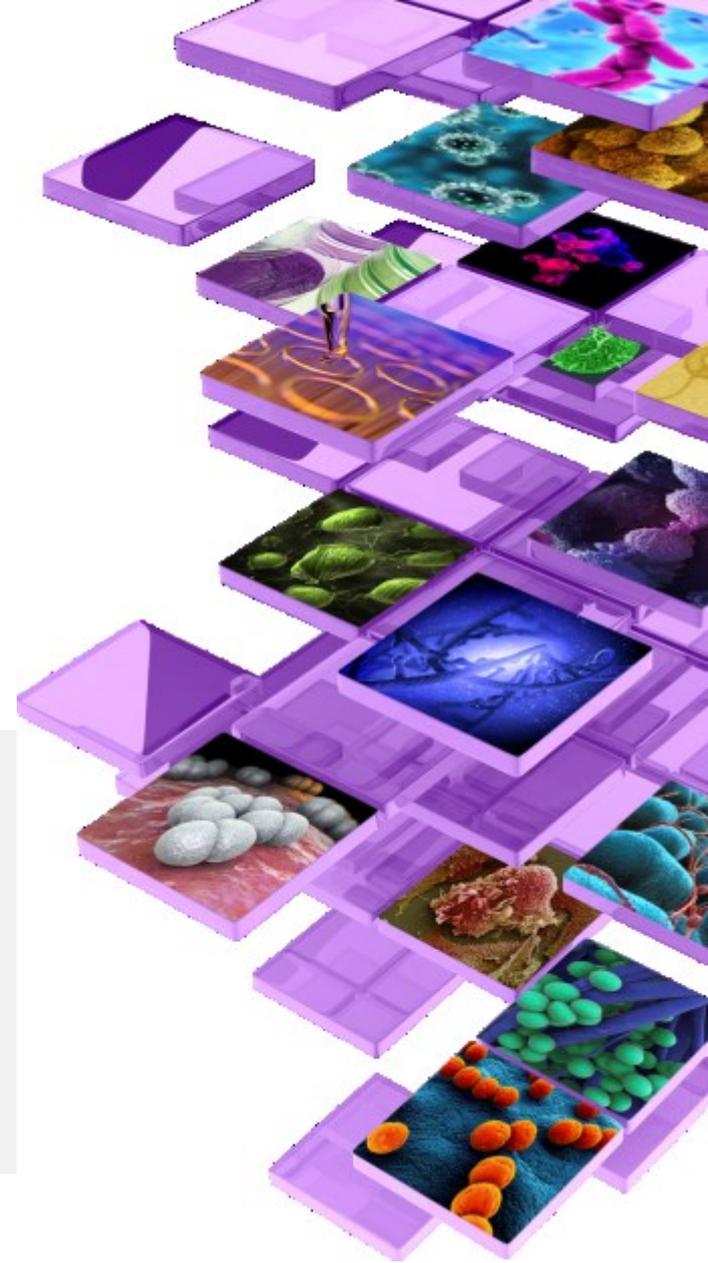


ATCC® Quantitated Nucleic Acids – Empowering Molecular-based Assay Development

Cynthia Long
Product Line Business Manager, ATCC

Fang Tian, Ph.D.
Lead Scientist, ATCC
September 22, 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



Agenda



- **Methods of nucleic acid quantitation**
- **Applications of quantitated nucleic acids**
- **Quantified genomic and synthetic microbial nucleic acids**
- Precision medicine and unmet needs in genetic tests
- ATCC human genomic DNAs as control materials
- ATCC quantified human genomic DNAs
 - Gene mutation allelic frequencies
 - Gene copy numbers

Utility for molecular testing



- Improved sensitivity
- Improved specificity
- Time to results
- Ability to multiplex

Analytical sensitivity, specificity, & interference

Analytical sensitivity

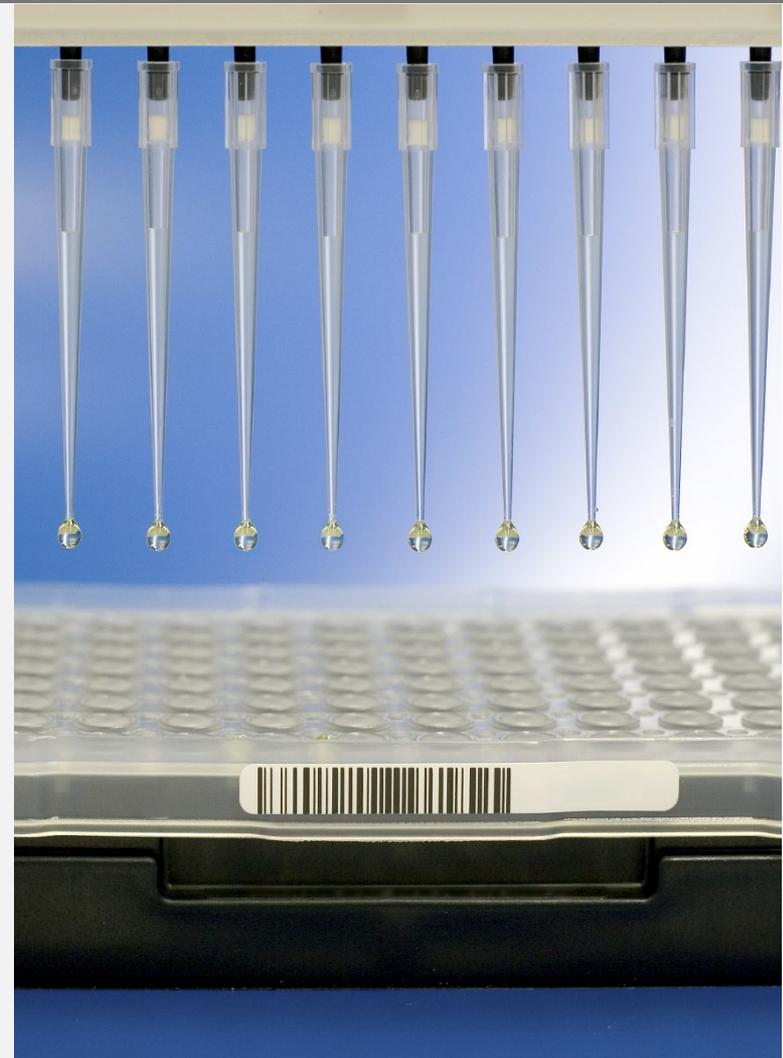
- Limit of detection (LoD)
- Reactivity

Analytical specificity

- Cross reactivity

Interference

- Clinical specimens



Controls for nucleic acid-based assays



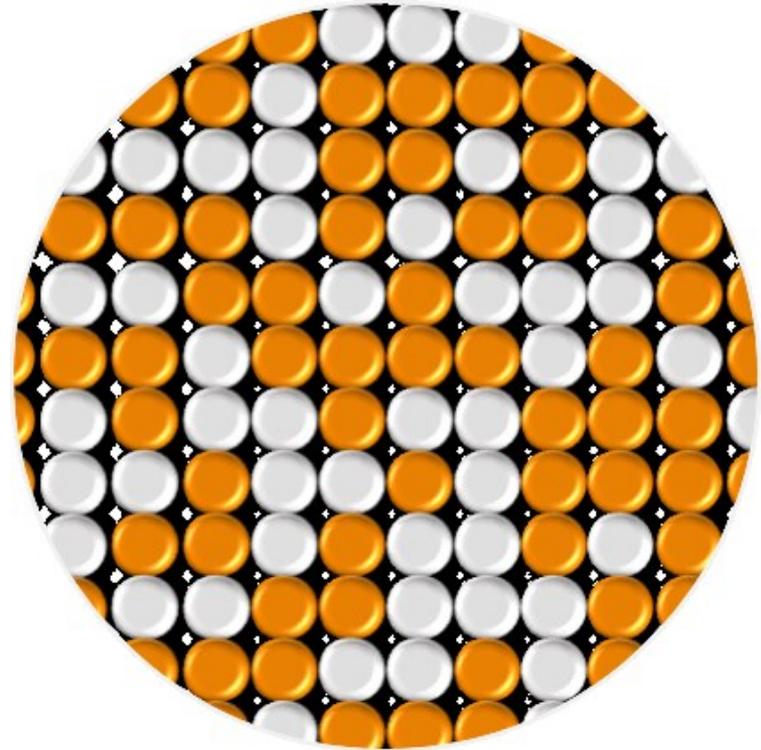
Used for verification of assay performance in your analytical and clinical studies:

- **Negative control:** To rule out contamination
- **Positive control:** To verify PCR reagents & instrument are running properly
- **Internal control:** To verify that a sample result is not caused by PCR inhibitors
- **External control:** To verify that lysis & extraction processes are functioning properly
- **Inclusivity/Exclusivity**

Advantages of digital PCR

Allows for absolute quantification of nucleic acids

- High precision and accuracy
- Target-specific quantification
- Copy number of individual genes
- Cost effective
- No need to generate cloned standards for a standard curve



Quantitative nucleic acid standards

Salient features

- Fully authenticated & characterized
- Quantitated by Droplet Digital™ PCR (ddPCR™)
- Compatible with several lab-developed and commercially available assays
- BSL-1 ready-to-use control

