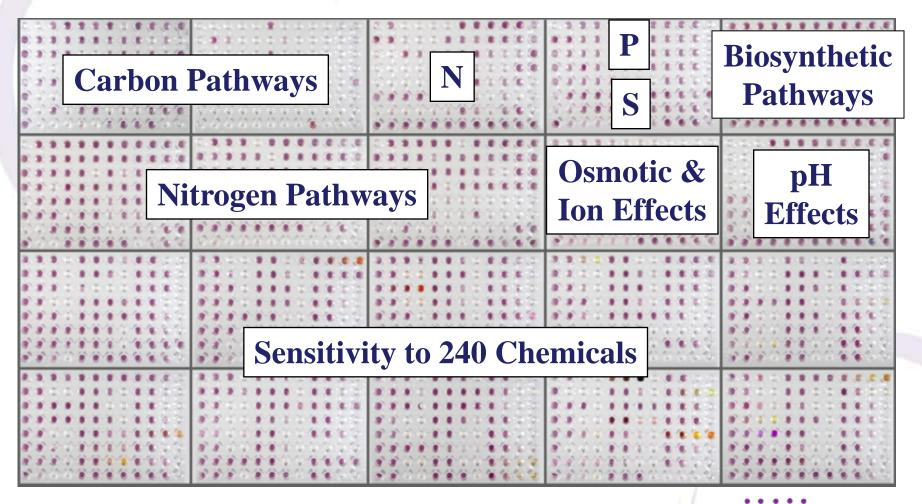
OmniLog[™] Incubator/Reader



incubation and recording of data in the OmniLog

Biolog

PM Platform - ~2,000 Phenotypic Assays



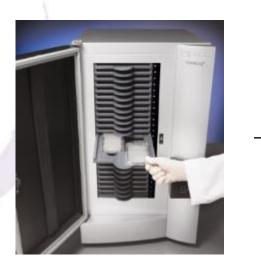
BIOLOG

PM Assays are Easy to Run



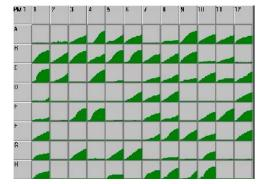
Assays Initiated by adding cells to wells

100 μ l per well





Holds 50 microplates at a set temperature and measures color formation at 15-minute intervals

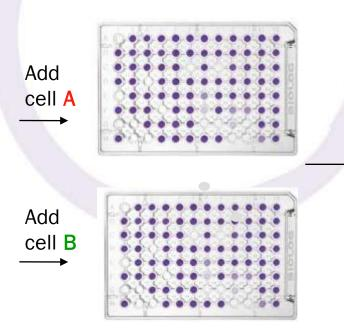


Kinetic assay readout for up to 5,000 wells

CVs typically < 10%

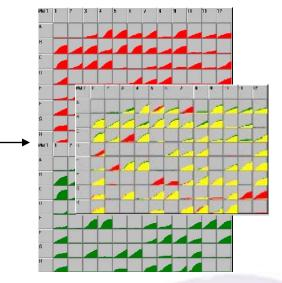
Biolog

PM Platform - Comparing Two Cell Lines



PM Pattern





OmniLog PM System

PM Kinetic Result

BIOLOG

PM Platform – Comparing Two Cell Lines

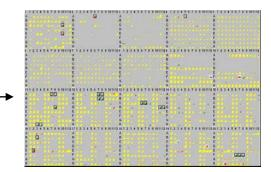
	: ÷. [300000000000		
****** **				
	· · · ·	004000440041	diliniiii f	
i valitadi		milia.	Minik.	
		·····	Silimily.	

PM Pattern

1 hr



OmniLog PM System



PM Kinetic Result

Automatic

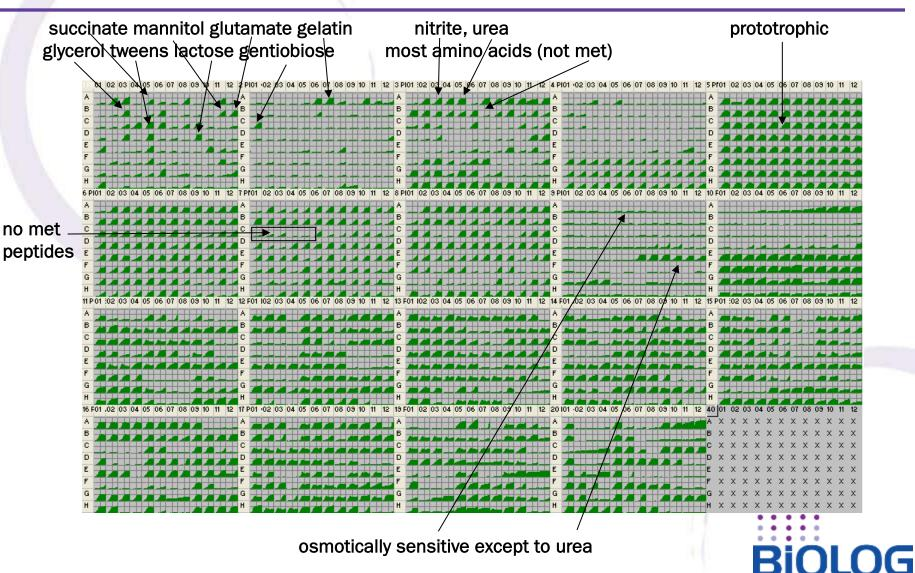
24-48 hr

BiOLOG

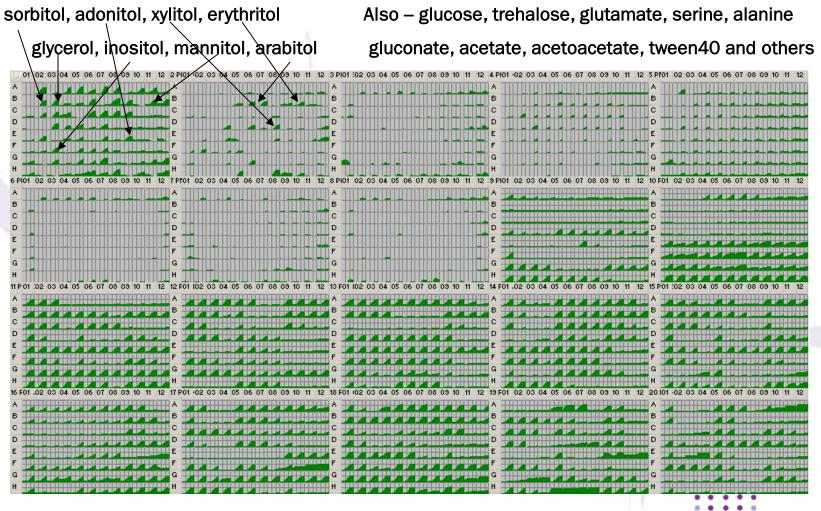
Examples of Phenomic Applications

- Profiling metabolism and chemical sensitivities of a cell
- Comparing properties of pathogenic vs non-pathogenic strains
- Determining the function(s) of a gene
- Analyzing environmental effects on cell phenotypes
- Optimizing production of a cell product e.g. a toxin

