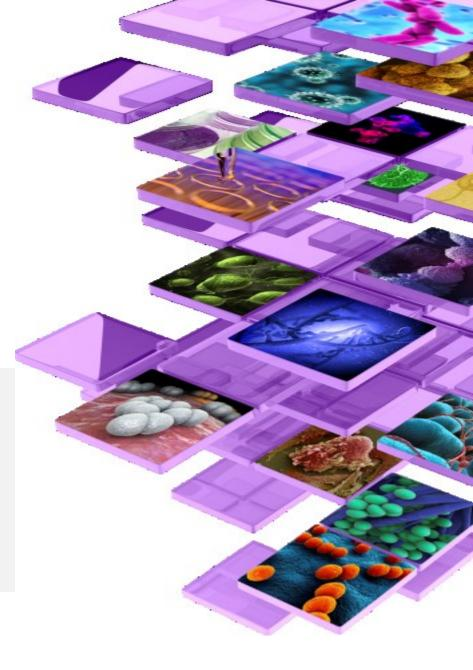
A Tale of 3 Mummies: A Microbiome Analysis of Life in the Peruvian Andes 1,000 Years Ago

Raul Cano, Ph.D. Professor Emeritus, California Polytechnic State University Director, Microbiome Research, ATCC-CTM April 14, 2016





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 one-third with advanced degrees



Established partner to global researchers and scientists





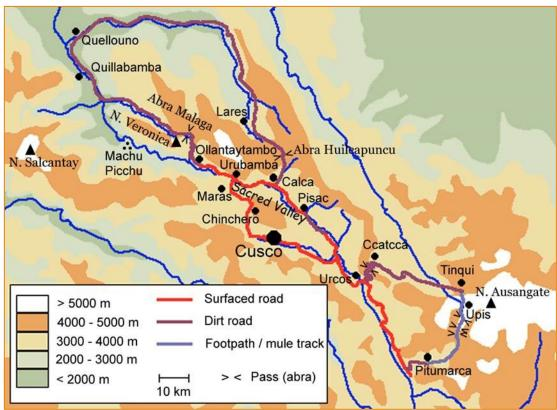




A Tale of 3 Mummies: A Microbiome Analysis of Life in the Peruvian Andes 1,000 years ago

Cuzco, Peru — Geography





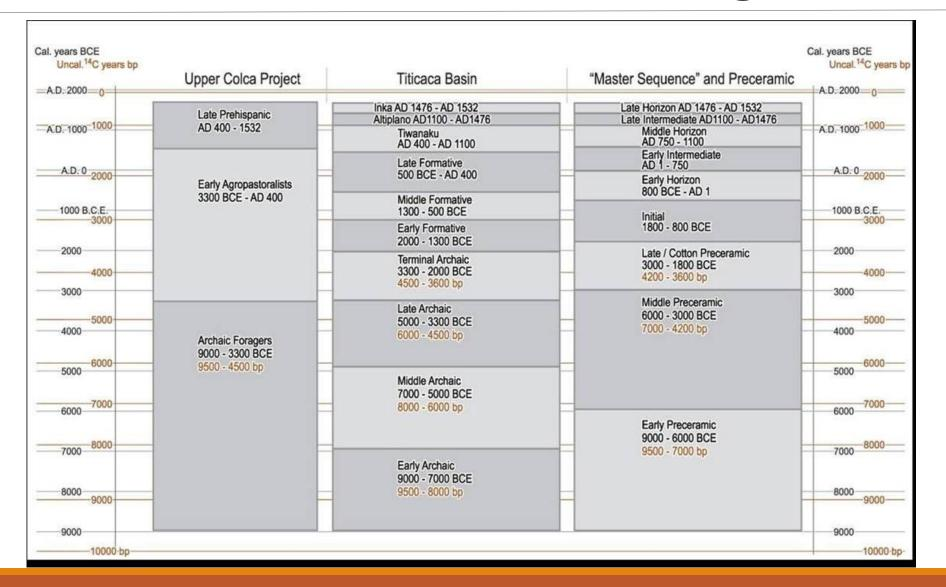


Collaborators and Funding

- Tasha Santiago-Rodriguez: Sample processing, data analysis, and MOST everything else
- Franco Rollo: My friend and gatekeeper
- Gino Fornaciari: Mummy Czar
- Stefania Luciani: Ancient DNA isolation
- Isolina Marota: Continue Franco Rollo's legacy (and lab)
- Gary Toranzos: Public health emphasis



Selected Andean Chronologies





Inca Accomplishments

- Excellent farmers, builders, and managers
- Roads and aqueducts
 - Built > 19,000 miles of roads (over mountains)
 - Built canals and aqueducts to carry water to dry areas
- Advancements in medicine and how to make them use of medicinal plants
- Arts and science
- Accomplished travelers



Inca Diet

- Crops cultivated across the Inca Empire included quinoa, maize, beans, grains, potatoes, sweet potatoes, peppers, tomatoes, peanuts, cashews, squash, cucumber, cotton, carob, and avocado
- Livestock was llama and alpaca herds; meat eaten as charqui
- Drank chicha a lot! (so l'm told)









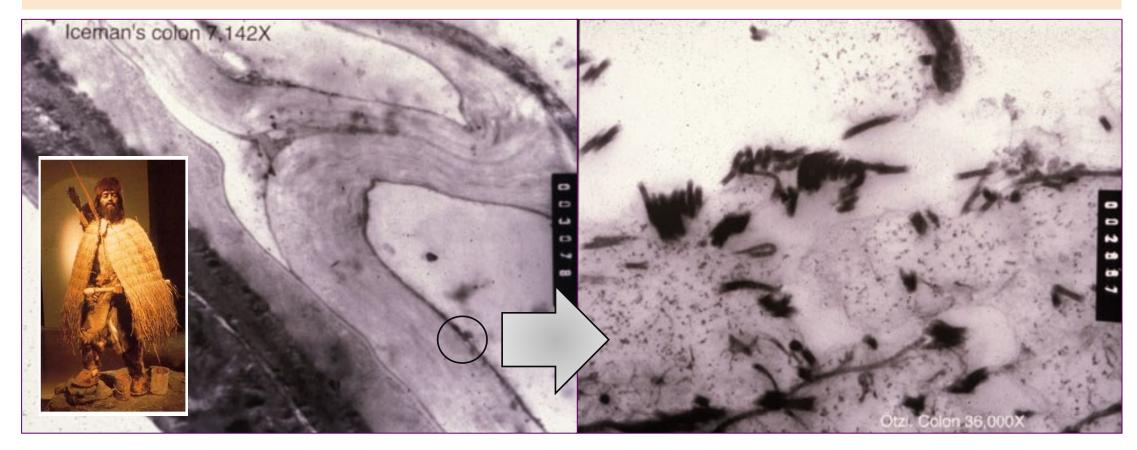
Questions

- Suitability for analysis
- Microbiome/metagenome structure
 - Microbial diversity
 - Metabolic diversity
 - Correlations?
- Medical and cultural aspects
 - AMR/MDR genotypes
 - Infectious diseases
 - Diet



What Got It All Started!