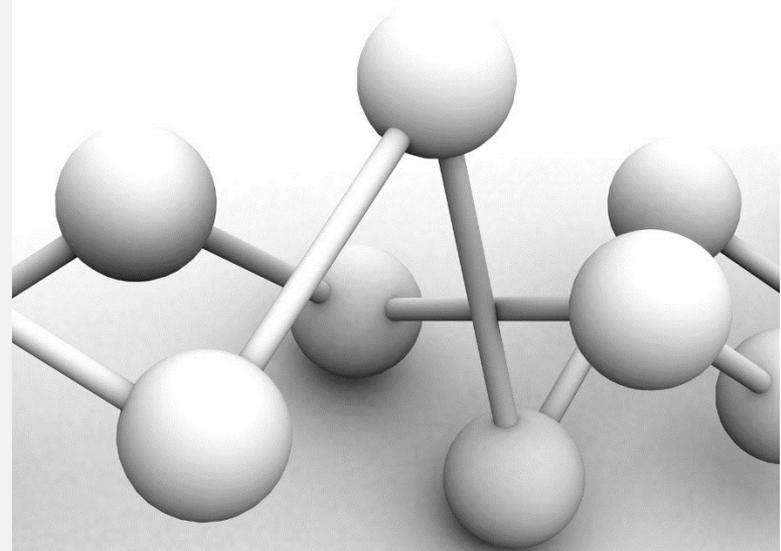
Applications

- Generation of a standard curve
- Positive control for molecular-based assays
- Independent standard for validation and verification studies
- Monitoring assay-to-assay and lot-to-lot variation
- New assay development
- Limit of detection studies

Quantitative nucleic acid standards

The collection encompasses over 60 quantitative preparations

- Synthetic nucleic acids
- Genomic nucleic acids
- Certified reference materials



ATCC Synthetic Molecular Standards

Can be used as a positive control for:

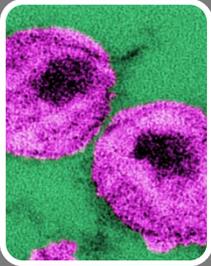
- Difficult-to-culture or unculturable strains
- Strains requiring BSL-3 containment
- Strains on the commerce control list

Advantages of synthetic nucleic acids:

- Eliminate the need to culture microorganisms
- Can be used in a BSL-1 facility
- No shipping restrictions
- Manufactured under ISO 13485:2003
- Quantified using ddPCR™
- Useful for monitoring assay-to-assay or lot-to-lot variation

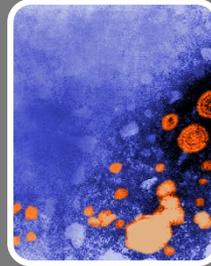


ATCC Synthetic Molecular Standards



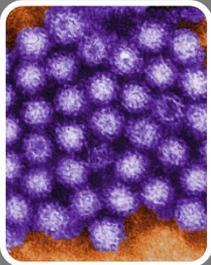
Blood-borne Disease

- BK virus
- Hepatitis B & C virus
- Epstein-Barr virus
- Human immunodeficiency virus



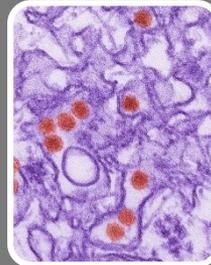
Sexually Transmitted Infections

- Hepatitis B & C virus
- Human immunodeficiency virus
- Human papillomavirus 16 & 18
- Mycoplasma genitalium*
- Treponema pallidum*



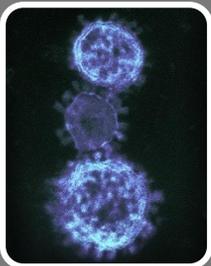
Enteric Disease

- Astrovirus
- Cyclospora cayentanensis*
- Norovirus GI & GII
- Sapovirus



Vector-borne Disease

- Chikungunya virus
- Dengue virus types 1-4
- Eastern equine encephalitis virus
- St. Louis encephalitis virus
- West Nile virus
- Zika virus

