

Simplifying assay development with molecular standards: Remove culturing from the equation

Kyle Young, B.S., M.B.A. Product Line Business Specialist, ATCC

Credible Leads to Incredible™



Agenda

- \checkmark What ATCC molecular standards provide
- ✓ Development of ATCC's synthetic molecular standards
- $\checkmark~$ The ATCC molecular standards portfolio
 - ✓ Including upcoming new products
- ✓ Validation data





About ATCC

- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's premier biological materials resource and standards development organization
 - 5,000 cell lines
 - 80,000 microorganisms
 - Genomic & synthetic nucleic acids
 - -Media/reagents

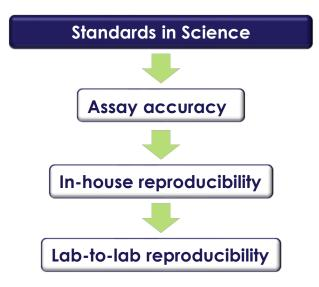
- ATCC collaborates with and supports the scientific community with industrystandard biological products and innovative solutions
- Growing portfolio of products and services
- Sales and distribution in 150 countries, 15 international distributors
- Talented team of 450+ employees, over one-third with advanced degrees



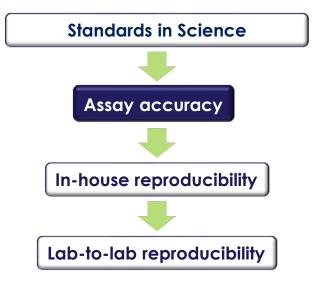


Molecular Standards

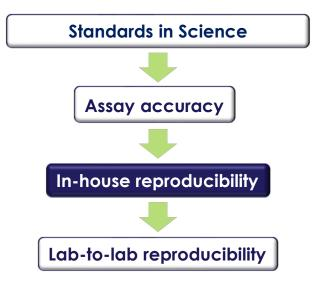




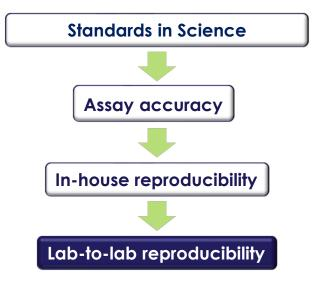














Project history

A culture collection provides good value for research.

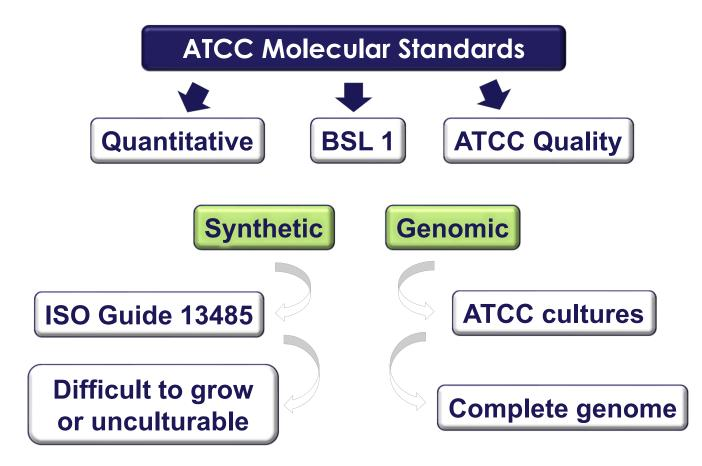
A standards organization with a culture collection provides even more!

In 2012, members of the ATCC R&D team realized there was more we could do for the scientific community, and began work on the first four molecular standards: three synthetic viruses and a quantitated bacterial DNA.

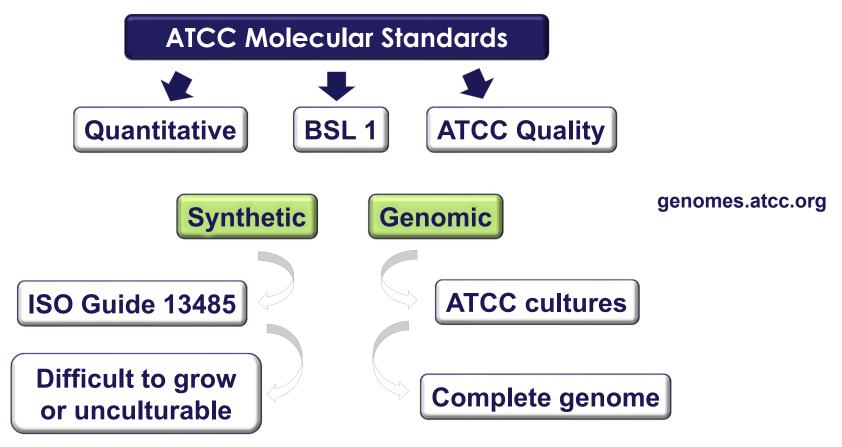
Those standards launched in May of 2013: VR-3198SD (West Nile virus RNA) VR-3199SD (Norovirus G1 RNA) VR-3299SD (Norovirus G2 RNA) 29212Q-FZ (Enterococcus faecalis)