Iceman colon, circa 1998

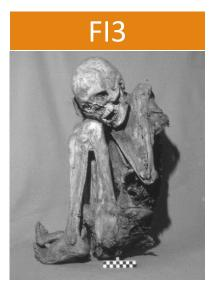




Cast of Characters



The Protagonists: Three Natural Mummies



14th Century male 25-30 years old Good preservation Leishmaniasis



11th Century female 18-23 years old Well preserved Cardiomegaly FI12



14th Century female 20-25 years old Good preservation Bronchopneumonia

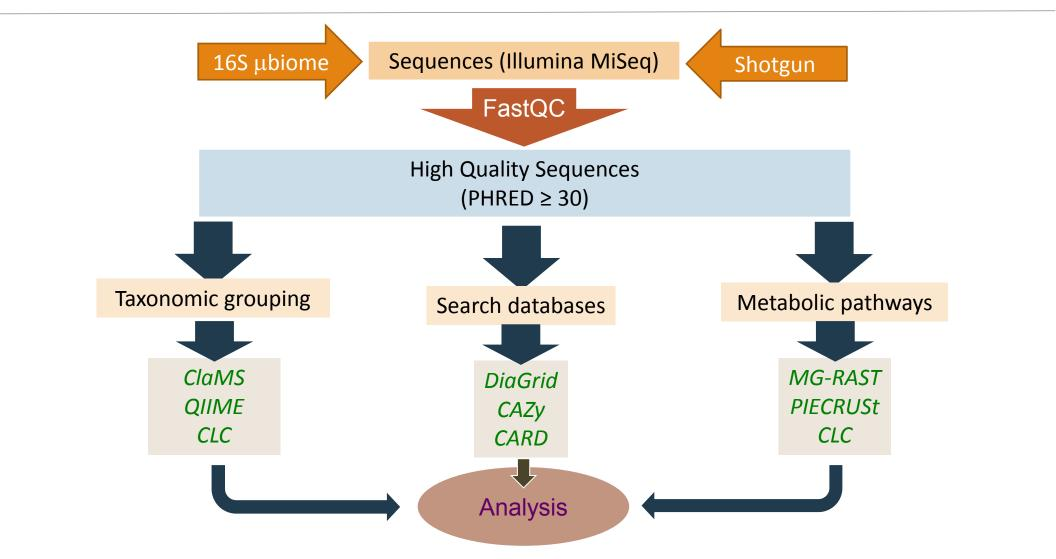


Supporting Cast

	Medici Family			Aragonese Family			
Sample ID	NASD3	NASD12	CM40	NASD22	NADS27	NASD29	
Personage	Francesco I	Ferdinando I	Gian Gastone	Ferrante I d'Aragona	Luigi Carafa	Unknown	
Lifespan	1541-1587	1549-1601	1671-1737	1431-1494	1511-1576	16th Century	
Comments	Grand Duke	Grand Duke	Gran Duke	King of Naples	Prince of Stigliano	?خ	
COD	Malaria	Tertian fever	Small Pox	Colon carcinoma	Natural causes	Cirrhosis	



Analysis Workflow



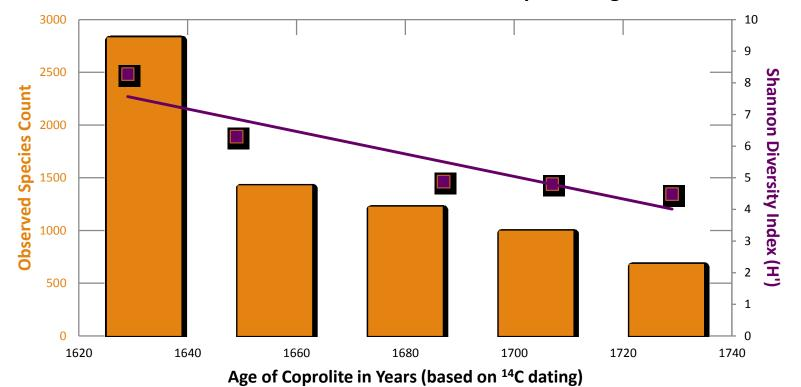


QC/QA of Sequences

ANCIENT DNA MAY BE DAMAGED – SEQUENCES OF POOR QUALITY



Genetic Information is Lost Over Time



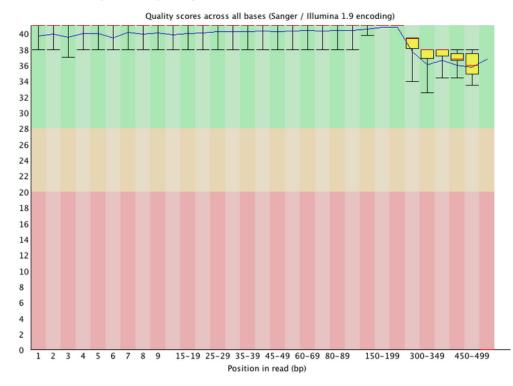
Information Loss as a Function of Coprolite Age