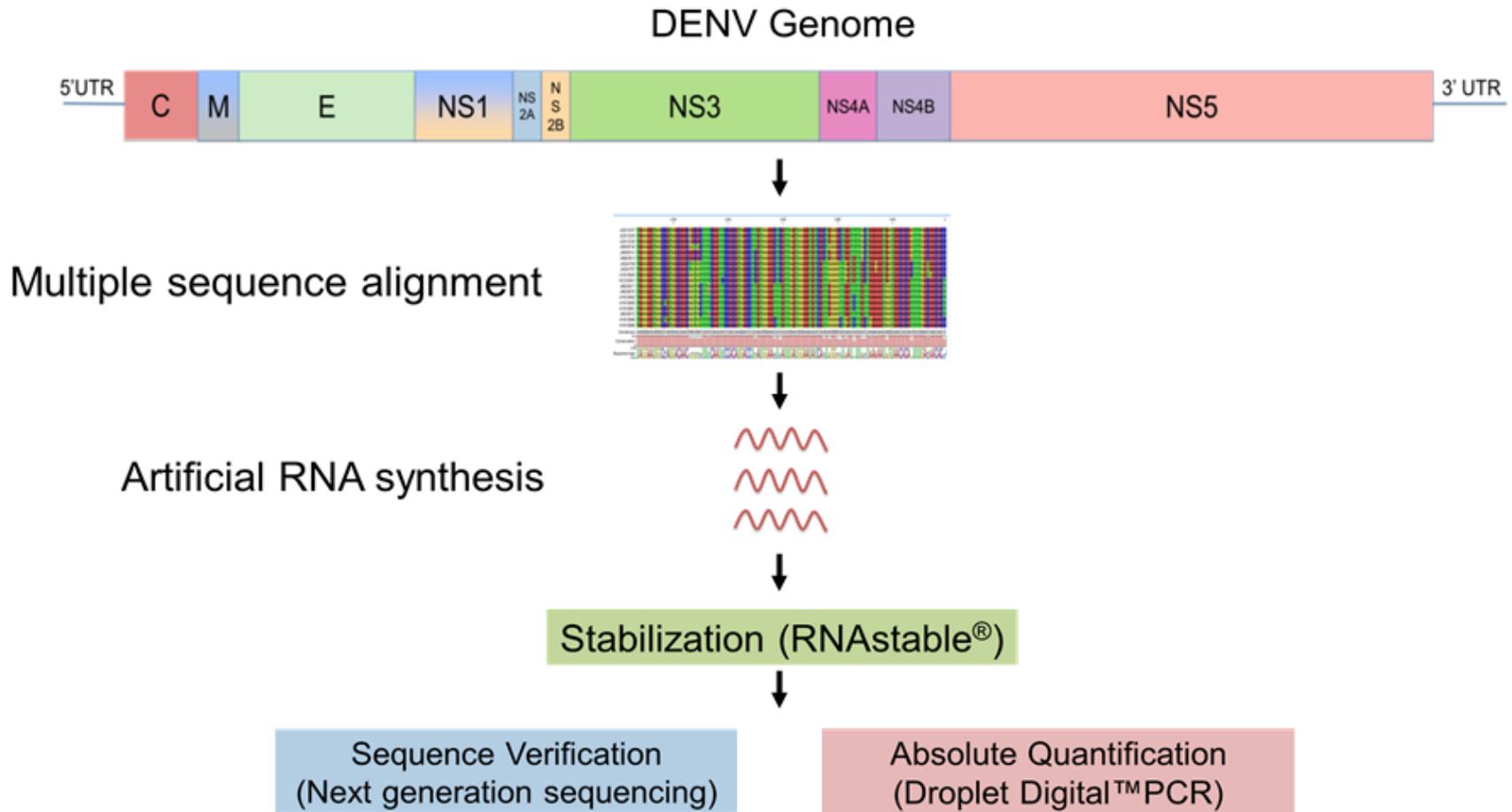


Respiratory Disease

- Human Bocavirus
- MERS-CoV
- Human metapneumovirus

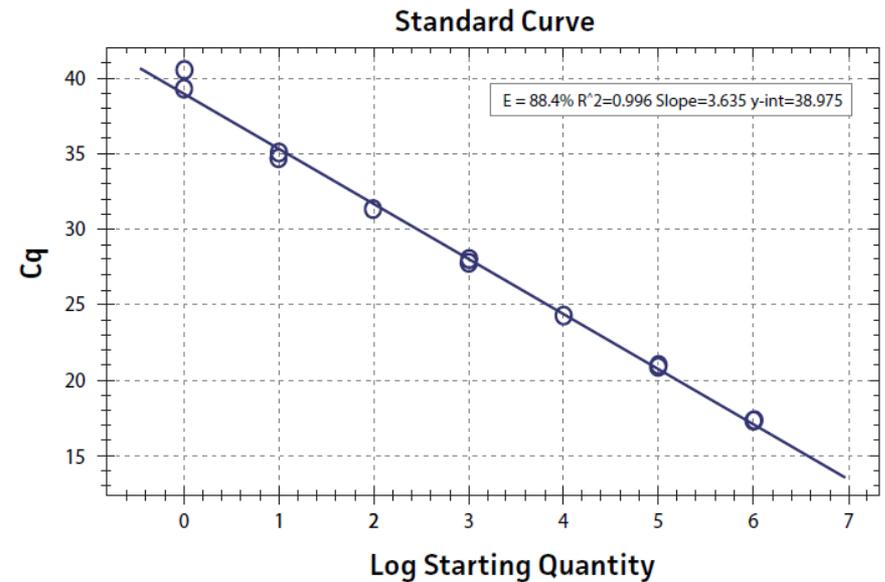
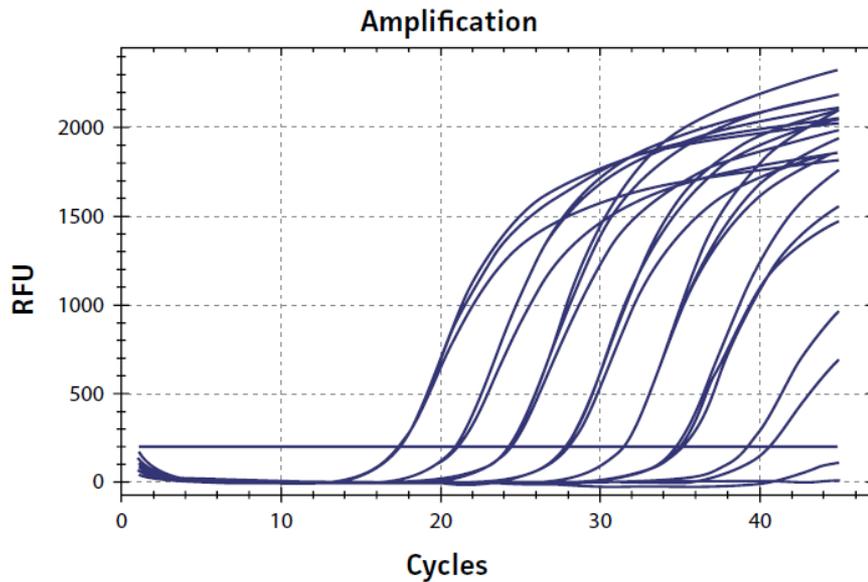
Available online at
www.atcc.org/synthetics

Synthetic Dengue virus RNA



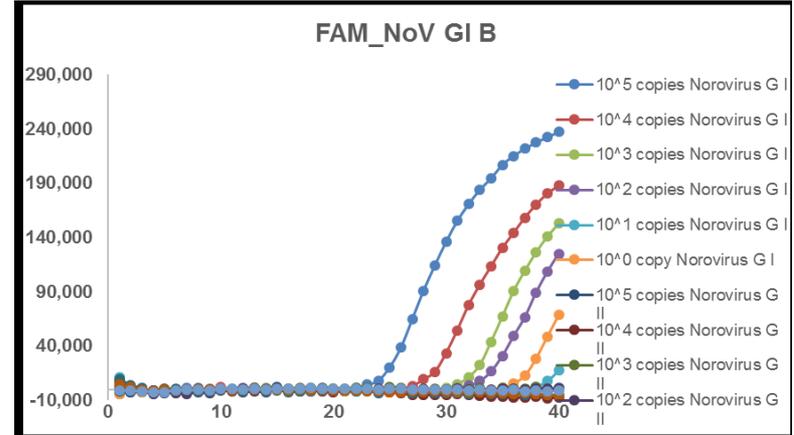
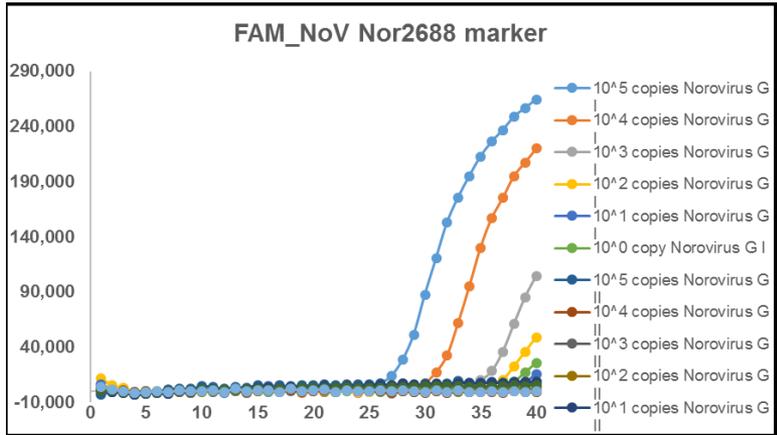
Synthetic Dengue virus RNA

Generation of standard curves using the DENV-4 Molecular Standard



Primer and Probe		DENV-1	DENV-2	DENV-3	DENV-4
CDC Assay	Slope	-3.244	-3.277	-3.315	-3.642
	R ²	0.990	0.996	0.987	0.996
Waggoner Assay	Slope	-3.536	-3.535	-3.705	-3.775
	R ²	0.991	0.997	0.989	0.996

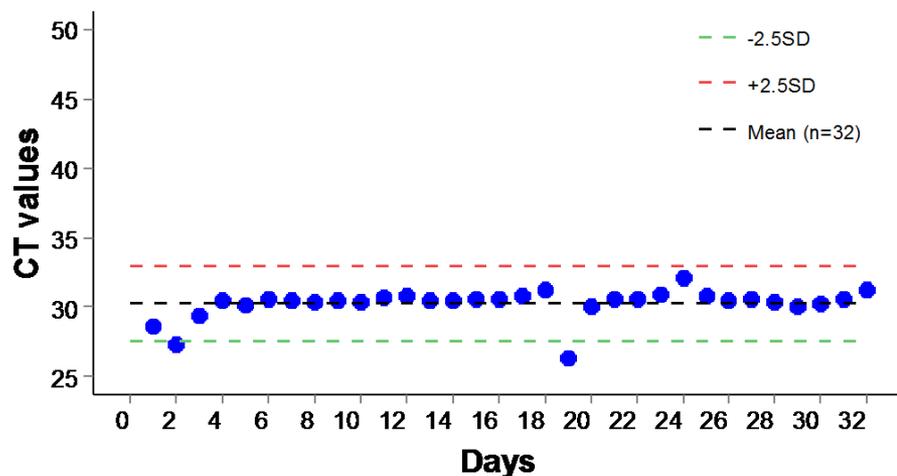
New assay development with Norovirus GI and GII



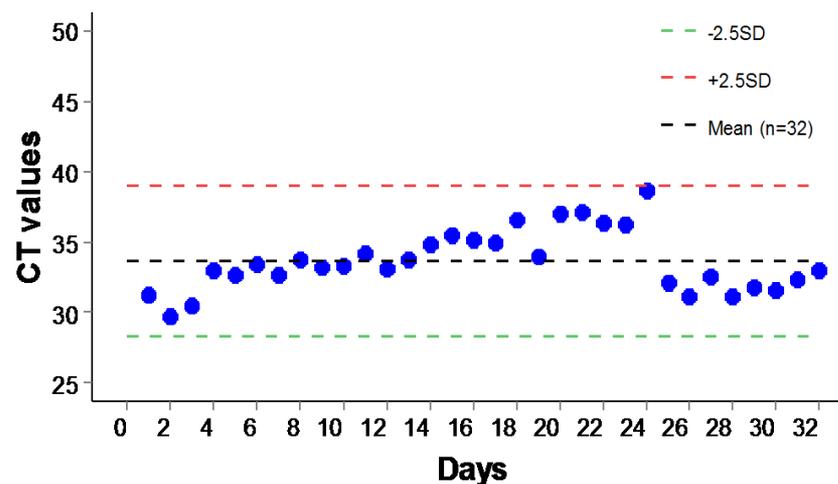
Primer	Probe Target	Limit of Detection	
		NoV GI Template	NoV GII Template
Set 1	FAM_NoV GI Nor2688	500 copies	-
	Cy5_NoV GII Nor21932	-	500 copies
Set 2	VIC_NoV GI A	5 copies	-
	FAM_NoV GI B	5 copies	-
	ROX_NoV GII C	-	5000 copies

Independent assay validation

NoV-GI control



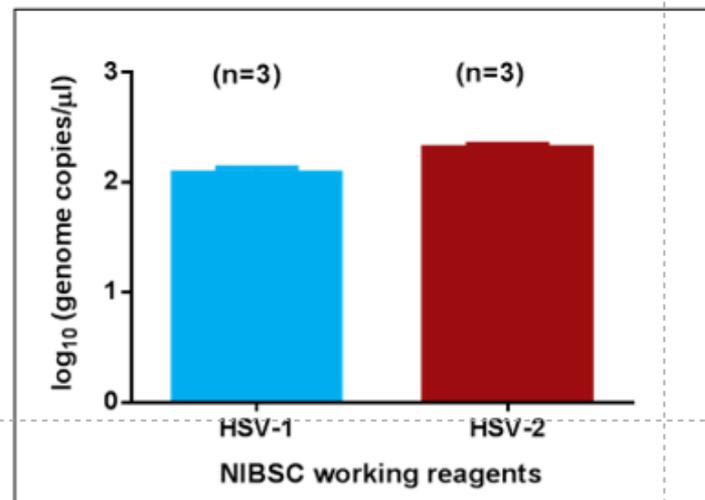
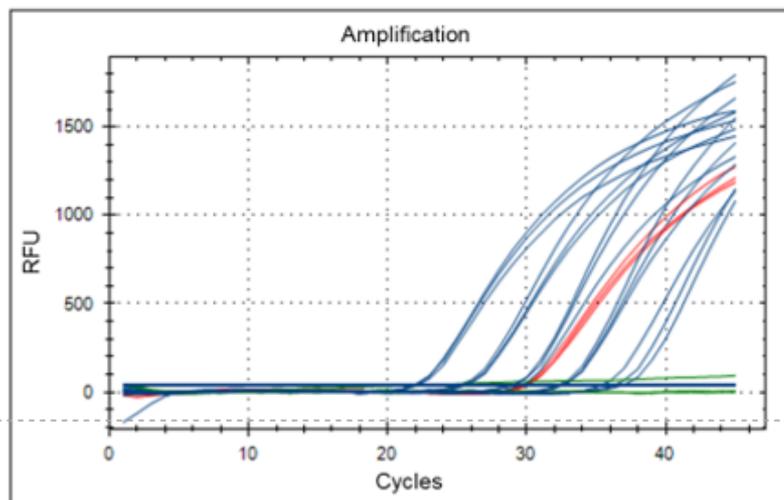
NoV-GII control