PM Analysis of Streptomyces coelicolor



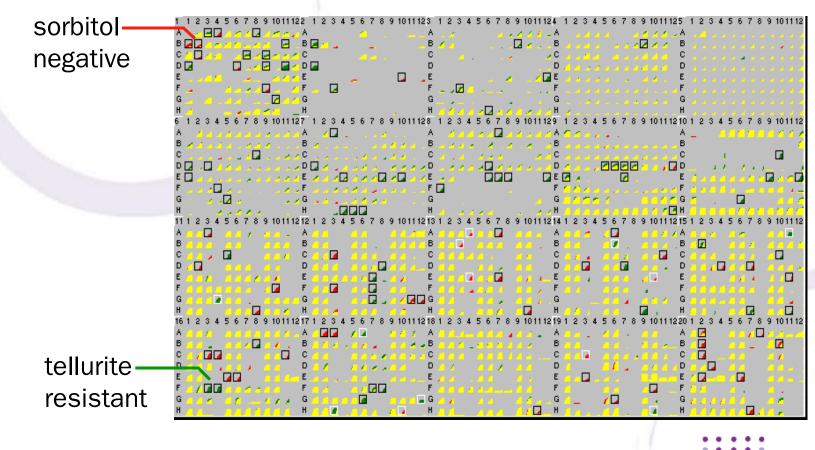
PM Analysis of Mycobacterium smegmatis MC²-155



BIOLOG

Comparing Two E. coli Strains:

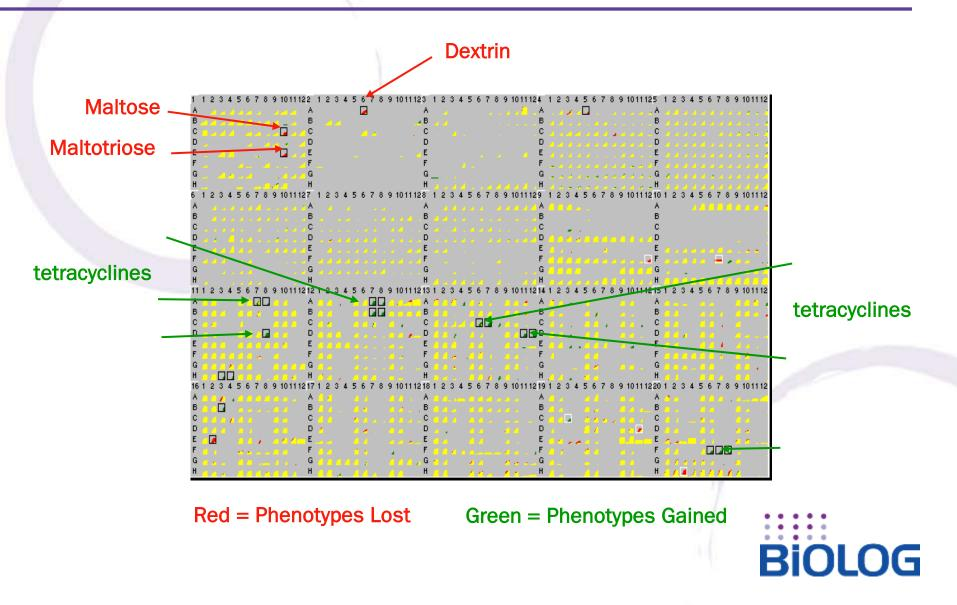
Pathogenic (0157) vs non-Pathogenic (MG1655) E. coli



BIOLOG

Comparing Wild Type vs Mutant:

E. coli malF::Tn10 vs MG1655



Comparing Psd.aeruginosa 8512 at 26 vs 36 C.

1 Pr01 02 03 04 05 06 07 08 09 10 11 12 2 Pl01 02 03 04 05 06 07 08 09 10 11 12 3 Pl01 02 03 04 05 06 07 08 09 10 11 12 4 Pl01 02 03 04 05 06 07 08 09 10 11 12 5 Pl01 02 03 04 05 06 07 08 09 10 11 12
6 PIO1 02 03 04 05 06 07 08 09 10 11 12 7 PIO1 02 03 04 05 06 07 08 09 10 11 12 8 PIO1 02 03 04 05 06 07 08 09 10 11 12 9 PIO1 02 03 04 05 06 07 08 09 10 11 12 10 FOI 02 03 04 05 06 07 08 09 10 11 12
· ANCHARANAACA MANAAAAAAAAA MAAAAAAAAAAAAAAAA TITA
" " " " " " " " " " " " " " " " " " " "

Biolog

Culture Conditions Inducing Toxin Synthesis

	Tme(dass)	The identian	6 61
	A STATE OF	State State	Карияна Совтт Анитика
	A COMPANY AND		Rithe
	The second		Ktosta Une 0
20 8	NUMBER		Bure: 91 8 Leterne 40 5
5 19		CONTRACTOR OF	L-Angloine 00 T
82 2	A DATE STREET,	No. 1	Lógetrán 80 % Lógetrán 10
100	0.00		L FRANKO AC 1
		The second second	L-Glutenine Gijelne
	1 TO BERRIES	CONTRACTOR OF	Lifeblacine Lisoleacine
		A CALIFORNIA	LLALONS LLANNE
	ALC: NOT	the state	E Misthonine E Promplake en
	T = PROUND	Thurs	L.Fraine L.Gette
	Contraction of the		Litheonine
	Sector Sector		L-Trystophan L-Trystophan
	1200		Dividine Dividine
			O Aspengine D-Aspeni o Asia
	100		D-Gildamic 2d 5
	13255	1988	Dilystro DiSette
	. TRANSPORT		D'Adrie L'Otrutine
	1.1.1	Me and a	Chonesente L'Orielline
			CARLENDER A LAND REPORT AND
	A MAGNINGS	a an an a	NP 10 May 4 - G som c Add L-Pyrogius nic Add L (vogius nic Add L (vogius nic Add
	100	建 田中人的时	Moliny land to .
		August -	MARSon in
		NUMBER OF A	E tylanice Lite werke
-	1 Contractor	1	Etylandanine Ozosoine
			Agnethe
		C A STATE	B-Phoryx Belonine
	STOLENSING.	人。 一個語	Kustanice Formar de
	A PARA AND		Chic are twitch
		100 A.	DL-Lacamète D-C activement D-Calentraumère
			O-Venicovinie
	BENETATION		Manager Construction Network Distances on the
	- Composition	A BOULDER	Ni Asiarya Tattaning ang ing Adamina
	1000000	Section Participation	Adenosine
	A STATE		Cyddinu Cyddina
	1000 COLOR	Contractory of the local division of the loc	Oueroveia
			Taymine Taymin ee
	1 Street and	公司 11日日	Lined Linder
	- ANNA		inceine Xenthee
	STATE IN STATE		Xandrosino. Una Add
		The local division of	Mean
			Alexilor Perahanic Acid
	机器理的	建 成 1 字	CL-ro-Amino-N-Curynic Add AAmino-N-Eurork Add
		16 60 5	eAniro-N-Capito Adi. E. Li u vinisto Gabrylo Adia
		Paris and	6-Amino-N-valorie Aced Exemino-re-salance Aced
	and the second	E-EASTREE	No-Net-
	CAL ASS	S. S. BELS	Ha-Gis Ala-Gis
	A DEC	Contraction of the second	Alter Try Alter His
		all a	Neri-m. Neri-m.
	- La Castalia	WHER !!	Gradian Gradia
	ALL SUPPLY		G pilder
	the second second	and the second	Wet-Me