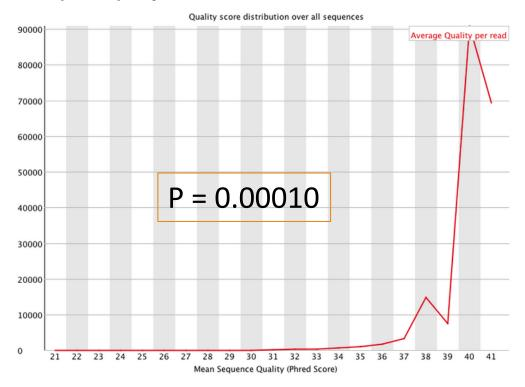
FastQC Results

Assurance of good NGS results

Per base sequence quality



Per sequence quality scores

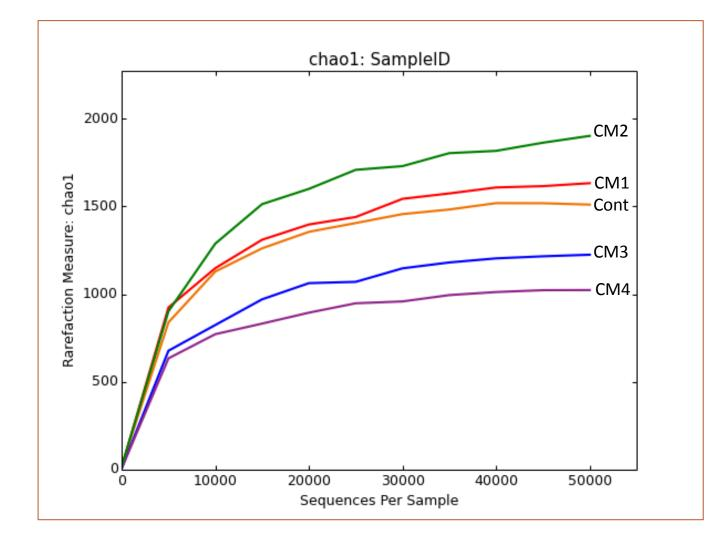




Sampling Depth

Rarefaction Curves

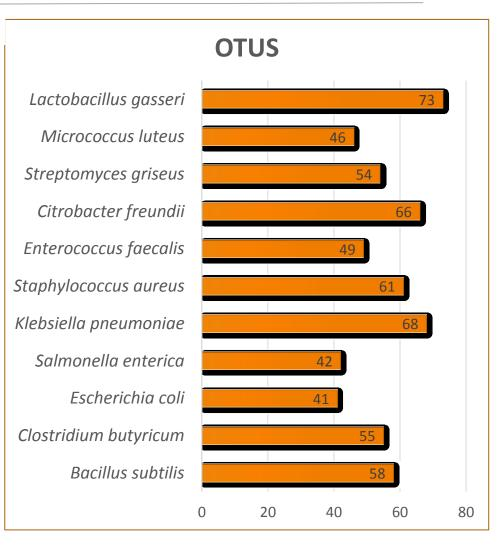
Assurance of sufficient sample size





Mock Community for Assessing DNA Extraction

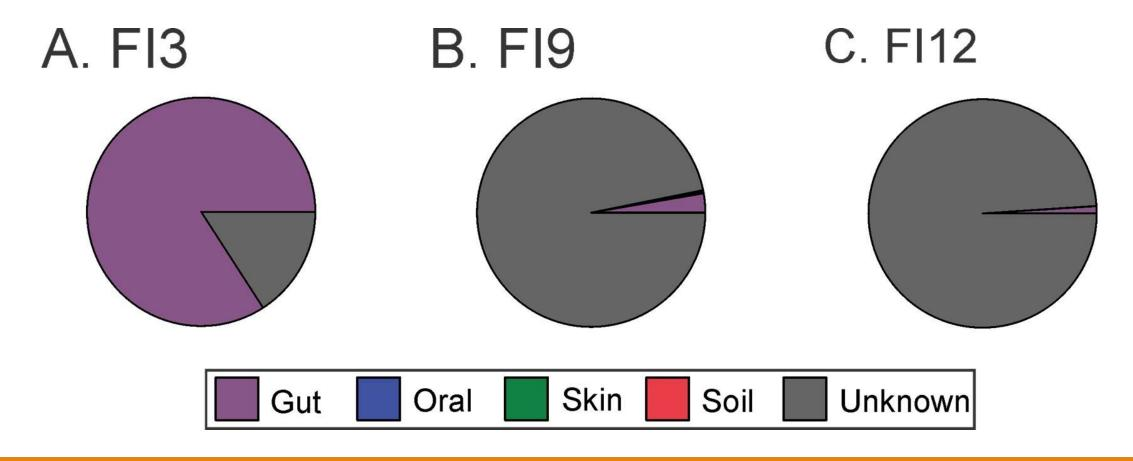
Taxon	ATCC [®] No.	Inoculum CFU/mL
Bacillus subtilis	6501™	1.00E+06
Clostridium butyricum	3627™	1.00E+06
Escherichia coli	11775™	1.00E+06
Salmonella enterica	8388™	1.00E+06
Klebsiella pneumoniae	13883™	1.00E+06
Staphylococcus aureus	12600™	1.00E+06
Enterococcus faecalis	19433™	1.00E+06
Citrobacter freundii	8090™	1.00E+06
Streptomyces griseus	23345™	1.00E+06
Micrococcus luteus	4698™	1.00E+06
Lactobacillus gasseri	33323™	1.00E+06





Source Tracker Results (from QIIME 1.9.0)

Assessing Environmental Contamination





Mummy FI9



Peruvian Mummy from the Museum of Anthropology

- A Peruvian natural mummy of a young woman 20±3 years old from Cuzco (Peru), housed in the Museum of Anthropology and Ethnology of Florence, was autopsied
- The funerary equipment is typical of the Andean highlands, Late Intermediate Period (1000-1476 A.D.)





After preliminary X-rays showing good preservation of the body, the mummy was then autopsied through the posterior thorax to avoid damage