ATCC Synthetic Molecular Standards	Average CT	Standard Deviation (SD)	Co-efficient of Variation (CV)	2.5 SD Range
NoV GI	30.27	1.09	3.6%	27-34
NoV GII	33.69	2.15	6.4%	29-39

Utility of quantitated molecular standards as controls

Quantification of NIBSC working reagents for HSV-1 and HSV-2 using ATCC quantitative molecular standards



National Institute of Biological Standards and Control
(NIBSC) working reagents for HSV-1 and HSV-2

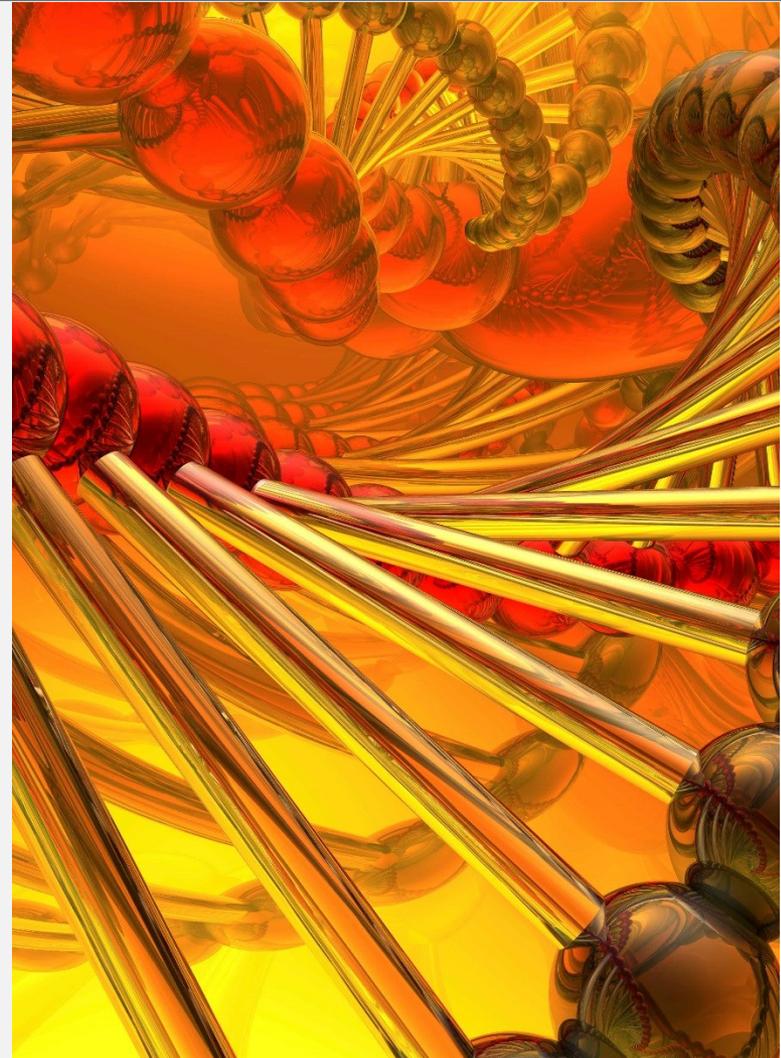
CRM Quantitated Mycoplasma Genomic DNA

Key features:

- Certified reference material
- Quantitated - Genome copy number is based on the quantification of the 16S rRNA gene from nine averaged samples using ddPCR™ (1.0x10⁶-1.0x10⁷ genome copies/μL)
- Extracted from titered mycoplasma reference materials

Use as a quantitative external control for:

- Inclusivity/exclusivity testing
- Establishing limits of detection (LoD)
- Verification or comparison of test methods
- Other molecular applications



Agenda



- Methods of nucleic acid quantitation
- Applications of quantitated nucleic acids
- Quantified genomic and synthetic microbial nucleic acids
- **Precision medicine and unmet needs in genetic tests**
- **ATCC human genomic DNAs as control materials**
- **ATCC quantified human genomic DNAs**
 - Gene mutation allelic frequencies
 - Gene copy numbers

Precision medicine

Precision Medicine Initiative®

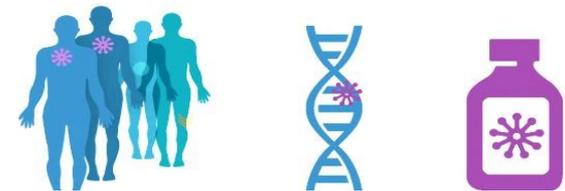
- Prevention and treatment strategies that take individual variability into account
 - Large-scale biologic databases
 - Molecular approach
 - Near-term focus on cancers
 - Longer term aim for a wide range of disease



The website of the White House (www.whitehouse.gov)

NATIONAL CANCER INSTITUTE PRECISION MEDICINE IN CANCER TREATMENT

Discovering unique therapies that treat an individual's cancer based on the specific genetic abnormalities of that person's tumor.

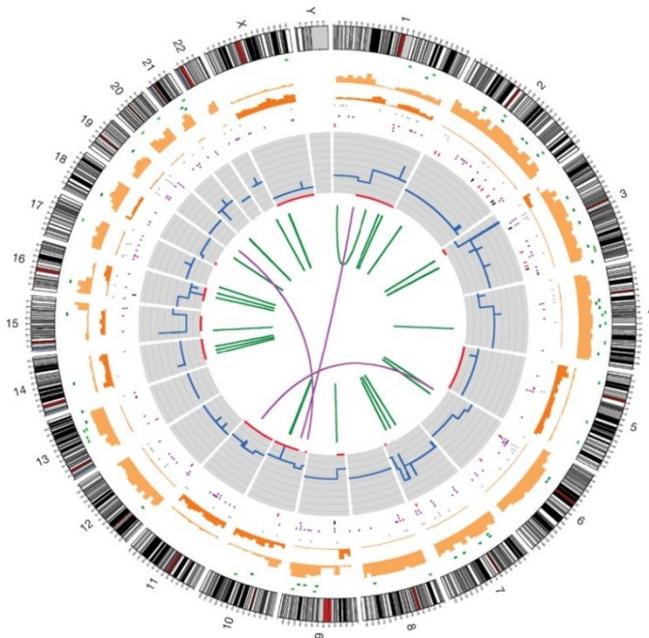


The website of the National Cancer Institute (www.cancer.gov)

Genetics of cancer

Genetic changes and cancer

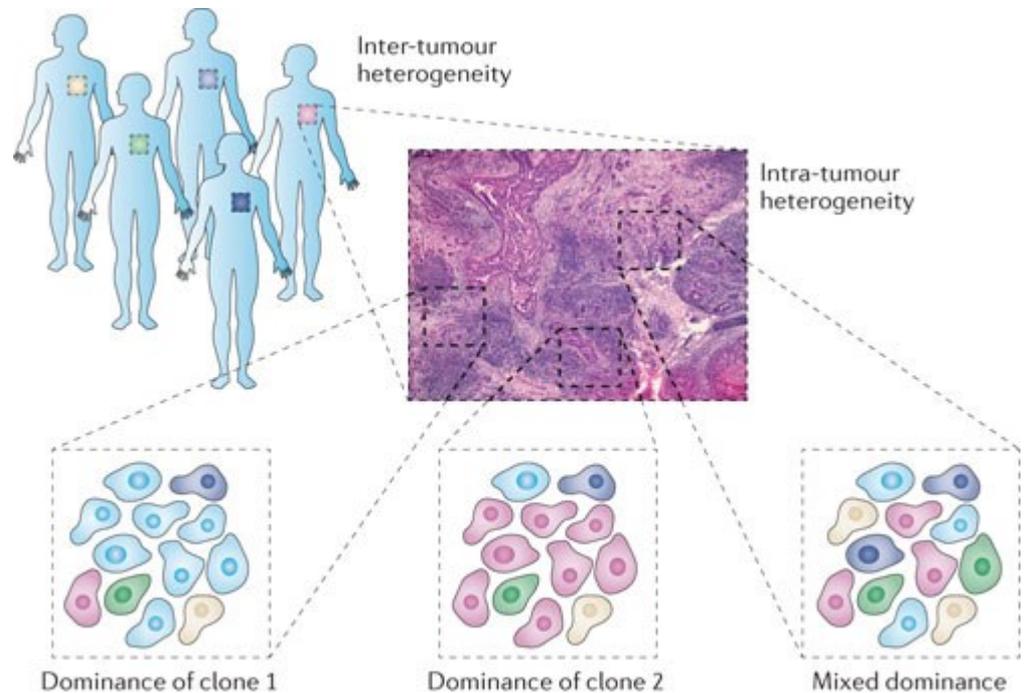
The catalogue of somatic mutations from a human cancer genome



Nature 463(7278): 191-196, 2010.