Culture conditions inducing synthesis of a trichothecene mycotoxin in the wheat pathogen, *Fusarium graminearum*. A special strain was constructed with a toxin gene promoter fused to GFP.

Induction was highest with arginine, putrescine, agmatine, and guanine as nitrogen sources.

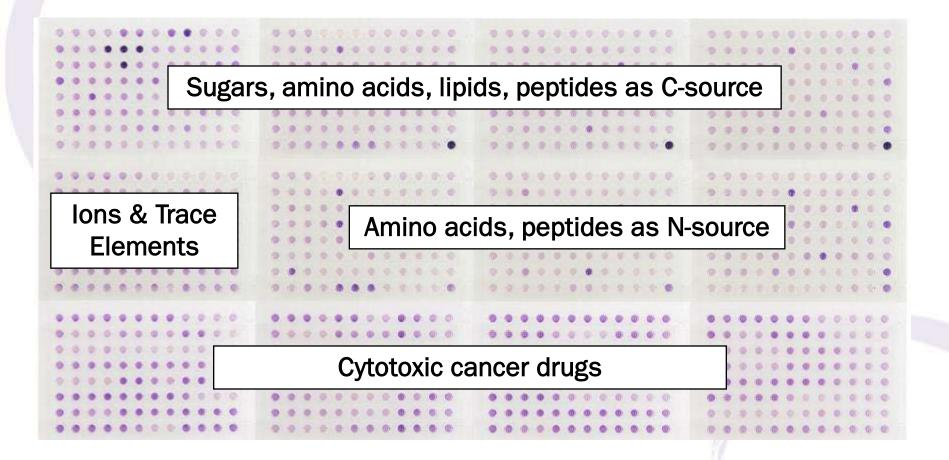
D. Gardiner et al – Fungal Gen & Biol (2009)

Phenotyping Human Cells:

Analyzing Human Cell Metabolism, Human Genes, Human Disorders, and Effect of Drugs



~1500 Assays and Culture Media for Mammalian Cells



Plus 3 panels with hormones, cytokines and other bioactives



Harry Eagle, JBC 1958

The Utilization of Carbohydrates by Human Cell Cultures

HARRY EASTR, STANLEY BANBAN, MINA LEVY, AND HENRY O. SCHOLZE

From the Section on Experimental Therapeutics, Laboratory of Infectious Discuss, Natural Institute of Allergy and Infectious Discusses, National Institutes of Health, Public Health Service, United States Department of Health, Education, and Welfere, Bethesda, Maryland

(Received for publication, April 21, 1958)

TABLE 1

Ability of a variety of carbohydrotes" and related arcpounds in substitute for in-planars in cell relaters The compounds were tested at varying levels (0.5 to 50 mW) in

a ghreine i an geiser wee "remplate" g<u>owih modum upsiemwaled with</u> dialyzed *serum*. The tetler in parentizers after eweb compound indicate the cell ince against which it was tested i

Company/a which did wat rapped grants	Concourds work[0] op abgin gravit, but ins abas with glasses	Congregade prim fride growth all rais conservable with that on glutour
Lasture (G, H. L, T)	Colleborn (I: G,	Treasure (C. G. D.
Sucrous (H. L. T.	$\Pi, \mathbf{K}, \mathbf{M}, \mathbf{T})$	H. M. TI
\mathbf{p} -Allow (G, T)	Melihinan (C, G, H, K, M, T)	Такеаана (С. С. Ц. Х. М. Т)
p. Algrage (G, T)	Sorbital (all)	7, 01. 1)
p-Galaritow (L)‡	·····	p.Fructose (all)
D-Galacturonic actd	D-Thion (G, T)	p-Galactore (G. G.
(H, T)	o-Xylner (all)	13, E, M, T)‡
to-Glumnic antid (h.,	o-Ribon (G, H,	s-Glucoss (all)
T;	1. II	c Glucase t PO, OI.
Methyl-a-n ghrenside		K. 5D
Mathyi dia glaziaide		E-Glucos-6-PO, (3),
a Closerosas acid		K.M)
(K, T) - Culue (K, T)		is-Mangues (all)
z-Gology (K, T) z-Almuzital (Alli		
L-Rhamness (N, T)		
1-Scrbese (all)		
L-Asabibuse (G)		
p-Arabinose (G)	6	8
p-Lyncer (C)		
p Ribnes (C. L. M.		
ገን <mark>ት</mark>		
Xylltel (G)		
p Brytheam (II, 1.)		
Arntabe (B. J. T)		
Citease (B, L, T)		
Pumarate (B, L, T)		
Givearal (E. L. T)		
Lastate (H, L)		
Malato (II, L. T)		
Ethosphoglycezate		
(H, K)		
Pyrovato (£1) Succlaste (£1, 3, T)		
automie (II, 4, 1)		ί.