Specifications

Synthetic Standards

Genomic Standards

Authentication	NGS to verify synthetic sequence	Authentication	Amplicon sequencing
Functionality & Identity	qPCR amplification, 3.32 cycles between Cq threshold	Integrity	High molecular weight DNA, by gel electrophoresis
Genome copy number by ddPCR™	1 x 10 ⁵ to 1 x 10 ⁶ genome copies/ μ L	Genome copy number by ddPCR™	1 x 10 ⁵ to 1 x 10 ⁶ genome copies/ μ L
Fill Volume	100 µL per vial	Fill Volume	100 µL per vial
Format	Frozen	Format	Frozen
Certifications	ISO 13485, ISO 9001, ISO 17025	Certifications	ISO 9001, ISO 17025



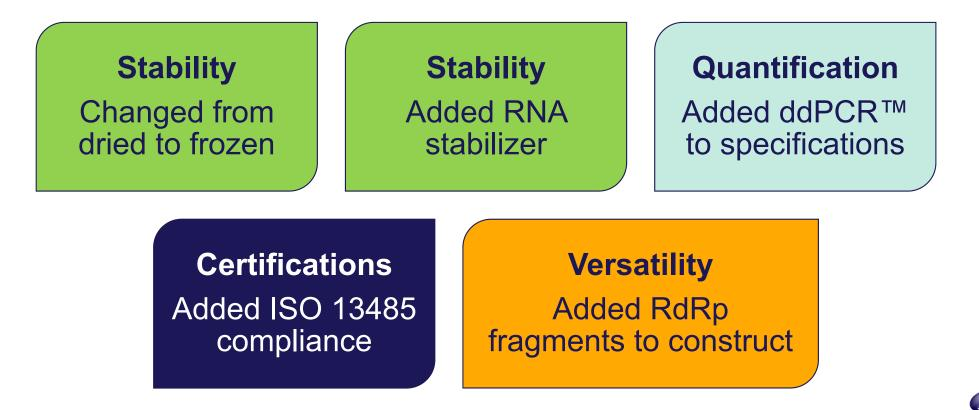


Synthetic Molecular Standard Design



Community response to the Norovirus standards

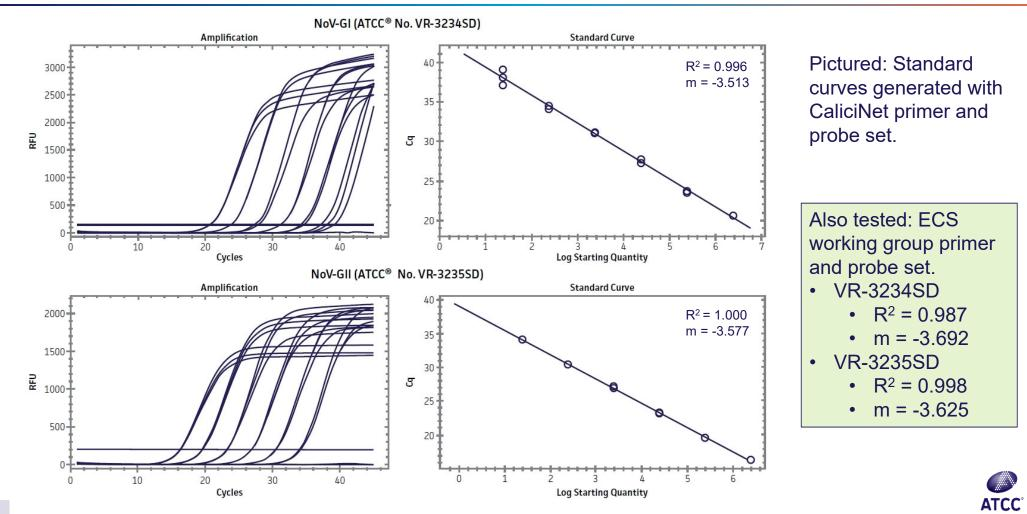
Interest in the standards was high, but feedback showed the synthetic constructs had room for improvement. ATCC modified the design and production processes, presenting the following changes at CVS in 2015.