Heterogeneity of cancer

Every cancer is unique

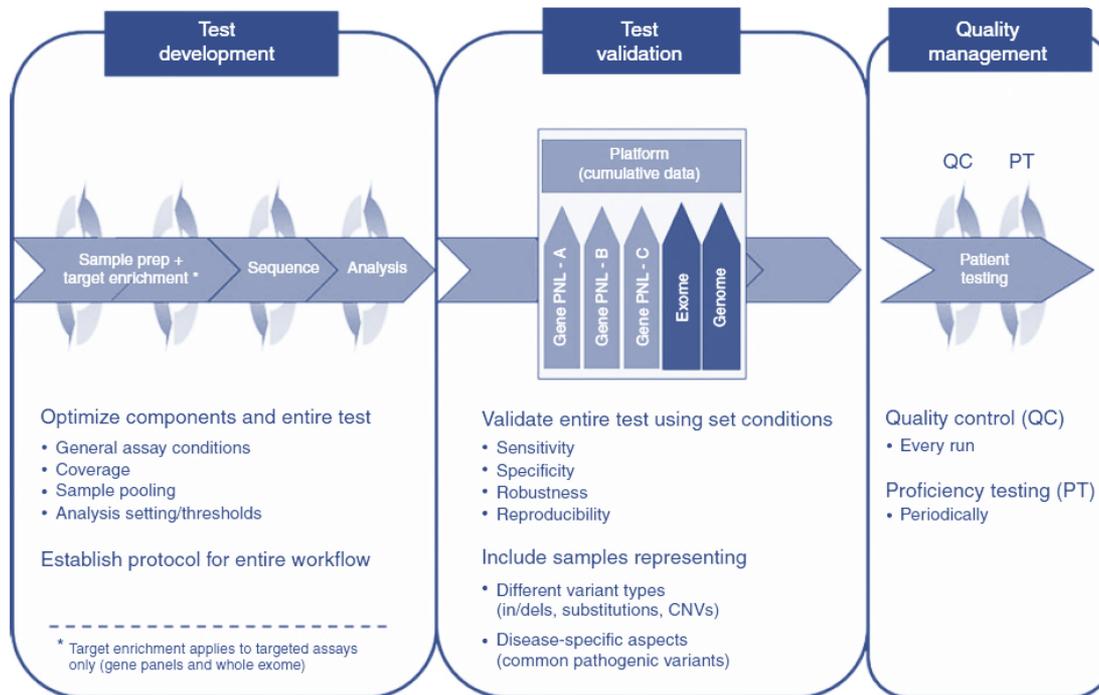


Nature Reviews Cancer 12: 323-334, 2012.

Nature Reviews | **Cancer**

ACMG guidelines for NGS

ACMG clinical laboratory standards for next-generation sequencing



Using reference materials

- Test validation
- Quality control
- Proficiency testing

A need for established, fully characterized, globally accepted reference materials

Genet Med 15(9): 733-747, 2013.

Unmet need - Reference materials

What is a reference material?

A material or substance, one or more of whose property values are sufficiently homogeneous and well established to be used for the calibration of a measuring system, the assessment of a measurement procedure, or for assigning values to materials (ISO 15195:2003)

A variety of reference materials, including:

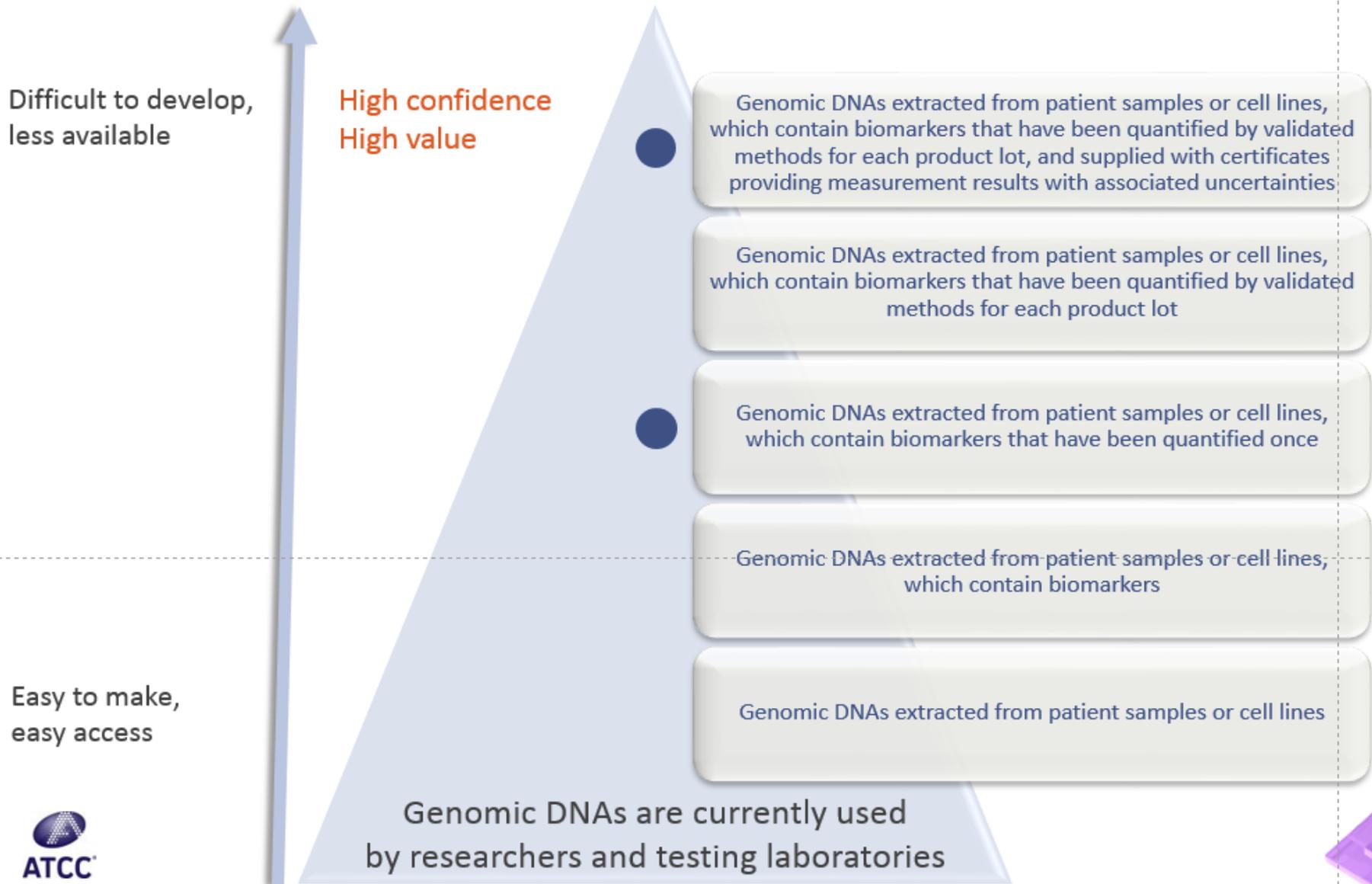
- Certified reference materials
- Standard reference materials
- Calibrators
- Characterized genomic nucleic acids

Reference material properties

- Qualitative
- Quantitative

A need for established, fully characterized, globally accepted reference materials

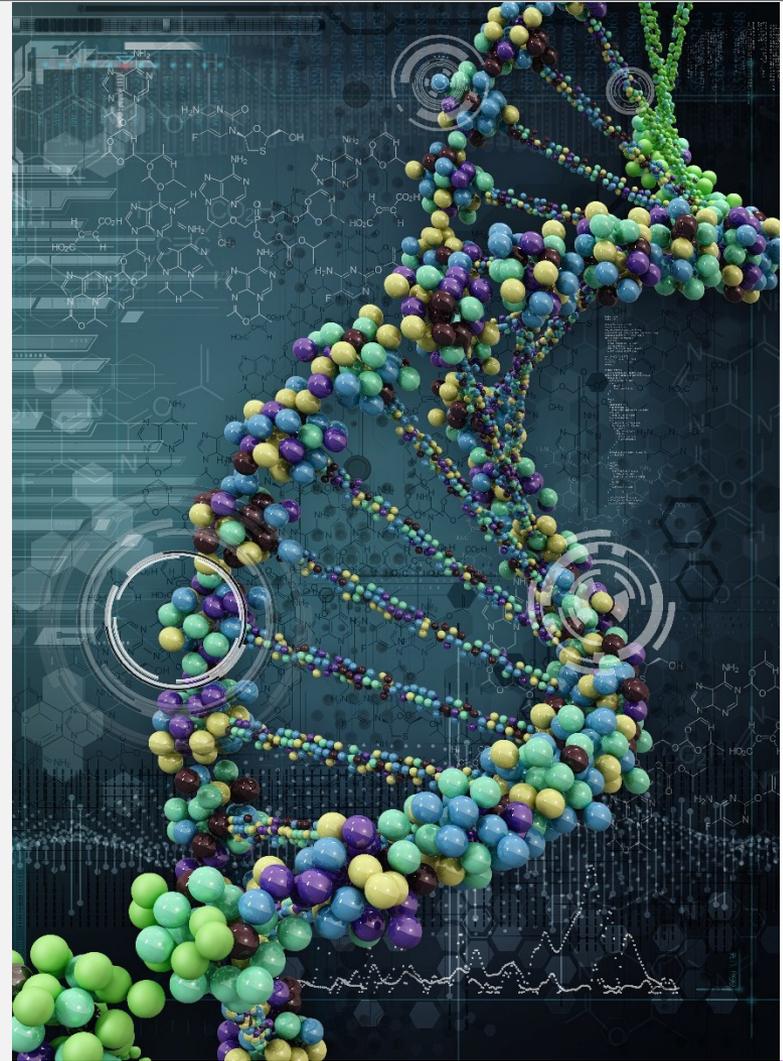
Example of various genomic DNA materials