The mummy in its shroud, in strict fetal position



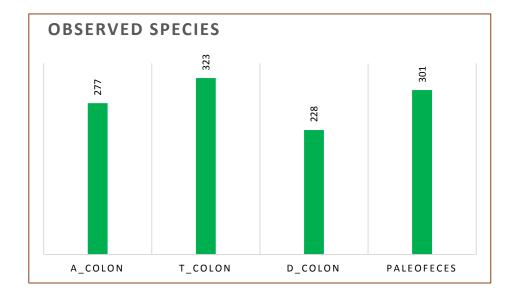
LL X-ray of the mummy

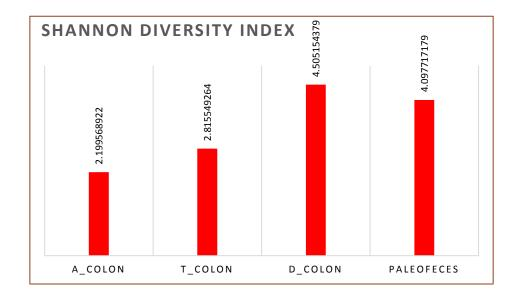


Autopsy by posterior breach



Alpha Diversity: FI9





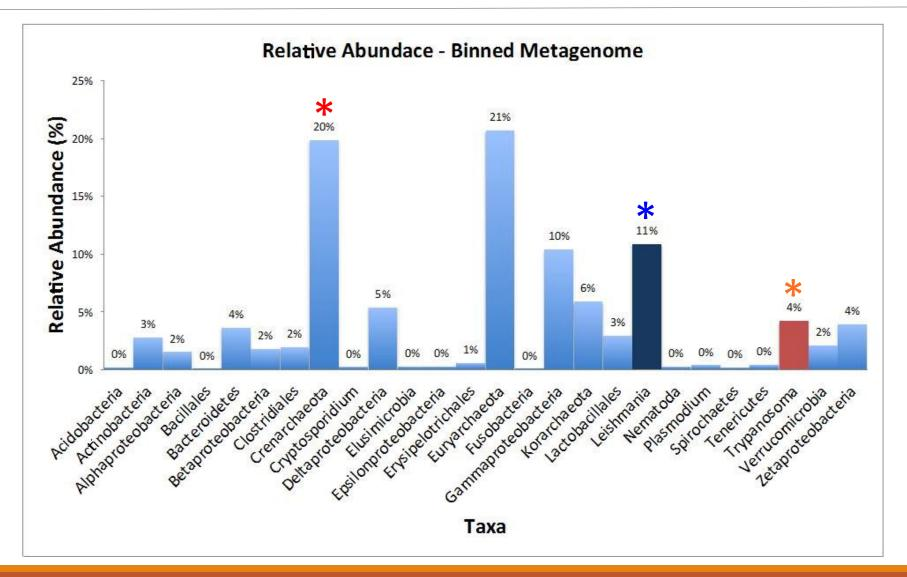


Taxonomic Diversity

Circa 1999: TRFL-P + clor	e and sequence	Circa 2014: NGS + QIIME			
BLAST ID	No. Clones	A. Coprolite	B. Descending colon		
<i>Clostridium botulinum</i> E	18	Turicibacter (89.2%)	Clostridium (81.4%)		
<i>Clostridium botulinum</i> G	2				
Clostridium perfringens	4	C. Transverse colon	D. Ascending colon Turicibacter (19.7%)		
Clostridium sp.	6				
Clostridium algidicarnis	1				
Eubacterium pectinii	2	Clostridium (96.2%)	Clostridium (68.3%)		



Metagenome Binning (ClaMS) – FI9





Paleopathology of Peruvian Mummies



Macroscopically, we (Gino) found a megavisceral syndrome in the form of <u>cardiomegaly</u>, megaesophagus, gastric ectasia, and megacolon, with enormous amounts of feces



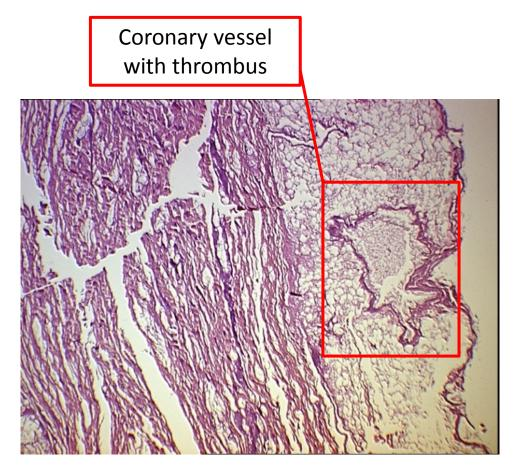
Megacolon with abundant feces



Cardiomegaly



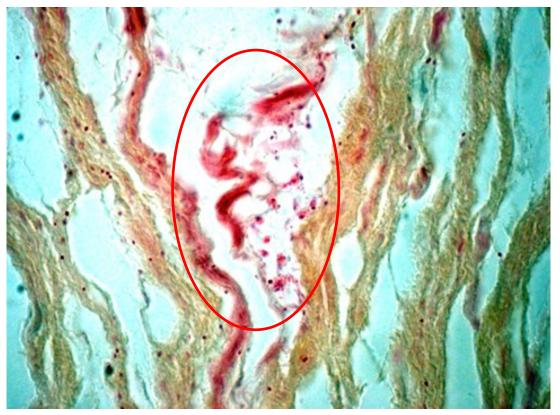
- Light microscopy showed massive fat substitution, fibrosis of the myocardium, and coronary thrombosis
- The mummy mega-visceral syndrome strongly suggested a case of Chagas' disease, chronic phase, caused by the protozoan parasite *Trypanosoma cruzi*



Fat substitution and fibrosis of myocardium (EE, 50X)



- Giemsa staining evidenced rare roundish intratissutal nests, about 15-20 µm large, in the myocardium, containing several ovular formations (1-2 µm) with small nuclei
- The findings correspond morphologically to intratissutal nests of amastigotes of *Trypanosoma cruzi*



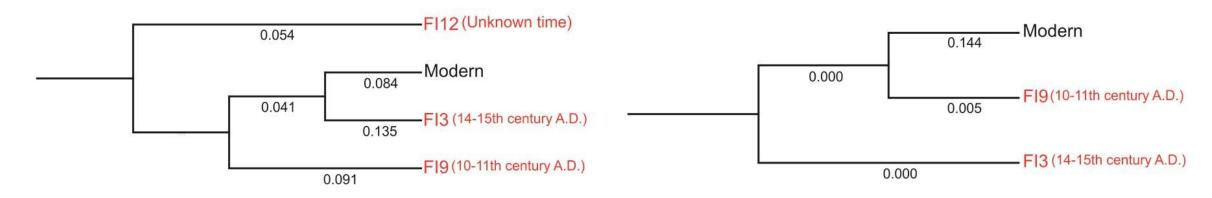
An intramyocardial nidus of amastigotes of *Trypanosoma cruzi* (Giemsa, 1000X)