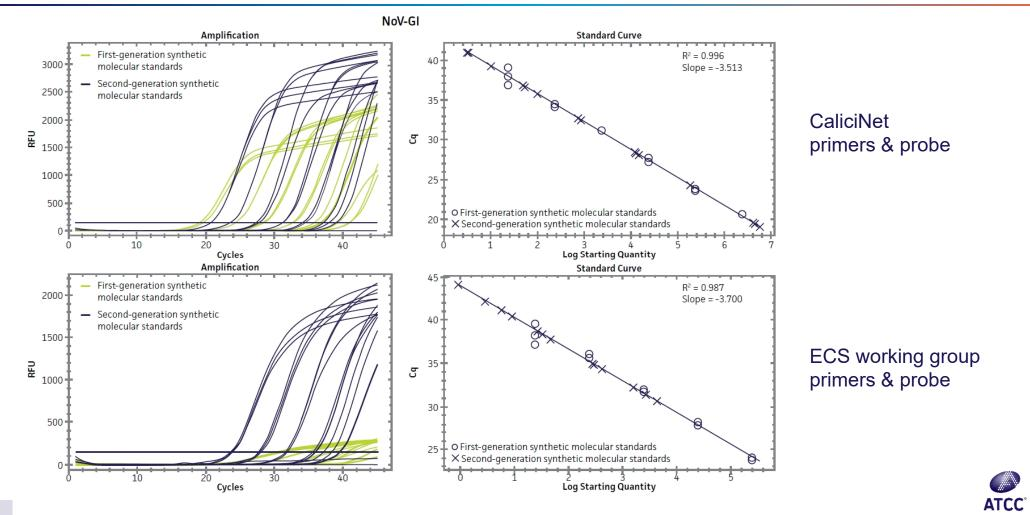
14

ATCC[°]

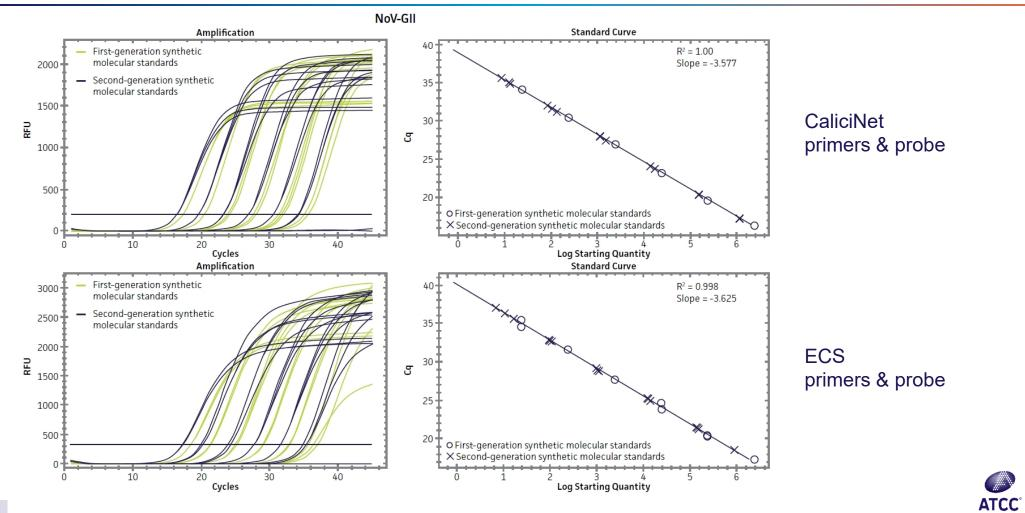
Validating the next generation of standards