ATCC human Genomic DNA (gDNA) control materials

Purified human gDNAs with oncology biomarkers

- Fully authenticated
- Short Tandem Repeat (STR) testing used to avoid contamination or misidentification
- Characterized genetic alterations
- Faithfully capture human cancer genome
- Reproducible results



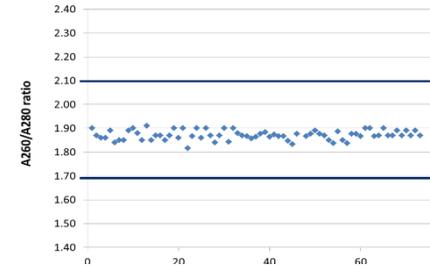
High quality gDNA for molecular tests

- Quantity
- Integrity
- Purity
- Identity
- Functional testing

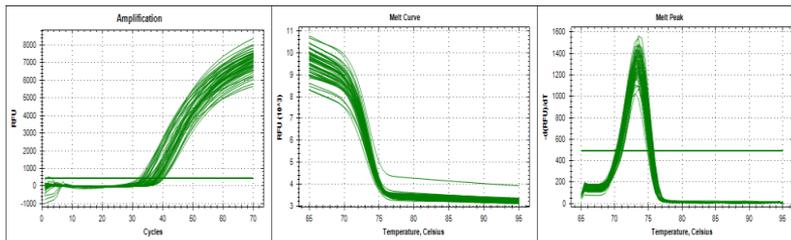
PicoGreen® dsDNA Quantitation



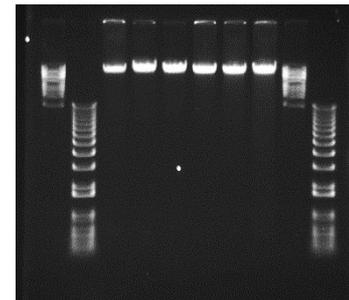
A260/280



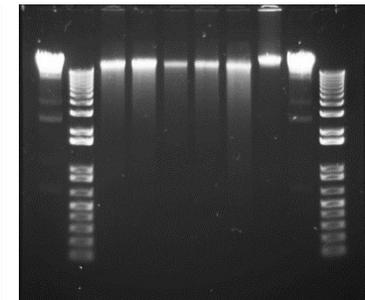
DNA tested in PCR based assay



Electrophoresis –uncut DNA



–DNA digestion



Point mutation validation

Example: RAS genetic alteration

ATCC® No.	Cell line name	Gene	AA Change	DNA Change	Zygoty	Coverage at Mutation Loci	% Zygoty
CRL-2177™	SW 1271	NRAS	p.Q61R	c.182A>G	Homozygous	26732	G = 99.8%
CRL-2273™	CHP-212	NRAS	p.Q61K	c.181C>A	Heterozygous	49859	C = 50.7, A = 49.1
CRL-7585™	Hs 852.T	NRAS	p.G12V	c.35G>T	Heterozygous	66411	G = 38.0, T = 61.8
CRL-9068™	NCI-H929	NRAS	p.G13D	c.38G>A	Heterozygous	21896	A = 53.9, G = 45.9
TIB-202™	THP-1	NRAS	p.G12D	c.35G>A	Heterozygous	60288	A = 70.1, G = 29.9
CRL-2547™	Panc 10.05	KRAS	p.G12D	c.35G>A	Heterozygous	42708	G = 52.7, A = 47.3
CRL-2549™	Panc 03.27	KRAS	p.G12V	c.35G>T	Heterozygous	58913	G = 47.0, T = 52.9
HTB-174™	NCI-H441	KRAS	p.G12V	c.35G>T	Heterozygous	87521	G = 52.8, T = 47.1
CL-187™	LS 180	KRAS	p.G12D	c.35G>A	Heterozygous	91234	G = 51.3, A = 48.6
CCL-225™	HCT-15	KRAS	p.G13D	c.38G>A	Heterozygous	49764	G = 52.1, A = 47.8

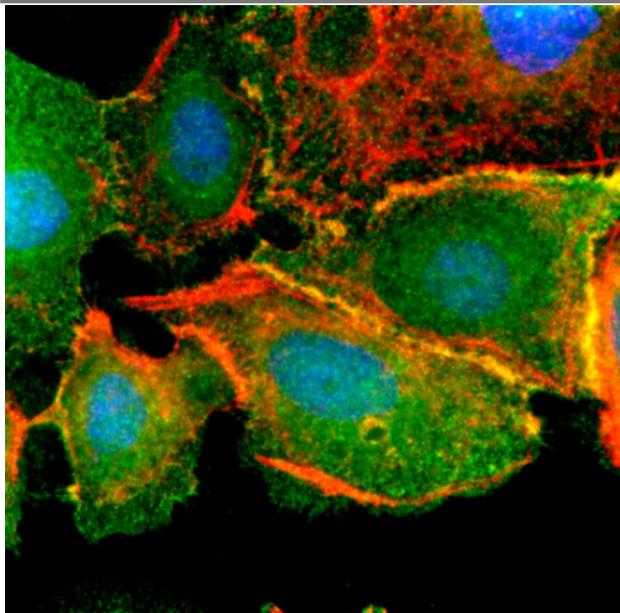
Point mutation validation

Example: EGFR genetic alteration

ATCC [®] No.	Cell line name	Gene	EGFR copy number variation	Measured CNV of EGFR	ERBB2 copy number variation	Measured CNV of ERBB2	Tumor source
CRL-2868™	HCC827	EGFR	Amplification	63.01	–	–	Lung
HTB-132™	MDA-MB-468	EGFR	Amplification	25.02	–	–	Breast
HTB-19™	BT-20	EGFR	Amplification	15.73	–	–	Breast
HTB-178™	NCI-H596	EGFR	Amplification	0.06	–	–	Lung
HTB-177™	NCI-H460	EGFR	–	–	–	–	Lung
CRL-5928™	NCI-H2170	ERBB2	–	–	Amplification	128.89	Lung
HTB-20™	BT-474	ERBB2	–	–	Amplification	29.70	Breast
HTB-27™	MDA-MB-361	ERBB2	–	–	Amplification	16.85	Breast

EGFR and HER2 are currently used as predictive markers of kinase inhibitor response in non-small cell lung cancer (NSCLC) and breast cancer therapy

ATCC quantitated human gDNA



- Purified from authenticated ATCC cell lines
- Contain oncology biomarkers
- Quantified gene mutation allelic frequencies, gene copy numbers
- Quantified by validated methods for each product lot
- Supplied with certificates providing measurement results with associated uncertainties for each product lot

Quantification of genetic variations in gDNAs

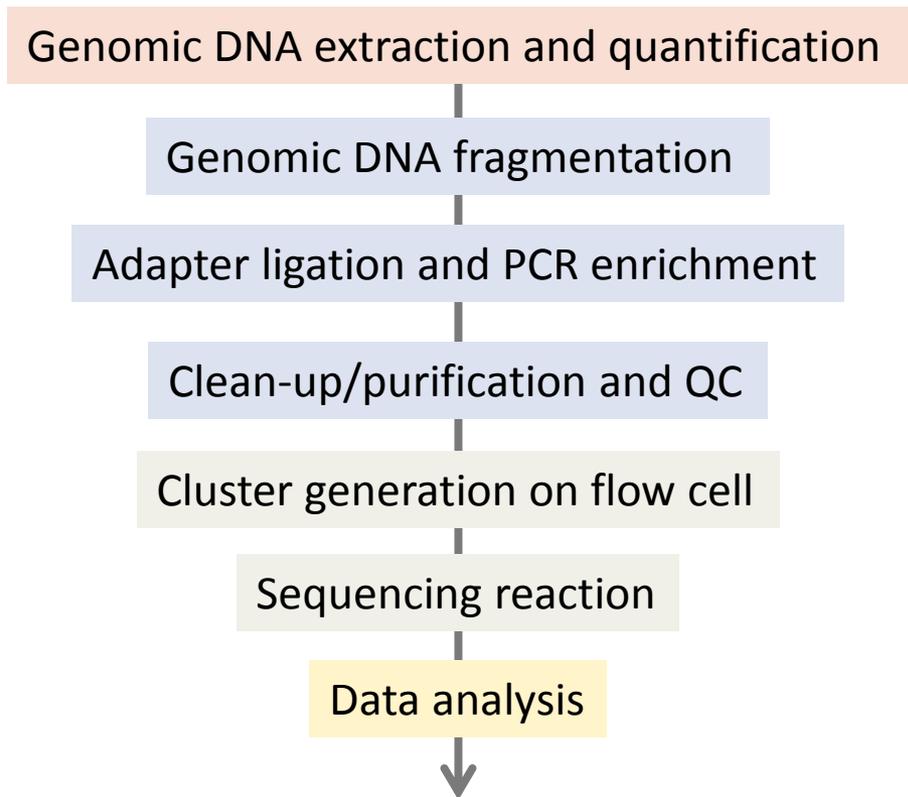
Coming soon: Quantitative gDNAs

ATCC® No.	Gene	Mutation
HTB-131DQ™	PIK3CA	H1074R
CRL-5908DQ™	EGFR	T790M; L858R
CRL-2868DQ™	EGFR	ELREA746del
CCL-231DQ™	EGFR	G719S
CRL-7898DQ™	BRAF	V600E
CCL-225DQ™	KRAS	G13D
CL-187DQ™	KRAS	G12D
CRL-2177DQ™	NRAS	Q61R
CCL-227DQ™	KRAS	G12V
	TP53	R273H
HTB-30DQ™	TP53	R175H
CRL-2158DQ™	TP53	R245S
CRL-1648DQ™	TP53	R248Q
HTB-122DQ™	TP53	R249S
HTB-111DQ™	pTEN	R130fs