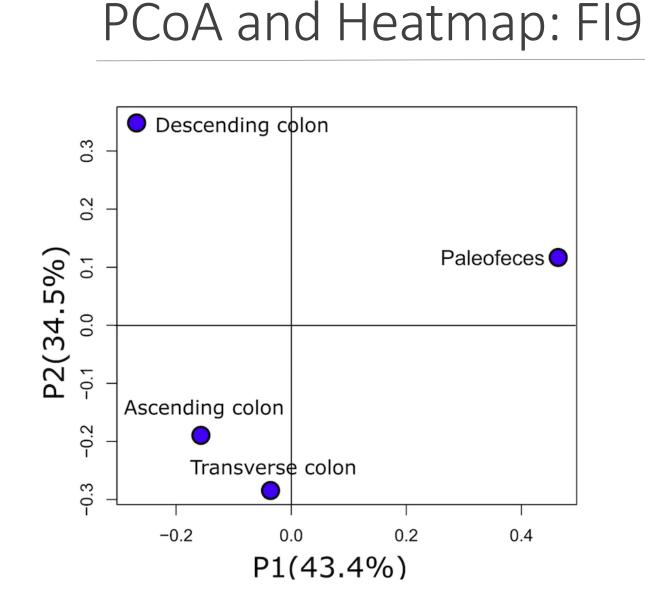
Phylogenetic Studies – Mixed Infections

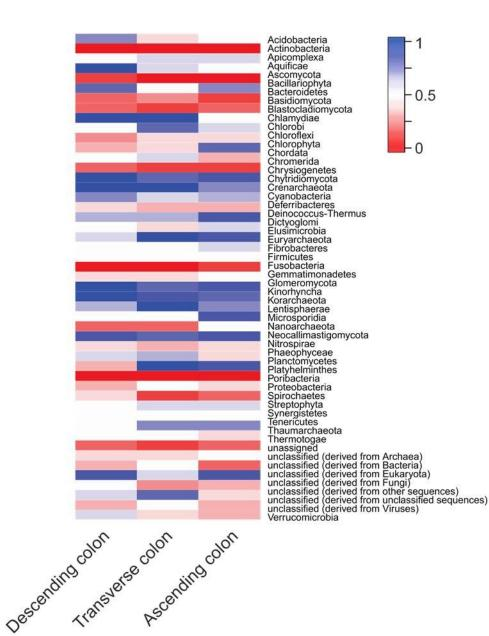
A. Trypanosoma cruzi

B. Leishmania donovani









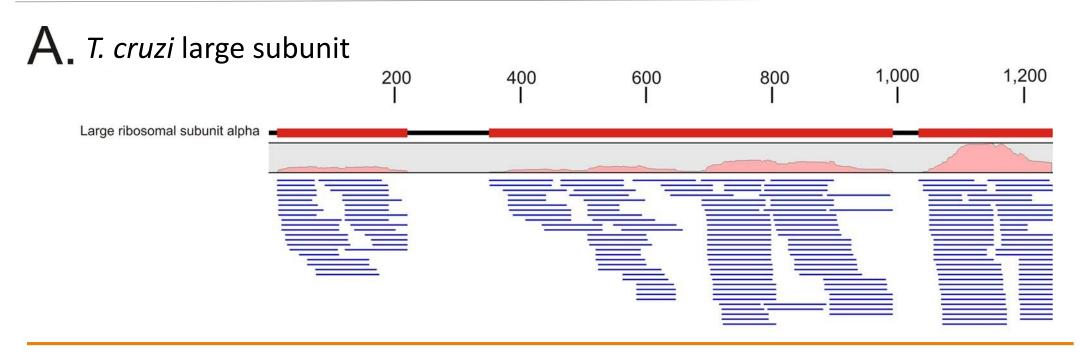


Principal Pathogens Detected

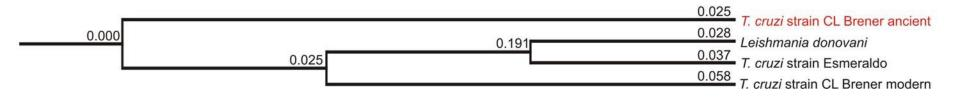
- Trypanosoma cruzi (Chagas disease)
- Leishmania donovani (Leishmaniasis)
- Clostridium
 - botulinum
 - difficile
- Human papillomavirus
 - HPV-21
 - HPV-49



Trypanosoma cruzi



B. Phylogenetic comparisons



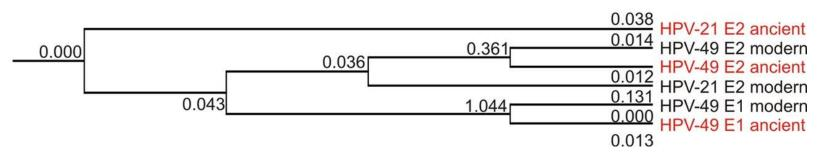


HPV – Sequence Homology and Phylogeny

A. <u>Sequence homology</u>

		1	2	3	4	5	6
HPV-21 E2 ancient	1	100.00	78.21	65.38	65.38	40.08	40.08
HPV-49 E2 modern	2	78.21	100.00	63.68	63.68	38.46	38.87
HPV-49 E2 ancient	3	65.38	63.68	100.00	97.44	34.82	35.22
HPV-21 E2 modern	4	65.38	63.68	97.44	100.00	35.63	36.03
HPV-49 E1 modern	5	40.08	38.46	34.82	35.63	100.00	98.76
HPV-49 E1 ancient	6	40.08	38.87	35.22	36.03	98.76	100.00

B. <u>Phylogenetic relationships</u>





Diet and the Microbiome



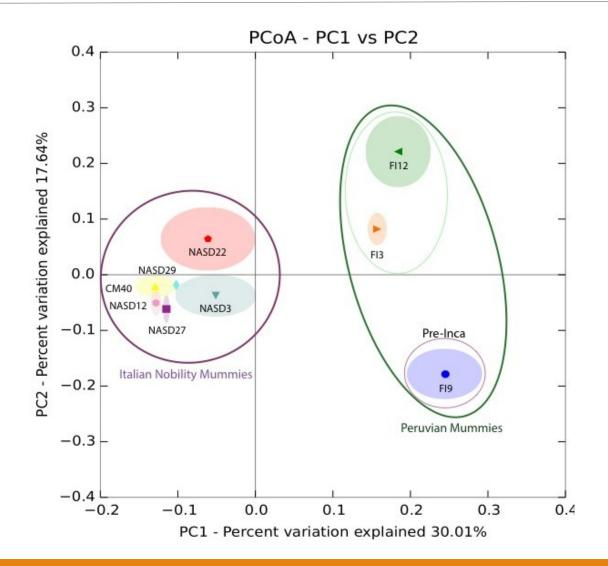
Significant "Hits" in Peruvian Mummies WGS