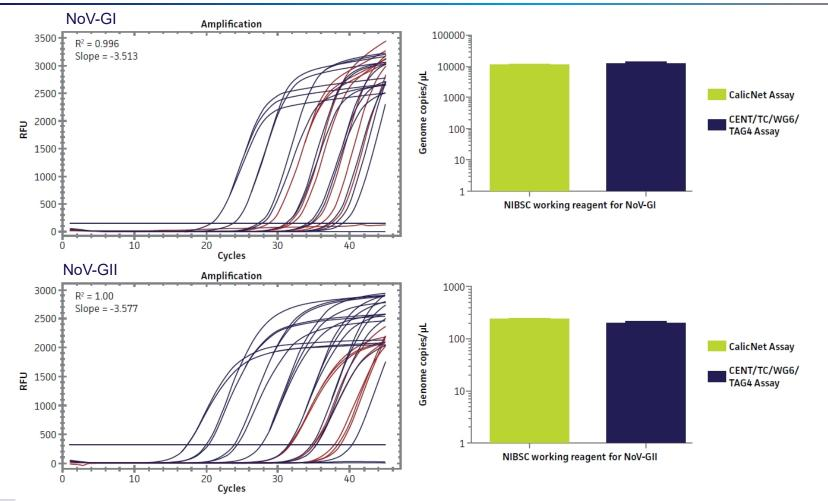
Old vs. new standards, part 1



Old vs. new standards, part 2

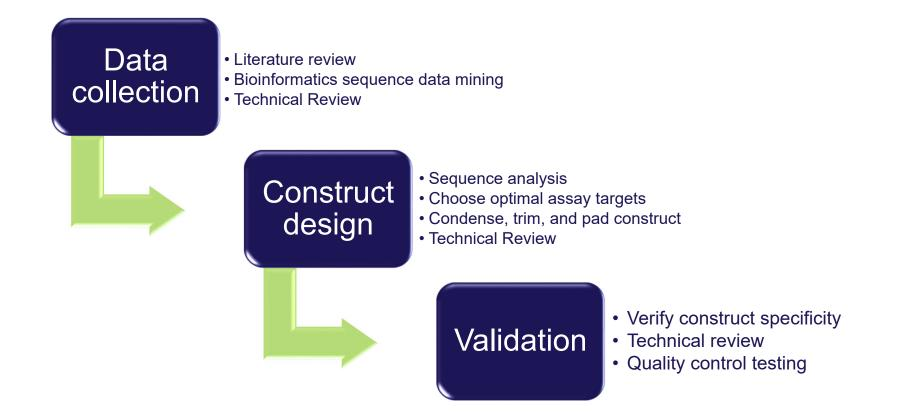


A test drive with NIBSC working reagents



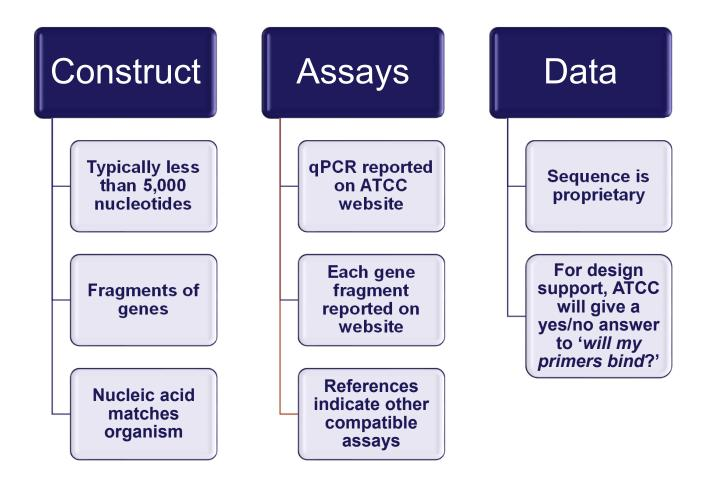


Design approach – synthetic standards





Design approach – synthetic standards







ATCC's Molecular Standards



Pathogen standards

- BK virus
- Hepatitis B virus
- Hepatitis C virus
- Epstein-Barr virus
- Human immunodeficiency virus 1
- Human T-cell leukemia virus
- Human cytomegalovirus
- Varicella-zoster virus
- Neisseria meningitides Plasmodium malariae
- Human parechovirus 3

Blood-borne disease

- Astrovirus
- Cyclospora cayetanensis
- Hepatitis A virus
- Hepatitis E virus
- Norovirus GI
- Norovirus GII
- Sapovirus
- Mycobacterium avium subsp.
- paratuberculosis
- Clostridiodes difficile
- · Salmonella enterica subsp. enterica serovar Typhimurium
- Cryptosporidium parvum
- •Human Enterovirus 71
- Rotavirus A
- Dientamoeba fragilis
 - Babesia canis
- Giardia lamblia

E. coli

- Murine norovirus
- Legionella pneumophila subsp.
- Pneumophila Human enterovirus 71 strain H Entamoeba histolvtica
- Gastro-Intestinal disease

- Human bocavirus
- Human coronavirus HKU1
- •Human coronavirus NL63
- •Human coronavirus 229E
- Middle East respiratory syndrome coronavirus
- •Human metapneumovirus
- Bordetella pertussis
- Mycobacterium africanum
- Mycobacterium bovis
- Mvcobacterium talmoniae
- Mvcobacterium microti
- Mycobacterium pinnipedii
- Mycobacterium tuberculosis
- Streptococcus pneumonia Human respiratory syncytial
- virus strain A2 Influenza B virus (BY) strain B/Wisconsin/1/2010BX-41A
- Influenza A virus (H3N2) strain A/Wisconsin/15/2009
- Influenza A virus (H1N1) strain A/PR/8/34
- Influenza B virus strain B/Florida/4/2006
- Betacoronavirus 1 strain OC43
- Human rhinovirus 77
- Bordetella pertussis
- Haemophilus influenza
- Streptococcus pneumoniae