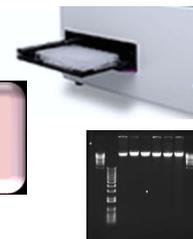
Use NGS to quantify mutations

Next-generation sequencing

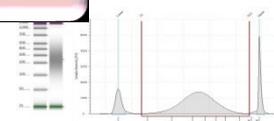
Workflow overview



High quality purified genomic DNA



Bioanalyser to ensure good quality library prep



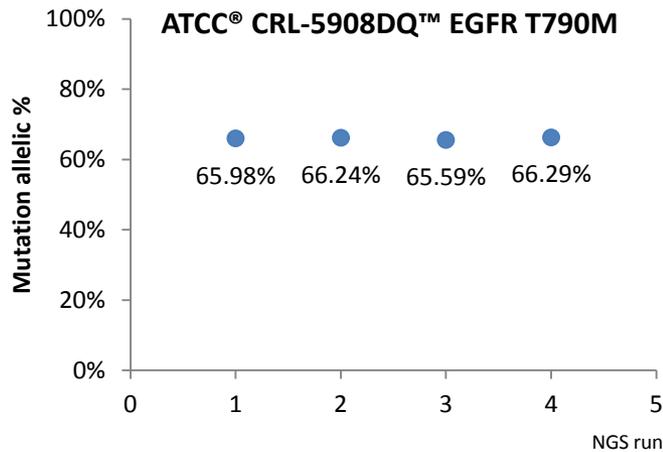
Deep sequencing to ensure the accuracy



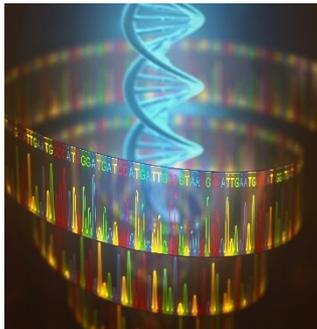
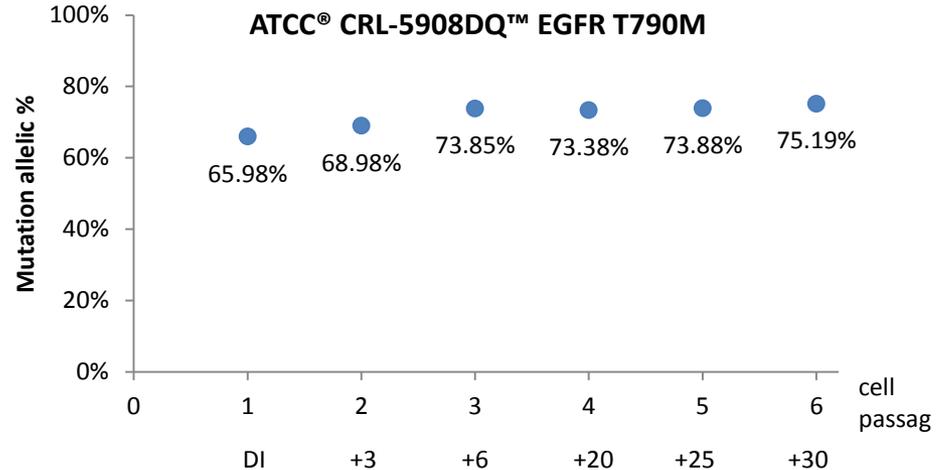
EGFR mutation quantification

Example of how cell passage number can affect mutation allelic frequency in some lines

Reproducibility in independent runs



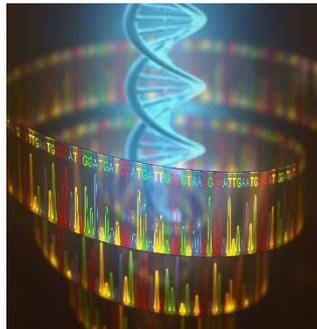
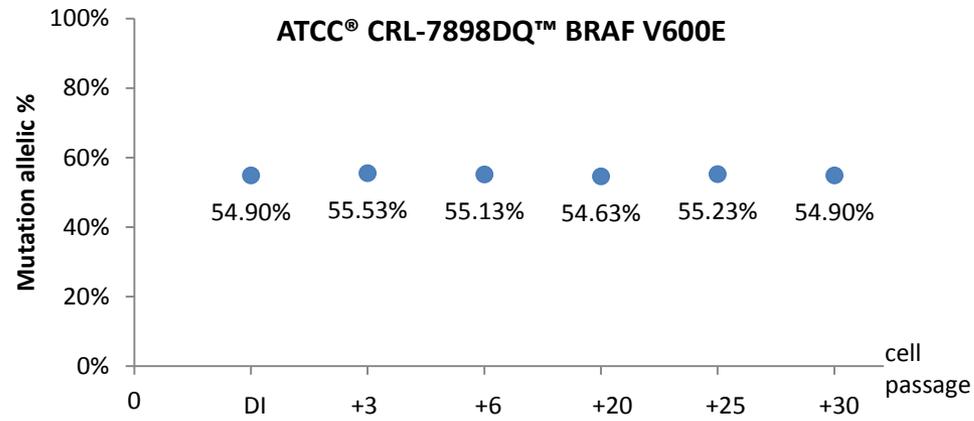
Mutation allelic frequency shifting



NGS
coverage >50,000X

NGS is performed on each lot to ensure biomarker quantification - your trusted control materials

BRAF mutation quantification



NGS
coverage >50,000X

The impact of passage number is cell line dependent, and some lines are relatively stable

Gene copy number quantification

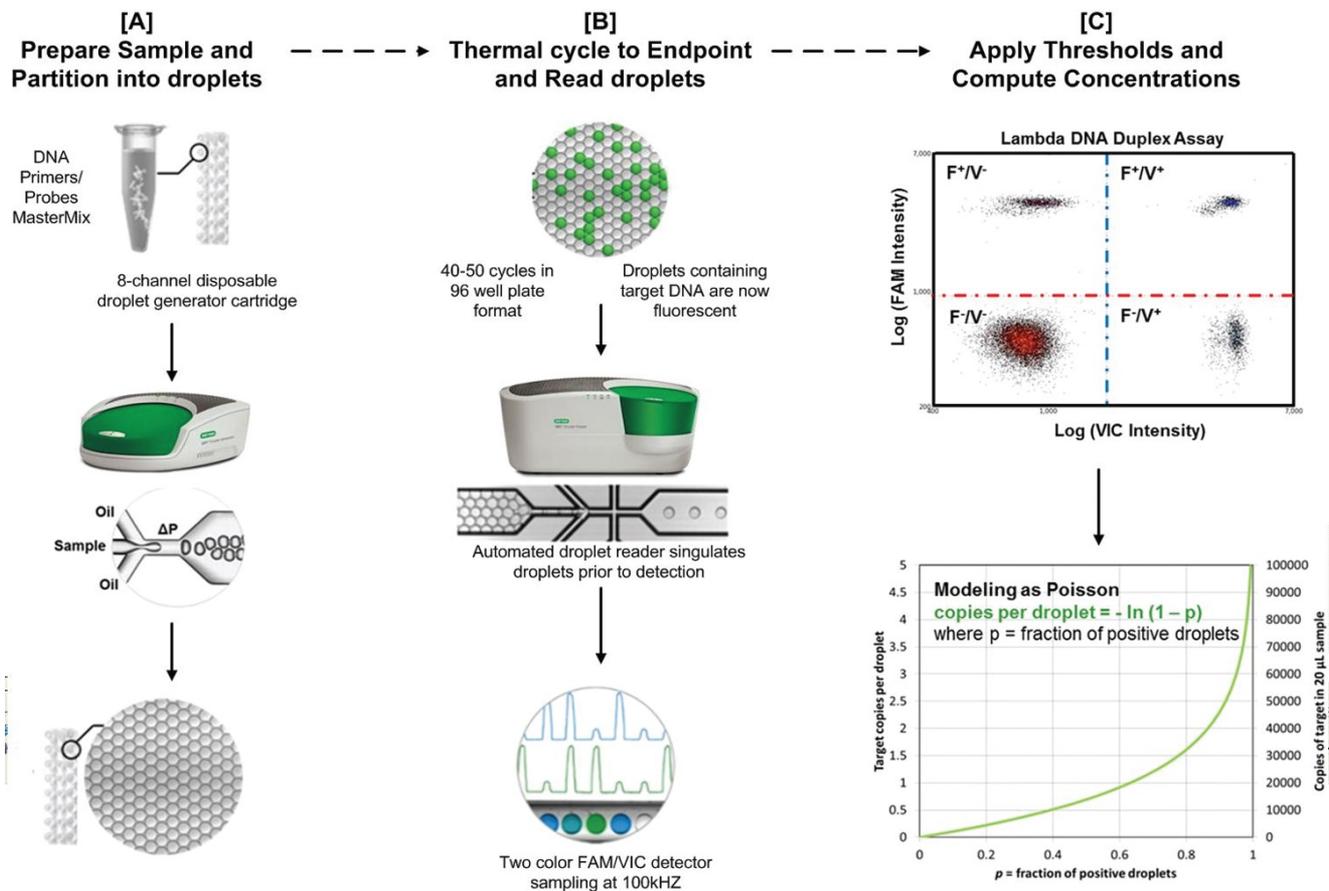
Gene amplifications have emerged as therapeutic targets and diagnostic biomarkers