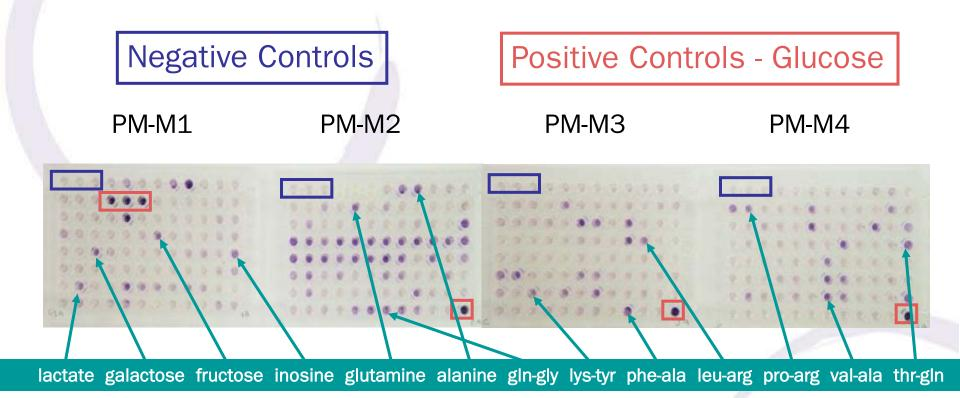
PM-M1 to M4: 367 Carbon-Energy Substrates for Cells

Negative Control	A2 Negative Control SUDSI	A3 Negative Control	A4 α-Cyclodextrin	A5 Dextrin	A6 Glycogen	A7 Maltitol	A8 Maltotriose	A9 D-Maltose	A10 D-Trehalose	A11 D-Cellobiose	A12 β-Gentiobiose
B1 D-Glucose-6- Phosphate	B2 α-D-Glucose-1- Phosphate	B3 L-Glucose	B4 α-D-Glucose	B5 α-D-Glucose	B6 α-D-Glucose	B7 3-O-Methyl-D- Glucose	B8 a-Methyl-D- Glucoside	B9 β-Methyl-D- Glucoside	B10 D-Salicin	B11 D-Sorbitol	B12 N-Acetyl-D- Glucosamine
C1 D-Glucosaminic Acid	C2 D-Glucuronic Acid	C3 Chon Sulfale)sacc		ës,	C11 Palatinose	C12 D-Turanose
)1)-Tagatose	D2 L-Sorbose	D3 L-Rhamnose	D4 L-Fucose	D-Fucose	D-Fructose-6- Phosphate	D-Fructose	Stachyose	D9 D-Raffinose	D10 D-Lactitol	D11 Lactulose	D12 α-D-Lactose
1 felibionic Acid	E2 D-Melibiose	E3 D-Galactose	E4 α-Methyl-D- Galactoside	E5 β-Methyl-D- Galactoside	E6 N-Acetyl- Neuraminic Acid	E7 Pectin	E8 Sedoheptulosan	E9 Thymidine		Ett Adenopine	E12 Inosine
1 .donitol	F2 L- Arabinose	F3 D-Arabinose	F4 β-Methyl-D- Xylopyranoside	F5 Xylitol	F6 Myo-Inositol	F7 Meso-Erythritol	F8 Propylene glycol	F9 Ethanolamine	F10 D,L-α-Glycerol- Phosphate	F11 Glycerol	F12 Citric Acid
i1 ricarballylic Icid	G2 D,L-Lactic Acid	G3 Methyl D-lactate	G4 Methyl pyrnyate	G8 Pyravic Acid	Gê α-Keto-Glutaric Ac <mark>a n C</mark>	G7 Succinamic Acid			G10 L-Malic Acid	G11 D-Malic Acid	G12 Meso-Tartaric Acid
H .cetoacetic Acid)	H2 y-Amino-N- Butyric Acid	H3 a-Keto-Buytric	H4 a-Hydroxy- Buty B cid	H5 D,L-β-Hydroxy- Hutvric Acid	H6 P-Hydroxy- Butyric \cid	H7 Butyric Acid	HB 2,3-Butanediol	H9 3-Hydroxy-2- But mone	H10 Propionic Acid	H11 Acetic Acid	H12 Hexanoic Acid

Each well has a different substrate

Biolog

PMs Assay Carbon/Energy Pathways in Cells



sugars, alcohols, acids

fatty acids, amino acids, and dipeptides



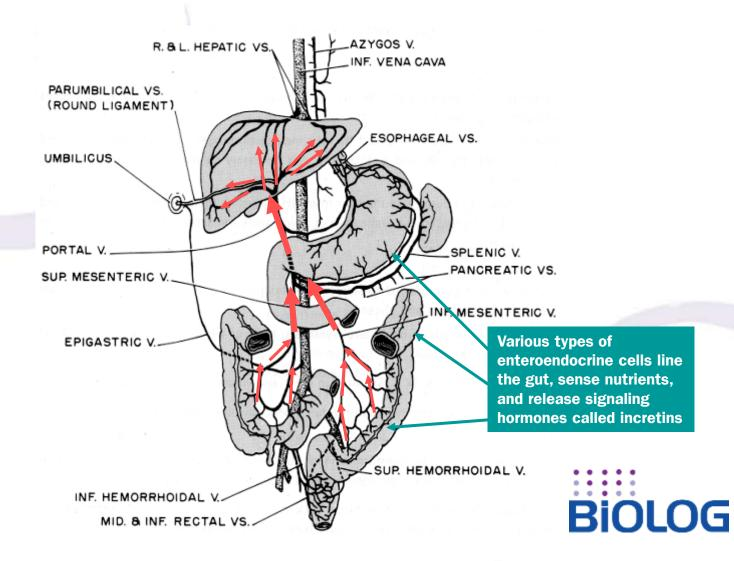
Each Cell has a Different Set of Energy Pathways

	PM-M1	PM-M2	PM-M3	PM-M4	
CCRF-CEM (lymphoid) maltotriose, maltose (A8-9) glucose (B4-6) mannose (C5)					
HL-60 (<mark>lymphoid</mark>)					
PC-3 (prostate) fructose (D7) uridine, adenosine, inosine (E10-12) pyruvate, succinamate, mono-methyl succinate (G5,7,9)					
A549 (lung) dextrin, glycogen (A5-6) darker wells in PM-M2, M3, and M4 correspond to glutamine and gln-peptides			•		
COLO 205 (colon) galactose (E3) lactate (G2) butyrate, propionate (H7,10)					
HepG2 (liver) darker wells in PM-M2, M3, and M4 correspond primarily to alanine and glutamine and ala, gln, and arg-peptides					
HepG2/C3A <mark>(liver</mark>)	**********		*********		
pyruvate (G5)			•••••••••••		