Respiratory disease

Hepatitis B virus

- Human immunodeficiency virus 1
- •Human papillomavirus 16 •Human papillomavirus 18
- •Human papillomavirus 31
- •Human T-cell leukemia virus 2
- Treponema pallidum Chlamydia trachomatis
- serovar I Chlamydia trachomatis
- serovar II
- Chlamydia trachomatis serovar III
- •Human herpesvirus 1
- •Human herpesvirus 2
- Neisseria gonorrhoeae
- •Human Herpesvirus 8
- •Human herpesvirus 7 •Human herpesvirus 6
- Mycoplasma genitalium

Sexually transmitted



Chikungunya virus

- •Dengue virus types 1-4 •Eastern equine
 - encephalitis virus
 - Plasmodium malariae
 - •St. Louis encephalitis virus
 - West Nile virus
 - Yellow fever virus
 - Zika virus
 - Borrelia burgdorferi
 - Plasmodium falciparum •Yellow fever virus
 - •Rift Valley Fever virus

Vector-borne disease



- Staphylococcus
- epidermidis

www.atcc.org/MolecularStandards

- Streptococcus pyogenes
- Candida albicans
- Pseudomonas aeruginosa

Epidermal & Nosocomial disease





Standards from the microbiome



Skin

- Staphylococcus aureus
- Candida parapsilosis
- Staphylococcus
 epidermidis
- Candida tropicalis



Gut

- Streptococcus agalactiae
- Faecalibacterium
 prausnitzii
- Klebsiella aerogenes
- Bacteroides fragilis
- Enterococcus faecium
- Citrobacter freundii
- Escherichia coli
- Klebsiella oxytoca
- Klebsiella pneumoniae
- Enterococcus faecalis
- Candida albicans
- Helicobacter pylori



Water safety standards

Item Number	Description
29212Q-FZ	Quantitative DNA from Enterococcus faecalis
8739DQ	Quantitative DNA from Escherichia coli
25923DQ	Quantitative DNA from Staphylococcus aureus subsp. aureus
6538DQ	Quantitative DNA from Staphylococcus aureus subsp. aureus
9027DQ	Quantitative DNA from Pseudomonas aeruginosa
13048DQ	Quantitative DNA from Klebsiella aerogenes
10231DQ	Quantitative DNA from Candida albicans
4617DQ	Quantitative DNA from Bordetella bronchiseptica
25285DQ	Quantitative DNA from Bacteroides fragilis



Food safety standards

		Designation	ATCC Item Number
Big 6 E. Coli	Campylobacter	Norovirus GI	VR-3234SD
	jejuni	Norovirus GII	VR-3235SD
		Murine norovirus	VR-2355SD
		<i>E. coli</i> 0145	BAA-2192DQ
Colmonollo	Norovirus	<i>E. coli</i> O45:H2	BAA-2193DQ
Salmonella enterica Norovirus		<i>E. coli</i> O26:H11	BAA-2196DQ
		<i>E. coli</i> O103:H11	BAA-2215DQ
	<i>E. coli</i> O121:H19	BAA-2219DQ	
		<i>E. coli</i> O111	BAA-2440DQ
Sapovirus		C. jejuni	700819DQ
		S. enterica	700720DQ
		Sapovirus	VR-3237SD



Recently available

Item Number	Description	Comments
VR-3265SD	Quantitative Synthetic Human herpesvirus 7 DNA	Implicated in numerous health complications, and a transplant concern.
30001DQ	Quantitative Genomic DNA from <i>Trichomonas</i> vaginalis	Causative agent of trichomoniasis.
VR-1360DQ	Quantitative Genomic DNA from <i>Chlamydophila</i> pneumoniae	A causative agent of pneumonia.
27853DQ	Quantitative Genomic DNA from <i>Pseudomonas</i> aeruginosa	Quality control strain for numerous applications. Opportunistic pathogen.
15311DQ	Quantitative Genomic DNA from <i>Bordetella</i> parapertussis	Causative agent of whooping cough.