ATCC® No.	Gene	Gene copy number variation	Parental cell line
CRL-2868DQ™	EGFR	amplification	HCC827
CRL-5928DQ™	ERBB2	amplification	NCI-H2170
CRL-5973DQ™	MET	amplification	SNU-5
CRL-5974DQ™	MYC	amplification	SNU-16

Use QX200 ddPCR™ (Bio-Rad) to perform quantification

- Absolute copy number of target gene
- Absolute copy number of housekeeping gene EIF2C1
- Relative copy number/ copy number variation

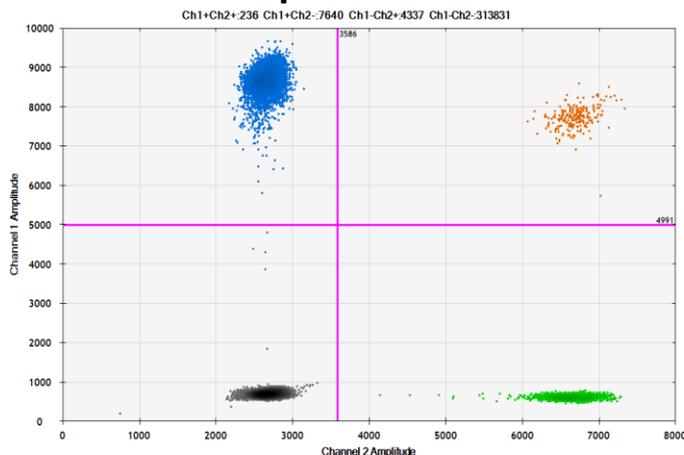
Overview of ddPCR™ workflow



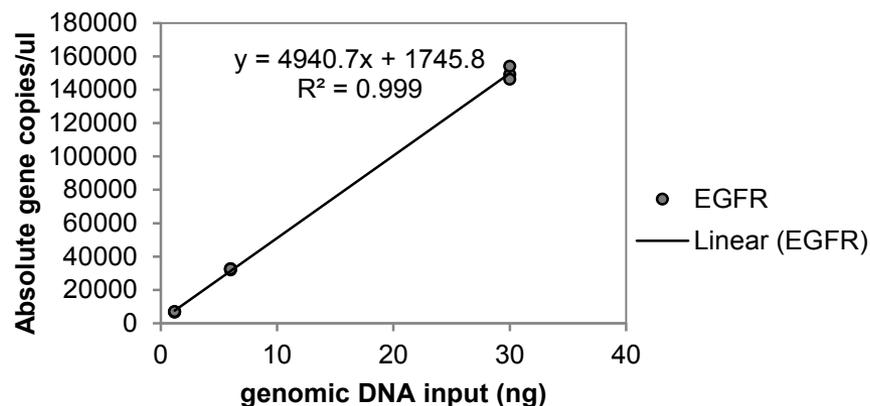
Anal Chem 84(2): 1003–1011, 2012.

Quantify EGFR amplification

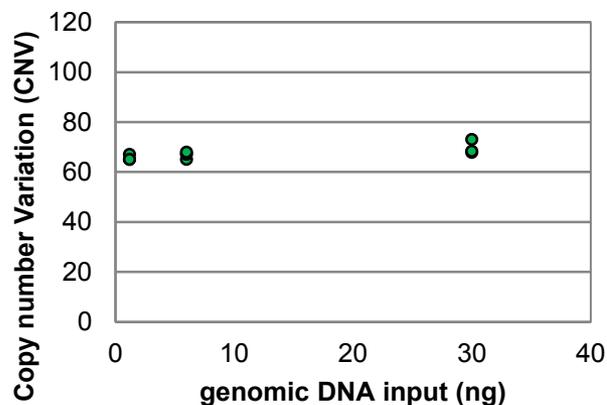
Multiplex ddPCR™



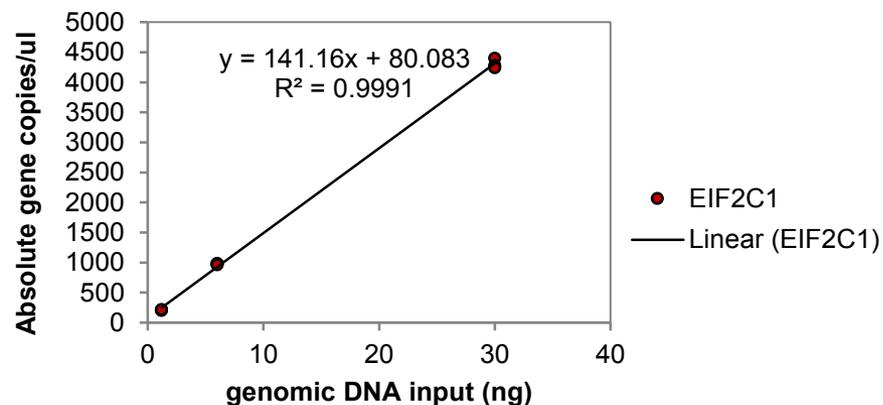
EGFR



CNV



EIF2C1

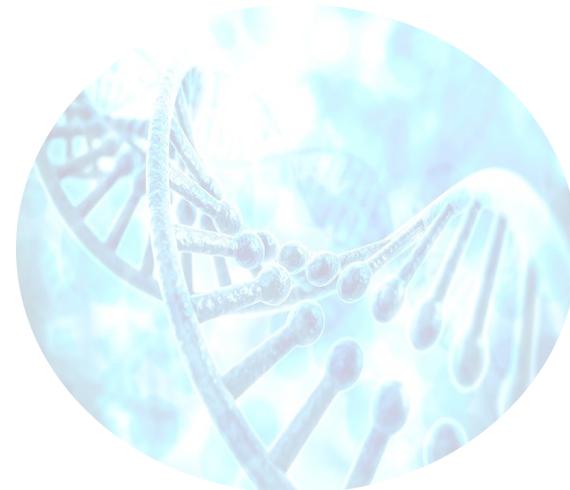


ddPCR™ is performed on each lot to ensure gene copy number quantification - your trusted control materials

Batch specific test results for each production lot

Example: ATCC® CRL-2868DQ™ quantitated human gDNA

- **Lot:** 63788713
- **Gene:** EGFR
- **Mutation:** ELREA746del
- **Mutation percentage:** 100%
- **EGFR absolute gene copy number:** 8217.83 copies/ng
- **EGFR relative gene copy number:** 65.83 copies



CoA report result – NGS (Coverage > 10,000X)

NGS result uncertainty is equal or smaller than $\pm 5\%$. The reported uncertainty represents uncertainty expressed at approximately the 99% confidence level using a coverage factor of $k=3$.

CoA report result – ddPCR™ (Average of nine data points)

ddPCR™ uncertainty is equal or smaller than $\pm 25\%$. The reported uncertainty represents uncertainty expressed at approximately the 99% confidence level using a coverage factor of $k=3$.

Conclusion

- Next-generation sequencing and other molecular tests have been used in clinical diagnostics, which is facilitating advances in disease prediction and therapeutic decision making
- Globally accepted, well-established reference materials are needed to ensure the reliability and reproducibility of diagnostic test results
- Quantitative gDNAs with known mutation allelic frequency and gene copy number provide a reliable and sustainable alternative to variable patient tissue derived controls in oncology molecular diagnostic assays



Acknowledgements

- Brian Chase
- John Foulke
- Michael Geimer
- Michael Jackson
- Stephen King
- Karin Kingdig
- Maria Mayda
- Lysa-Anne Volpe
- Melisa Wilson



Disclaimers

© 2016 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise. PicoGreen® and RiboGreen® are registered trademarks of Invitrogen. Bio-Rad®, Droplet Digital™, and QX200™ are trademarks of Bio-Rad Laboratories, Inc. RNAsable® is a registered trademark of Biomātrica, Inc. Precision Medicine Initiative is a registered trademark of the US Department of Health and Human Services.

ATCC nucleic acids are for research use only



Thank you for joining today!

Register for more ATCC “*Excellence in Research*” webinars, or watch recorded webinars, at www.atcc.org/webinars.

- **September 29, 2016**
12:00 PM ET
Keeping Cells Happy – Topics in Cell Health Maintenance and Viability
Steve Budd, M.S., *Product Line Business Specialist*, ATCC
- **November 3, 2016**
12:00 PM ET
The Biology of Anaerobic Bacteria and Predominant Propagation Practices
Allison Faust, *Senior Biologist*, ATCC
Nancy Krueger, *Senior Biologist*, ATCC