Adapted from Bochner et. al. PLoS ONE (2011) 6:e18147

BIOLOG

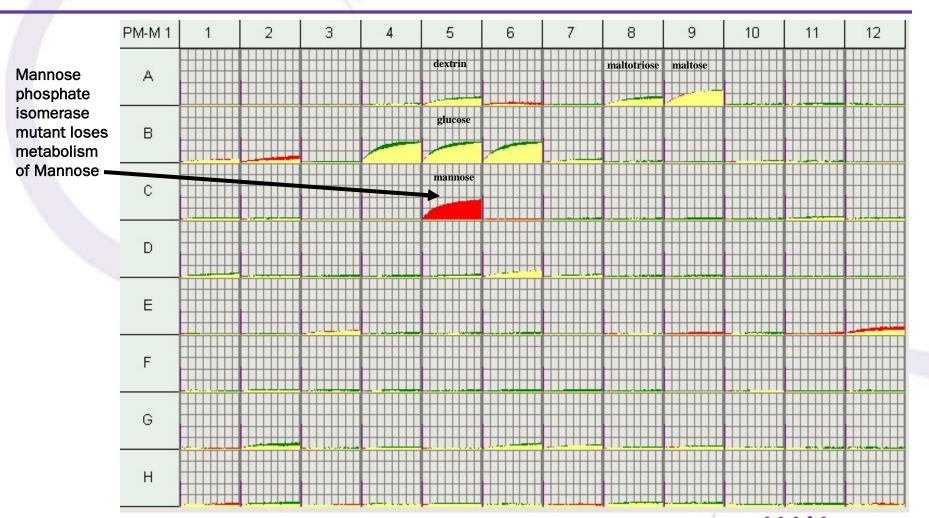
Physiology of Nutrient Absorption



Analyzing Mutations in Human Cells

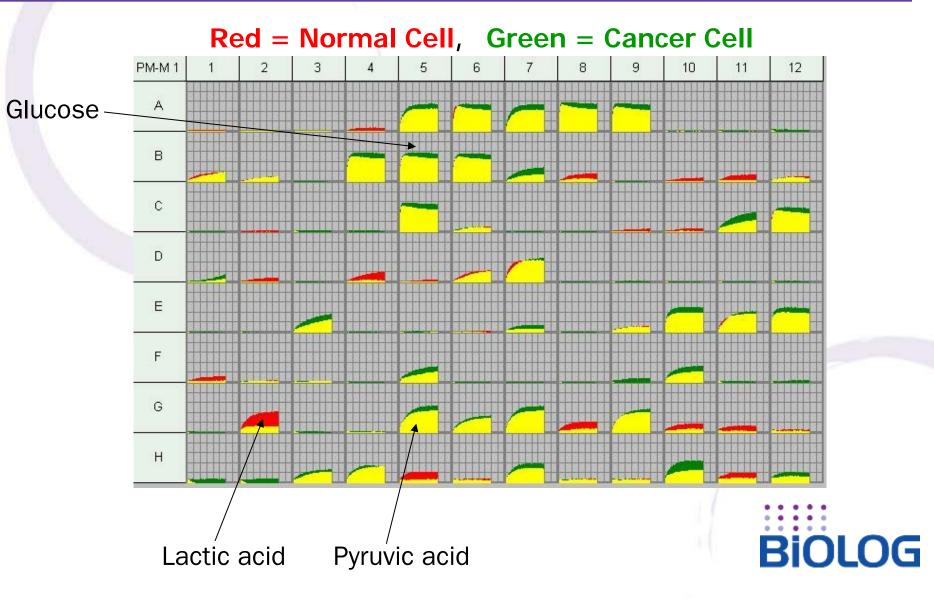


Comparison of HAP1 vs HAP1_MPI_124-16 on Biolog Redox Dye MB reduction (Average of 4) after dispensed into PM-M1 plates at 20,000/well in IF-M1 + 0.3 mM Gln + 5% FCS + PS medium and incubated for 20 h.

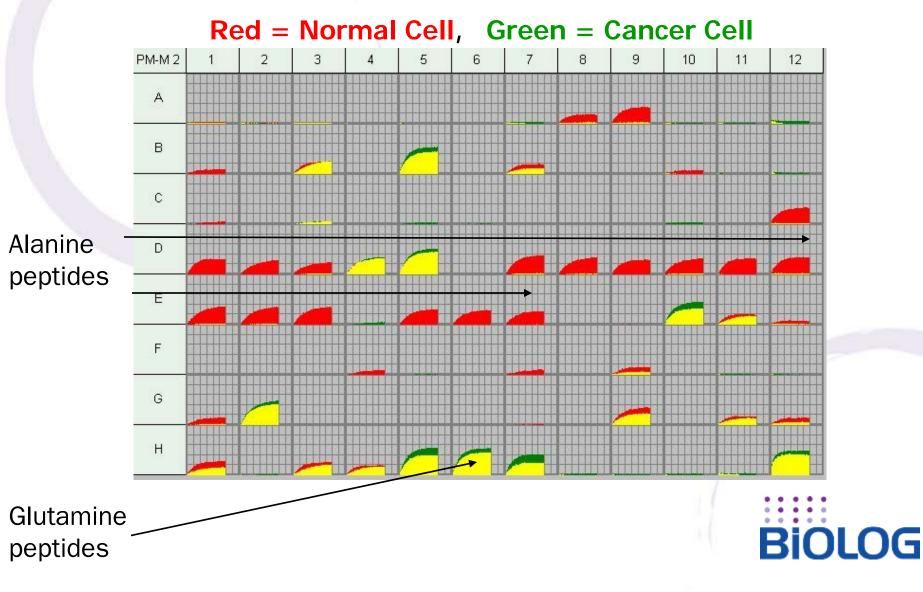


HAP1 and HAP1_MPI_MPI_124-16 (Mannose Phosphate Isomerase), both grown in IMDM + 10% FCS + PS, were dispensed into a PM-M1 plate at 20,000 cells/well in 50 uL IF-M1 + 0.3 mM Gln + **SOLOG** FCS + PS, incubated 20 h at 37°C under 5% CO₂-95% air before adding 10 uL/well Dye MB, plates sealed with tape and placed in a 37°C OmniLog. Dye reduction over ten hour minus background

Metabolic Changes in MCF10a vs PI3K Clone CL1



Metabolic Changes in MCF10a vs PI3K Clone CL1



Assay medium: IF-M1 +5% horse serum +0.3mM Gln +1xP/S+ 0.1 ug/mL cholera toxin +10 ug/mL insulin+0.5 ug/mL hydrocortisone +0.2 ng/m EGF

20141208-12_PM-M Tox2_PI3K Inhibitor Titration_68h in CO2_24h in OL with Dye MA

Cells: 20k/well

BYL 719 (nM)

MCF10a / PIK3CA (H1047R/+) MCF10a (WT) BYL 719 (nM) BYL 719 (nM) 0 0.5 1 2 3.9 7.8 16 31 63 125 250 500 0 0.5 1 2 3.9 7.8 16 31 63 125 250 500 0 0.5 1 2 3.9 7.8 16 31 63 125 250 500 0 0.5 1 2 3.9 7.8 16 31 63 125 250 500 0 0.5 1 2 3.9 7.8 16 31 63 125 250 500 None Glycogen Incatate <t

BIOLOG

Part Two:

Quality Control of Cells



Who Needs to Perform Cell Line QC?