Soon available

Coming soon, in 2019

Item Number	Description
VR-1493DQ	Quantitative Genomic DNA from Human herpesvirus 1
VR-94DQ	Quantitative Genomic RNA from Human respiratory syncytial virus
22905DQ	Quantitative Genomic DNA from Proteus vulgaris
VR-1826DQ	Quantitative Genomic RNA from Enterovirus D68
MYA-2876DQ	Quantitative Genomic DNA from Candida albicans
VR-3268SD	Quantitative Synthetic Lassa virus RNA
VR-3269SD	Quantitative Synthetic Nipah virus RNA
VR-3266SD	Quantitative Synthetic Human immunodeficiency virus 2 RNA
VR-3274SD	Quantitative Synthetic West Nile virus
MYA-4941DQ	Quantitative Genomic DNA from Saccharomyces cerevisiae

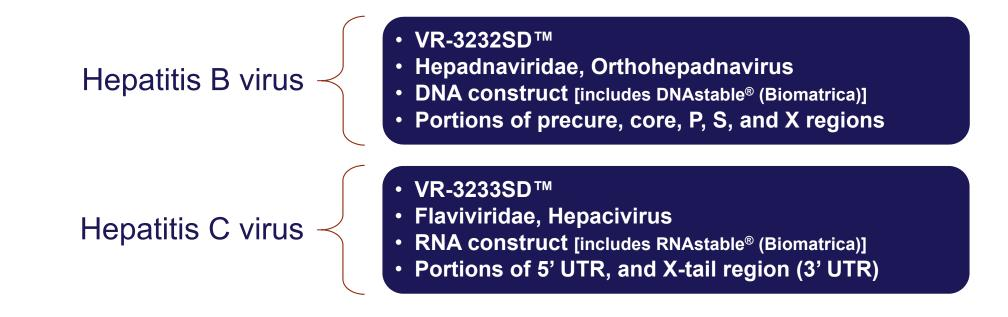




Validation of Molecular Standards



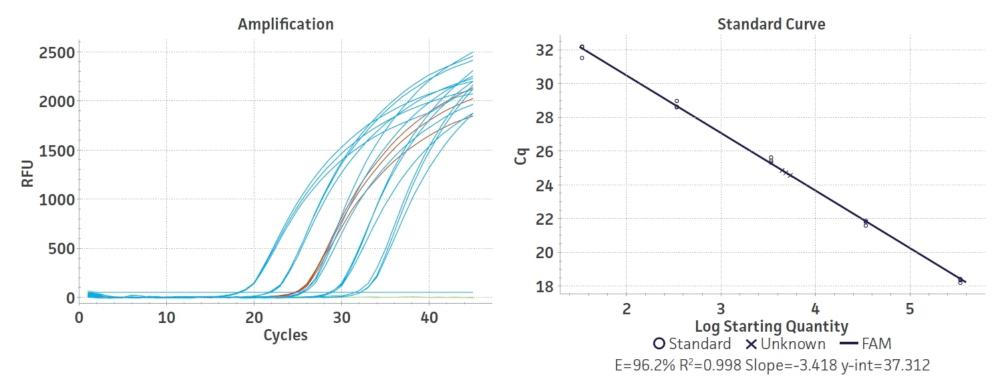
Validation of synthetic standards for hepatitis viruses



ATCC has also designed synthetic constructs for Hepatitis A virus (VR-3257SD[™]) and Hepatitis E virus (VR-3258SD[™]), and ATCC maintains a number of Hepatitis A viral stocks in its collection.



Hepatitis B virus

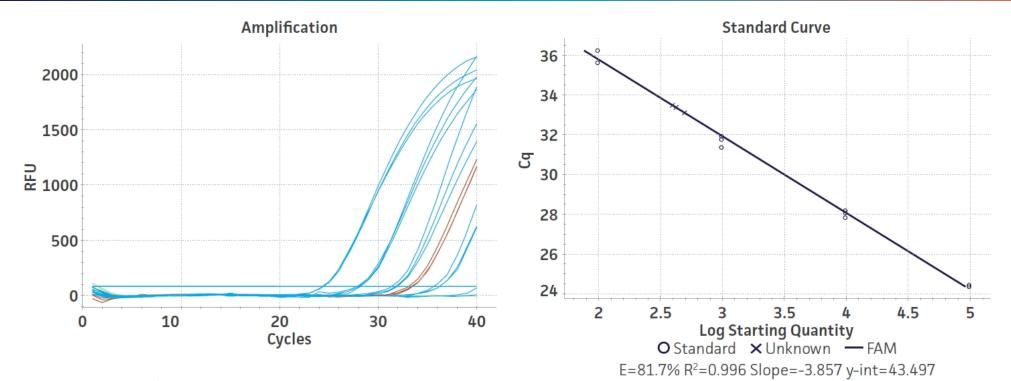


Blue = VR-3232SD[™] Red = NIBSC code 10/264 (3rd WHO international working reagent for HBV)

Sun S, et al. Development of a new duplex real-time polymerase chain reaction assay for hepatitis B viral DNA detection. Virol. J. 8: 227, 2011. PubMed: 21569595