Crop plant	Number of hits
Соса	517
Beans	1122
Potatoes	865
Peanuts	912
Pepper	877
Tomatoes	922
Corn	589
Rice	435
Avocado	911
Quinoa	339
Sweet potatoes	673

Animals	Number of hits
Alpaca	618
Llama	459
Chimney swift	155
Bony fishes	377

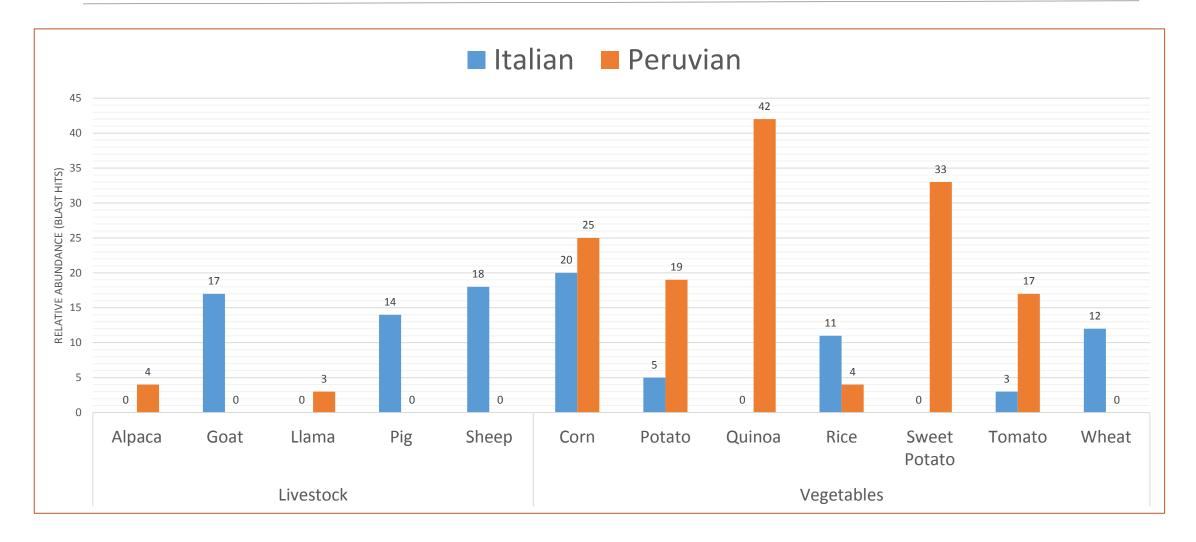


Gut Microbiome Comparisons: Two Diets



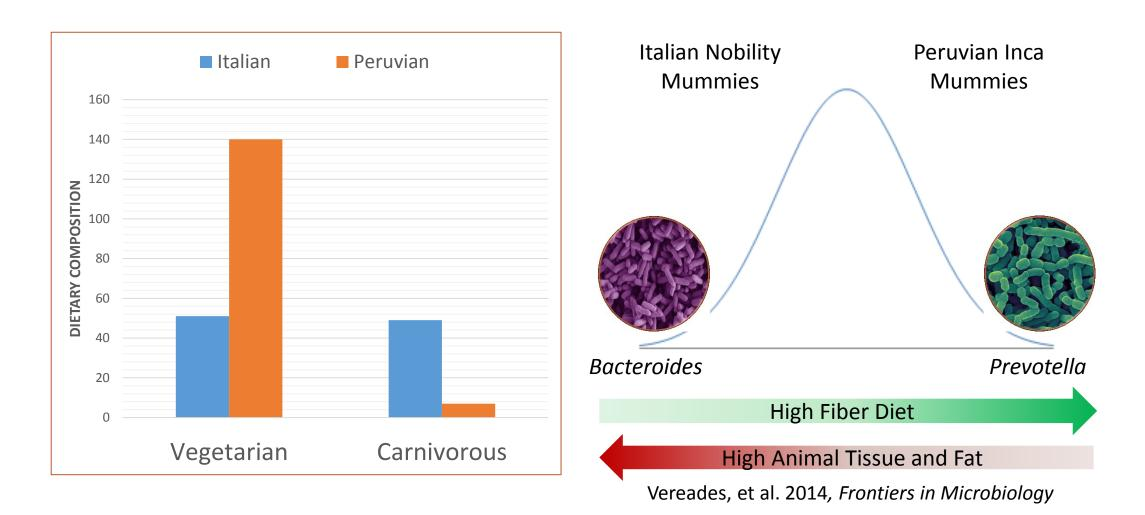


Principal Dietary Composition – Comparative Study



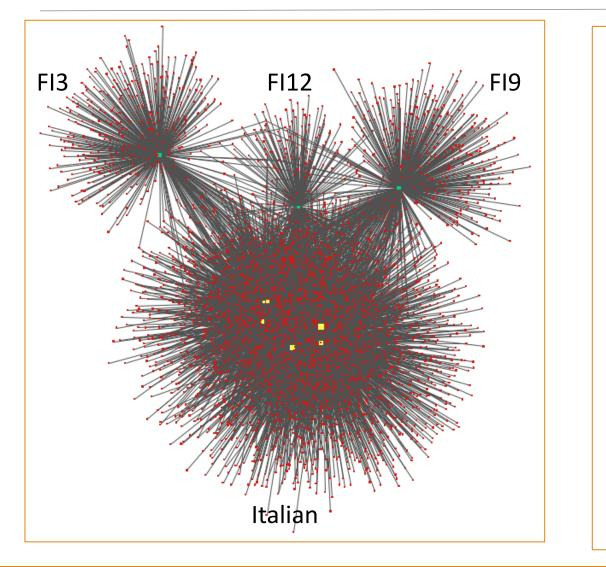


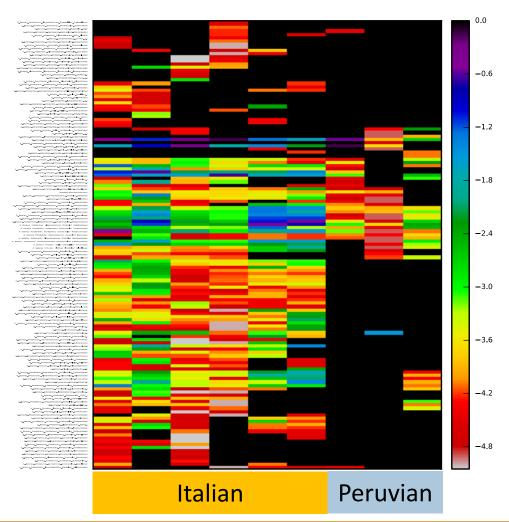
Dietary Impact on Microbial Prevalence





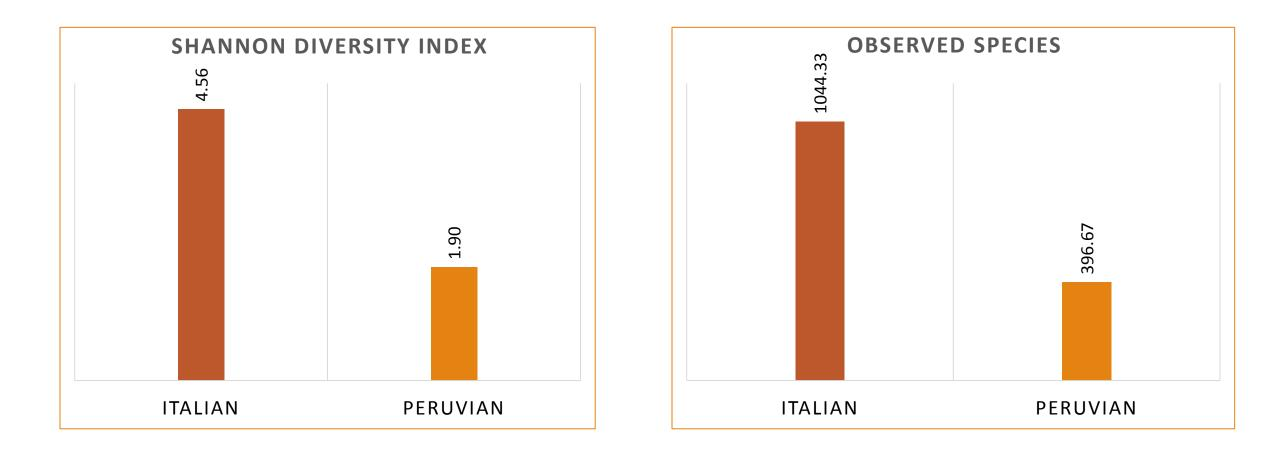
Dietary Impact on Microbial Prevalence ... Continued