- Culture collections and cell banks
- Bioprocess scientists banking cells (seed cultures) to inoculate fermentations
- Cell-based assay labs
- Anyone using cells in their research



How is Cell Line QC Currently Performed ?

For Microbial Cells

• A mixture of genetic and phenotypic tests

For Human and Animal Cells

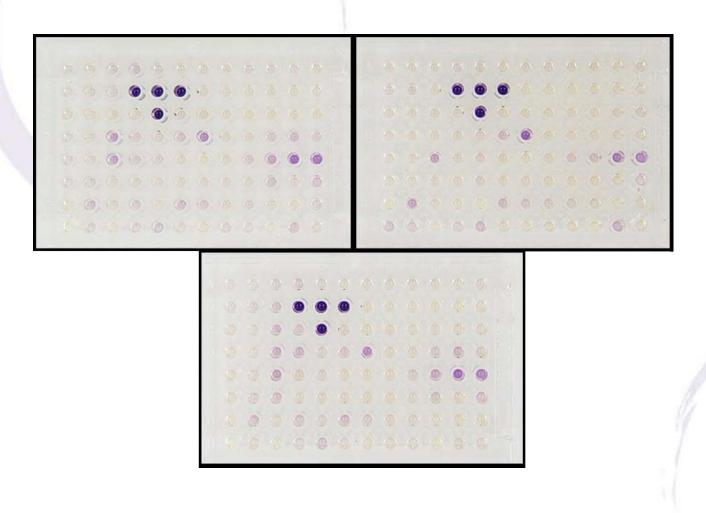
- STR profiling according to cell line authentication method ASN-0002 (genetic method)
- Many labs rely on judging cell morphology



Analyzing Metabolic and Other Phenotypes Provides a Complementary QC Method

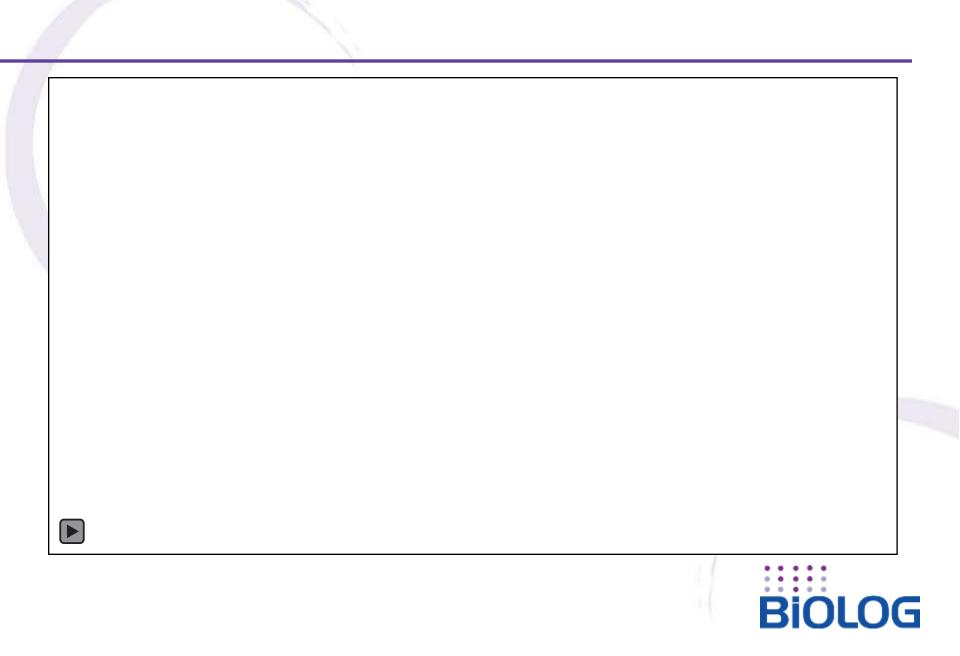
RiOl OG

PM Fingerprints are Highly Reproducible



CVs typically are less than 10%

Biolog



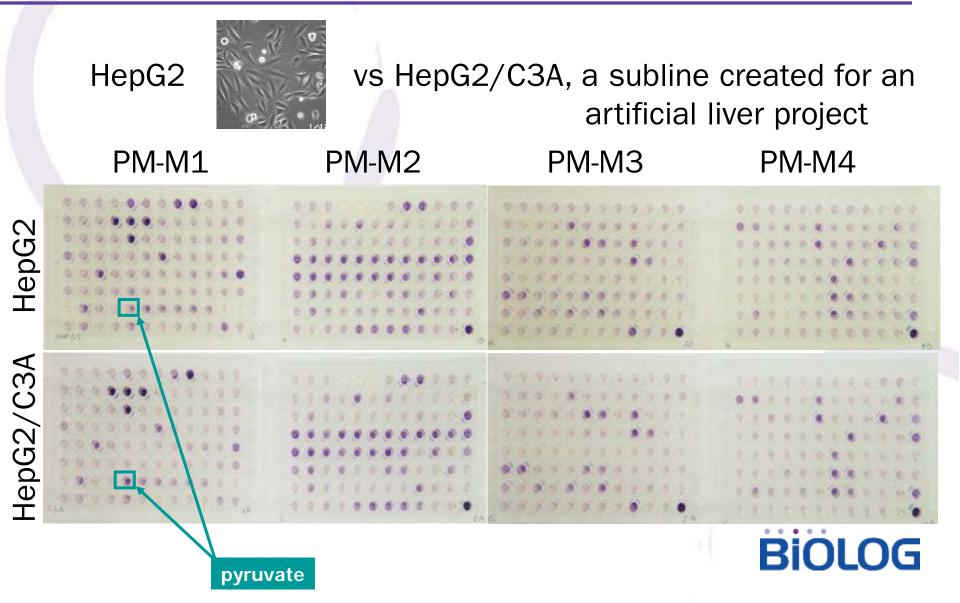
Cells Have Different Metabolic Phenotypes

	PM-M1	PM-M2	PM-M3	PM-M4	
CCRF-CEM (lymphoid) maltotriose, maltose (A8-9) glucose (B4-6) mannose (C5)					
HL-60 (<mark>lymphoid</mark>)		•			
PC-3 (prostate) fructose (D7) uridine, adenosine, inosine (E10-12) pyruvate, succinamate, mono-methyl succinate (G5,7,9)					
A549 (lung) dextrin, glycogen (A5-6) darker wells in PM-M2, M3, and M4 correspond to glutamine and gln-peptides	• • • • • • • • • • • • • • • • • • •	:	•		
COLO 205 (<mark>colon</mark>) galactose (E3) lactate (G2) butyrate, propionate (H7,10)					
HepG2 (liver) darker wells in PM-M2, M3, and M4 correspond primarily to alanine and glutamine and ala, gln, and arg-peptides					
HepG2/C3A <mark>(liver</mark>)	***		**********		
pyruvate (G5)			5 5 - 5 - 5 - 5 - 5 - 5 - 5 - 5 - 5 - 5		