Hepatitis C virus

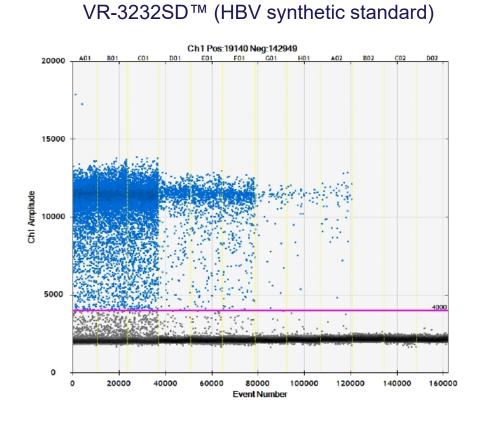


Blue = VR-3233SD[™] Red = NIBSC code 06/102 (4th WHO international standard for HCV)

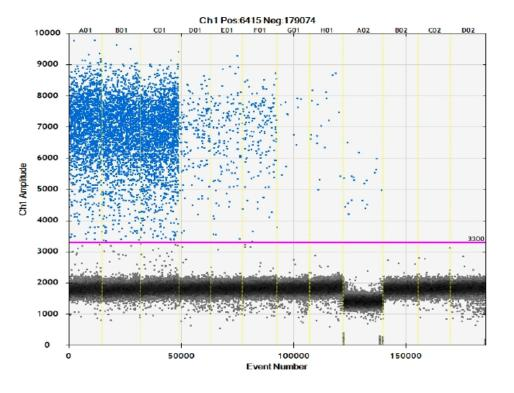
Lee SC, *et al.* Improved version 2.0 qualitative and quantitative AMPLICOR reverse transcription-PCR tests for hepatitis C virus RNA: calibration to international units, enhanced genotype reactivity, and performance characteristics. J. Clin. Microbiol. 38(11): 4171-4179, 2000. PubMed: 11060086

ATCC

Hepatitis viruses – ddPCR™

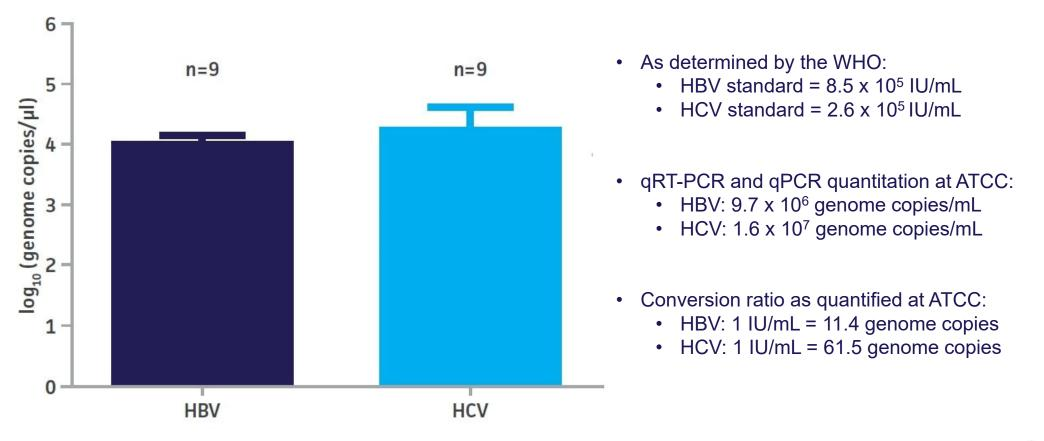


VR-3233SD™ (HCV synthetic standard)



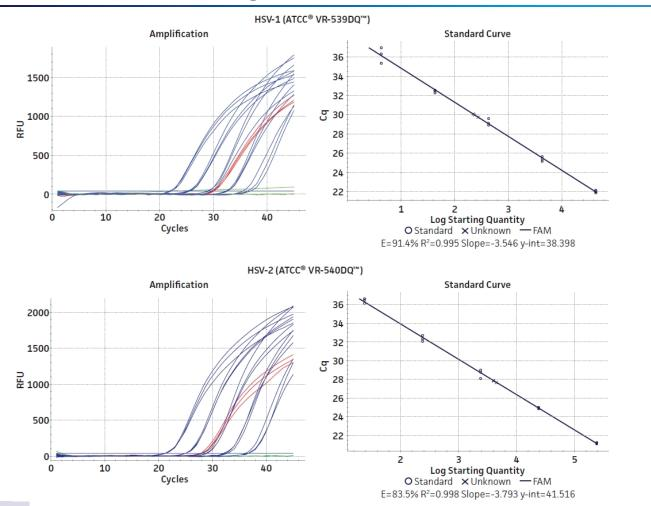
ATCC

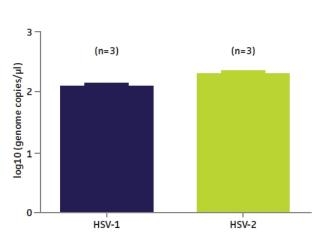
Quantitation of NIBSC Hepatitis standards