Alpha Diversity: Italian v. Peruvian Mummies

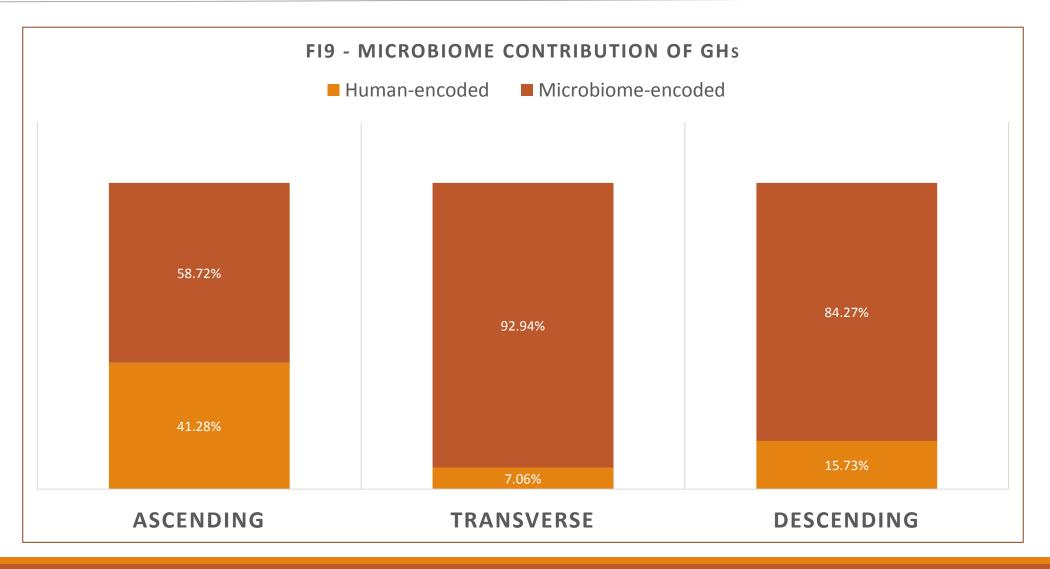




Carbohydrate Metabolism

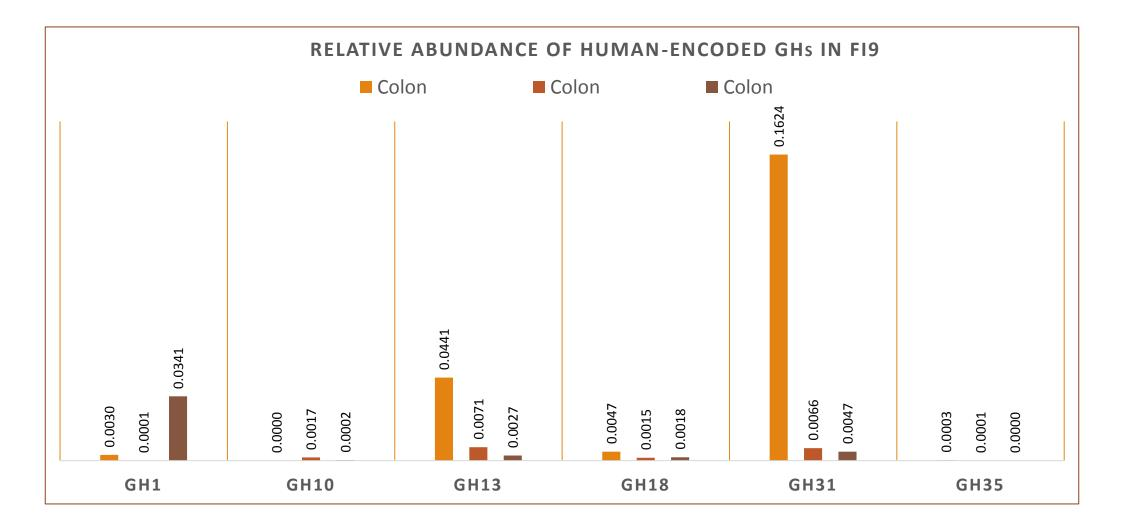


Human-Encoded GH: By Anatomical Site of FI9



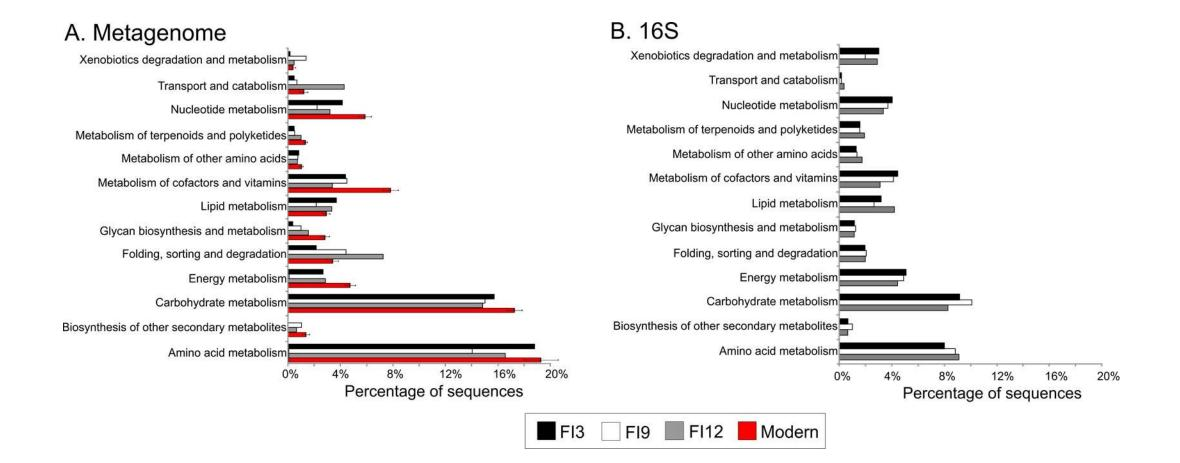


Human-Encoded GH: By Anatomical Site



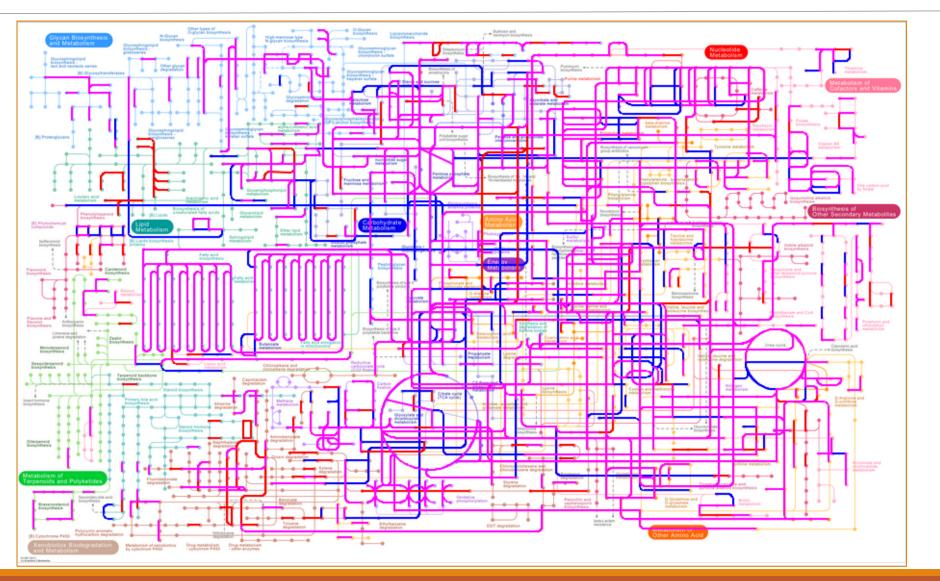


Metabolism (BlastX v PIECRUSt)



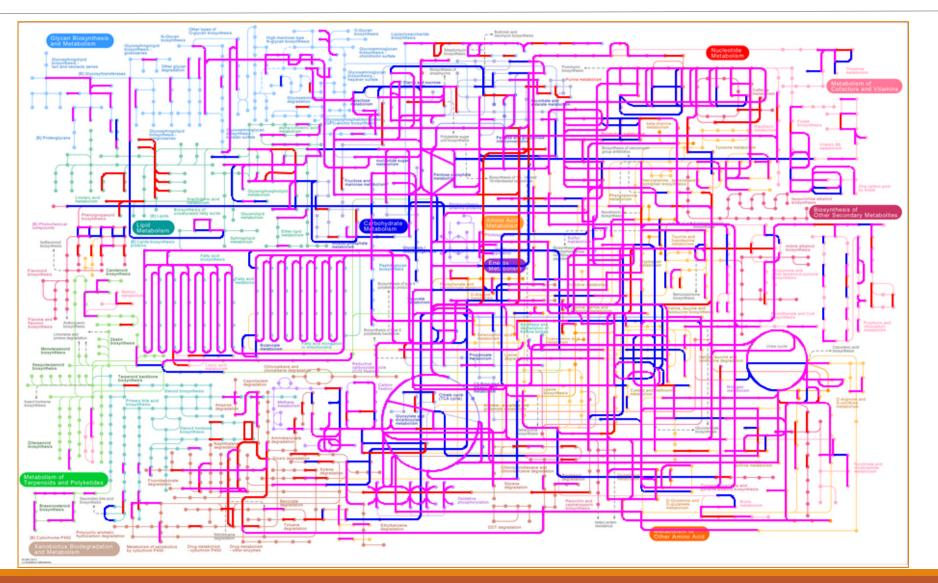


FI3 v. NASD27





NASD 22 v. NASD 27





Final Comments (Finally!)

- Ancient DNA must be suitable for analysis
 - Sample collection and processing
 - Taphonomy
 - Damage analysis recommended
 - Can help answer questions by providing a time reference
 - Mindful of information loss
 - Ask the correct and appropriate questions



Thank you for your interest

Questions?



Thank you for joining today!

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

April 21, 2016
 10:00 AM, 3:00 PM EST
 Steven Budd, M.S., M.B.A., *Product Line Business Specialist*, ATCC
 Best Practices in Cryopreservation

April 28, 2016

 10:00 AM, 3:00 PM EST
 Frank Simione, M.S., *Director, Standards, Standards Resource Organization*, ATCC
 The ATCC Story: A Ninety Year Celebration



Please email additional questions to: tech@atcc.org