Adapted from Bochner et. al. PLoS ONE (2011) 6:e18147

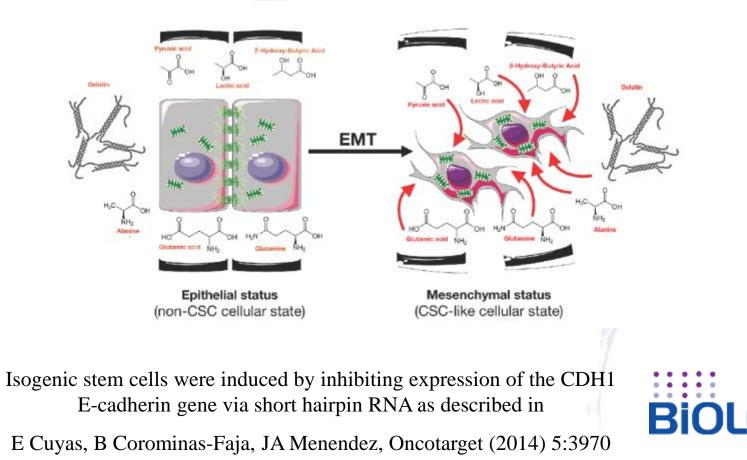
BIOLOG

PMs Distinguish Closely Related Sublines



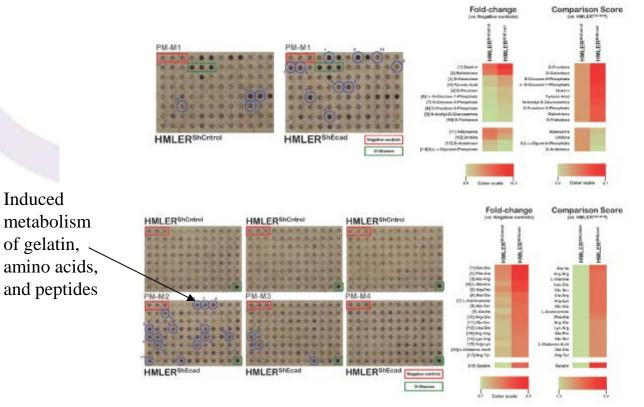
PM Comparison of Breast Cancer vs Cancer Stem Cell

Using the method of Robert Weinberg, cells can be induced to undergo an epithelial to mesenchymal transition (EMT) with stem cell-like properties



PM Comparison of Breast Cancer vs Cancer Stem Cell

"Most changes that occurred following the acquisition of a CS-like cellular state (28 out of 31, 90%) were increases in the ability to generate energy"

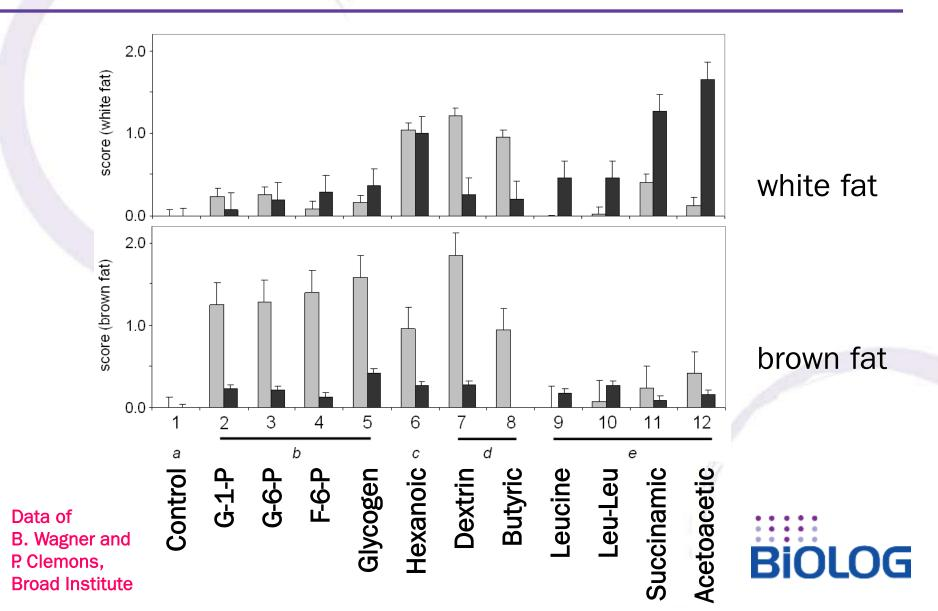


Isogenic stem cells were induced by inhibiting expression of the CDH1 E-cadherin gene via short hairpin RNA as described in

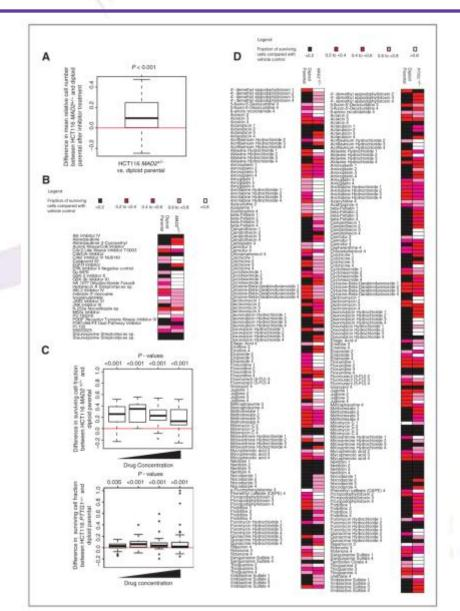
E Cuyas, B Corominas-Faja, JA Menendez, Oncotarget (2014) 5:3970

BiOl

Metabolic Differences in Preadipocytes and Adipocytes



Resistance to Anti-Cancer Drugs in CIN⁺ Cancers



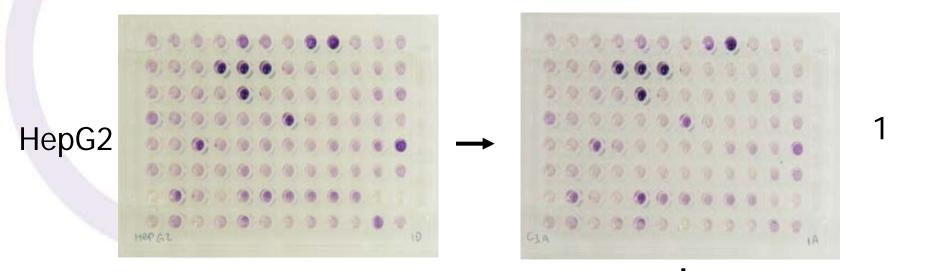
From Lee et. al. Cancer Res.(2011) 71:1858