Human herpes viruses

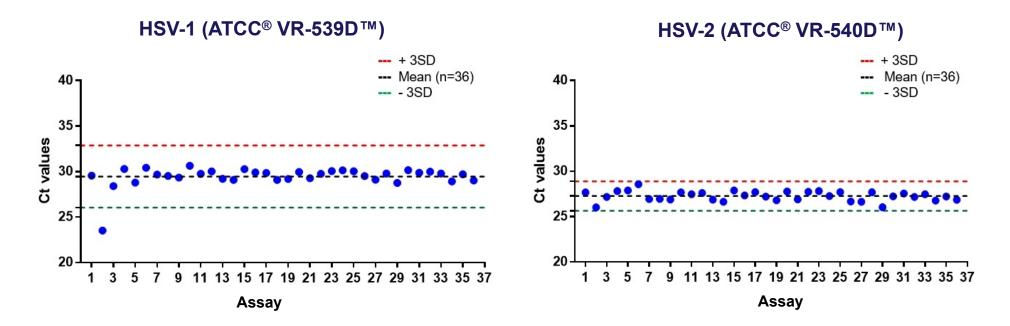




Ryncarz AJ, *et al.* Development of a high-throughput quantitative assay for detecting herpes simplex virus DNA in clinical samples. J Clin Microbiol 37(6): 1941-1947, 1999. Pubmed: 10325351



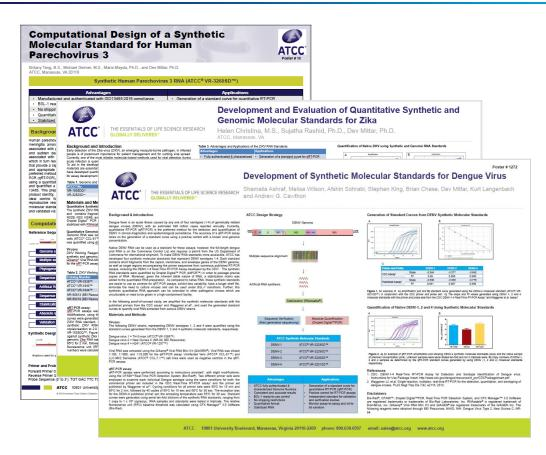
Human herpes viruses



	Average Ct	Standard Deviation	Coefficient of Variation
HSV-1 (ATCC [®] VR-539D™)	29.46	1.14	3.9%
HSV-2 (ATCC [®] VR-540D™)	27.27	0.54	2.0%



Other application data and posters



- Association of Molecular Pathology, 7-9 November, Baltimore
- American Society of Tropical Medicine and Hygiene, 20-24 November, National Harbor

https://www.atcc.org/Documents/Learning_Center/Research/Posters.aspx



Summary

- ATCC Molecular Standards are a consistent and reliable control for assay development and validation.
- Genomic standards eliminate the costs of growth and extraction.
- Synthetic standards provide controls for organisms that are difficult to culture.
- 159 standards currently in the portfolio.
 - 115 genomic standards
 - 44 synthetic standards
 - Standards for pathogens, microbiome, & food safety
 - Another 20 to 25 more to launch in 2019!





Thank you to the project team!

ATCC R&D, Technical Transfer, and Marketing Teams

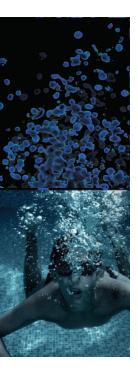
Cincinnati Children's Hospital, Department of Pathology, Donna Diorio

National Institute for Biological Standards and Control (NIBSC)

Stanford University Medical Center – Benjamin Pinsky







Questions?

Credible Leads to Incredible™



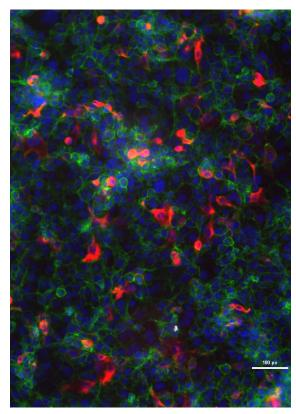
Cultivating collaboration to support global health

Go to www.atcc.org/molecularstandards for more information

Upcoming webinars:

- EMT/MET reporter cell lines: Elevating biological models of metastasis October 24, 12:00 ET
- On the edge of the bubble: Use of exosomes as reference materials in biomedical research October 31, 12:00 ET

www.atcc.org/webinars



ΔΤϹϹ

© 2019 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. ddPCR is a trademark of Bio-Rad Laboratories, Inc.