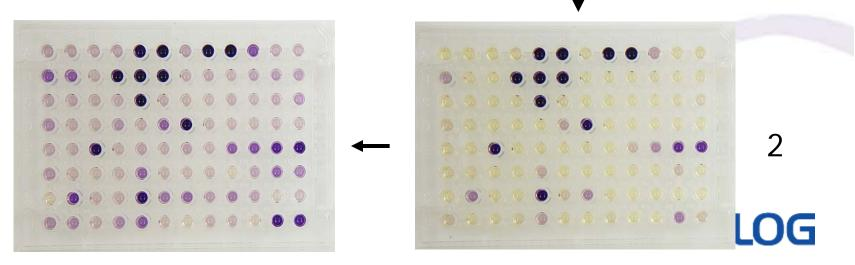


Examples of Cell Line Instability Detected by Metabolic Phenotype Analysis



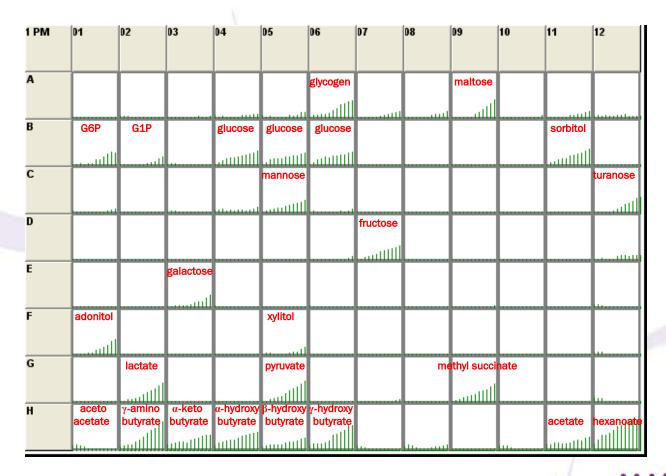
Modification of a Hepatocyte Cell Line





3

Detailed Metabolic Analysis of Primary Hepatocytes



Primary rat hepatocytes, 20,000 per well, incubated for 3 hours

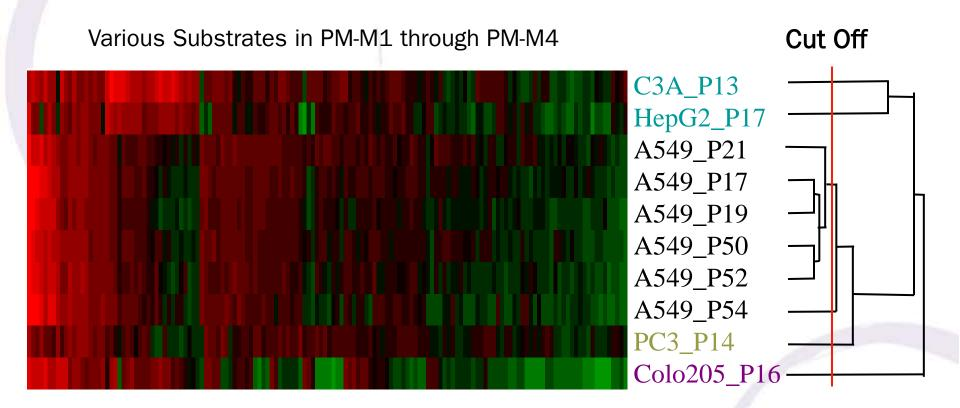


Hepatocyte Preps Show Lot-to-Lot Metabolic Differences

Lot 208 Lot 212 Lot 256 A в С D G

BIOLOG

PMs Detect Changes with Cell Passaging



- ·QA criteria for cell line stability can be experimentally determined
- Cross-contamination can be easily detected

Issues and Limitations of DNA-based QC

- Wide ranging cell stability from unstable stem cells to relatively stable cell lines that have been maintained in artificial culture for decades.
- DNA-based QC would not be a preferred method for characterizing and distinguishing
- cells of different tissue/organ type
- cells with point mutations
- ips reprogrammed cells, stem cells, differentiated cells
- cells with chromosome instability
- cells with epigenetic changes
- cells with metabolic changes

All of these can potentially be detected by analysis of metabolic and other phenotypes



Other Reasons for Performing Metabolic QC

Cell Metabolism based Bioprocess Production

- CHO cells producing recombinant proteins
- Hybridoma cells producing monoclonal antibodies
- Yeast producing special wines
- Bacteria producing antibiotics
- Bacterial cocktails producing special yogurts or probiotics for faecal transplants.

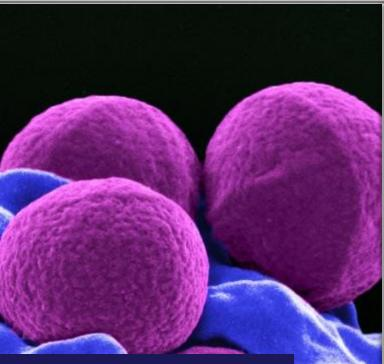
Cell-based Assays

• Primary liver cells used in tox assays

Summary

- There is no perfect solution or right answer to how much and what type of cell line QC should be performed. It is a judgement call based on assessment of how stable the cell is and how best to detect changes that it is prone to make.
- In general, it is prudent to perform a metabolic/phenotypic QC in addition to a genetic QC as these are complementary analyses.
- Biolog phenotypic assays provide an ideal metabolic/phenotypic platform that spans the spectrum of cells from bacteria to human cells.
- Biolog assays are very easy and inexpensive to perform and they can be read with any microplate reader or even by eye.
 However the best data is obtained by measuring metabolic rates using the OmniLog instrument.

The Importance of Authentication



Methicillin-resistant *Staphylococcus aureus* -Photo courtesy of NIAID

- The use of minimally cultured strains that have been propagated and preserved under the proper conditions can help prevent:
 - Genetic drift
 - Phenotypic variation
 - Changes in functional characteristics
 - Contamination
 - Misidentification



Thank you for joining today!

Register for more ATCC *"Excellence in Research"* webinars, or watch recorded webinars, at <u>www.atcc.org/webinars</u>.

November 12, 2015
10:00 AM, 3:00 PM EST
Bill Hirt, Ph.D., *Director of Accreditation*, ANAB
How Does ISO 17025 Accreditation Build
International Confidence?



Please email additional questions to: tech@atcc.org

