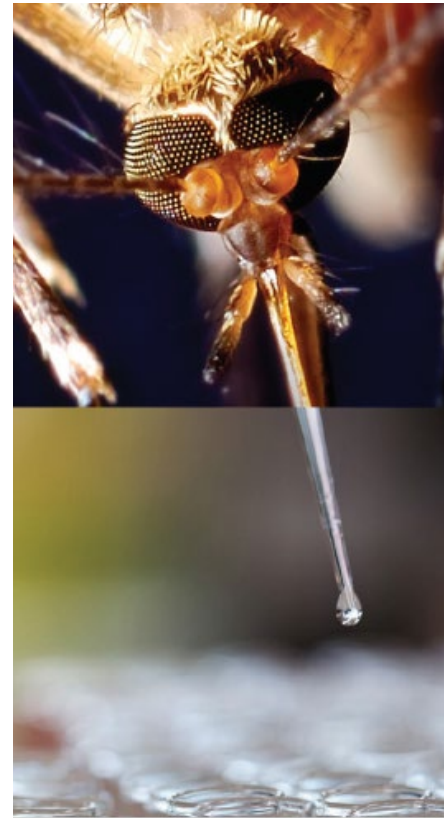
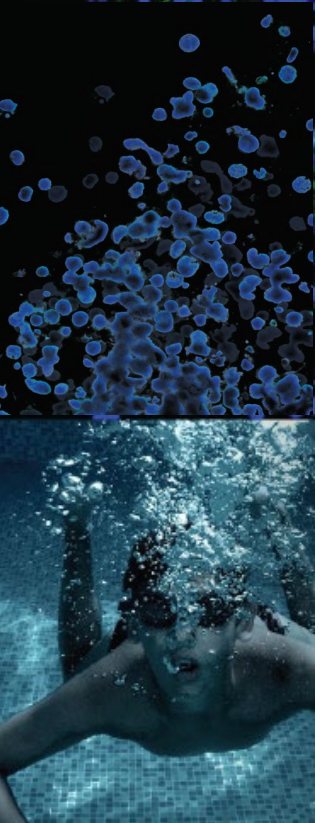




# Simplifying assay development with molecular standards: Remove culturing from the equation

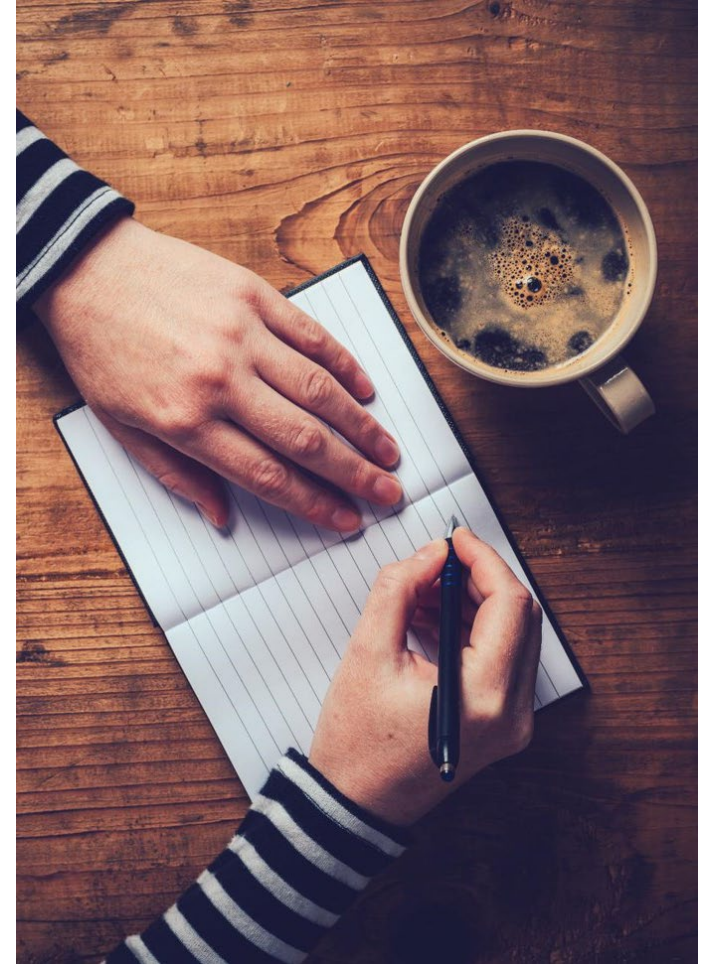
Kyle Young, BS, MBA  
Product Specialist, ATCC

Credible Leads to Incredible™



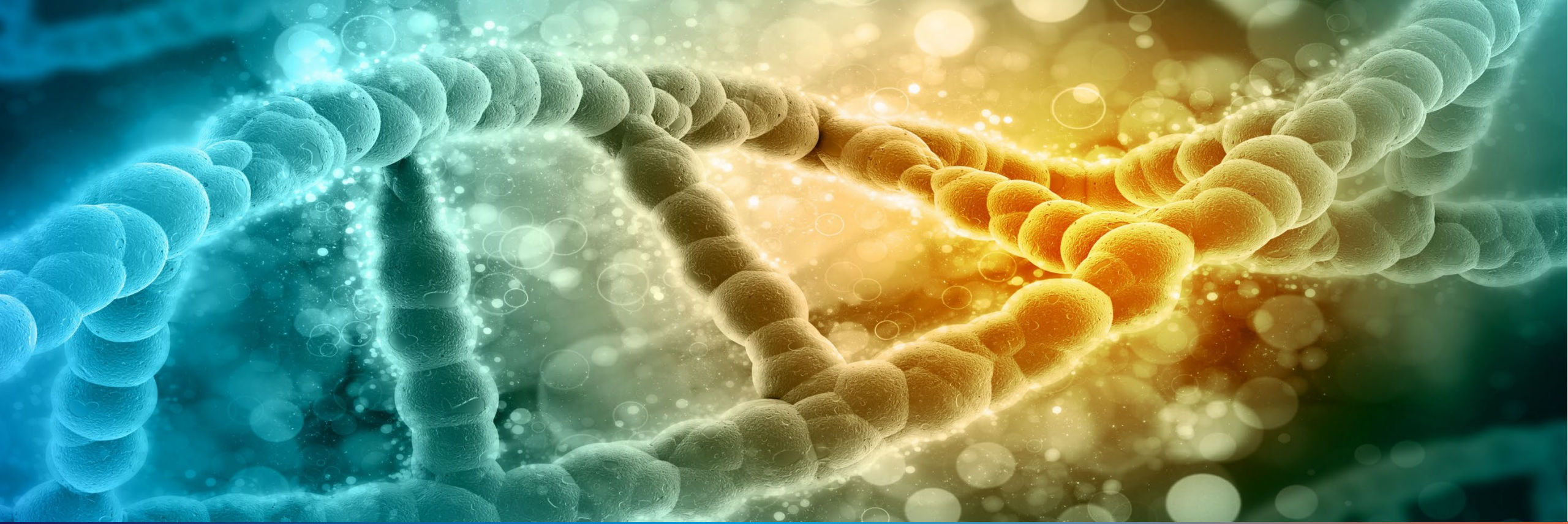
# Agenda

- ✓ How ATCC molecular standards accelerate assay development
- ✓ Development process for ATCC's synthetic molecular standards
- ✓ Validation data = materials you can trust



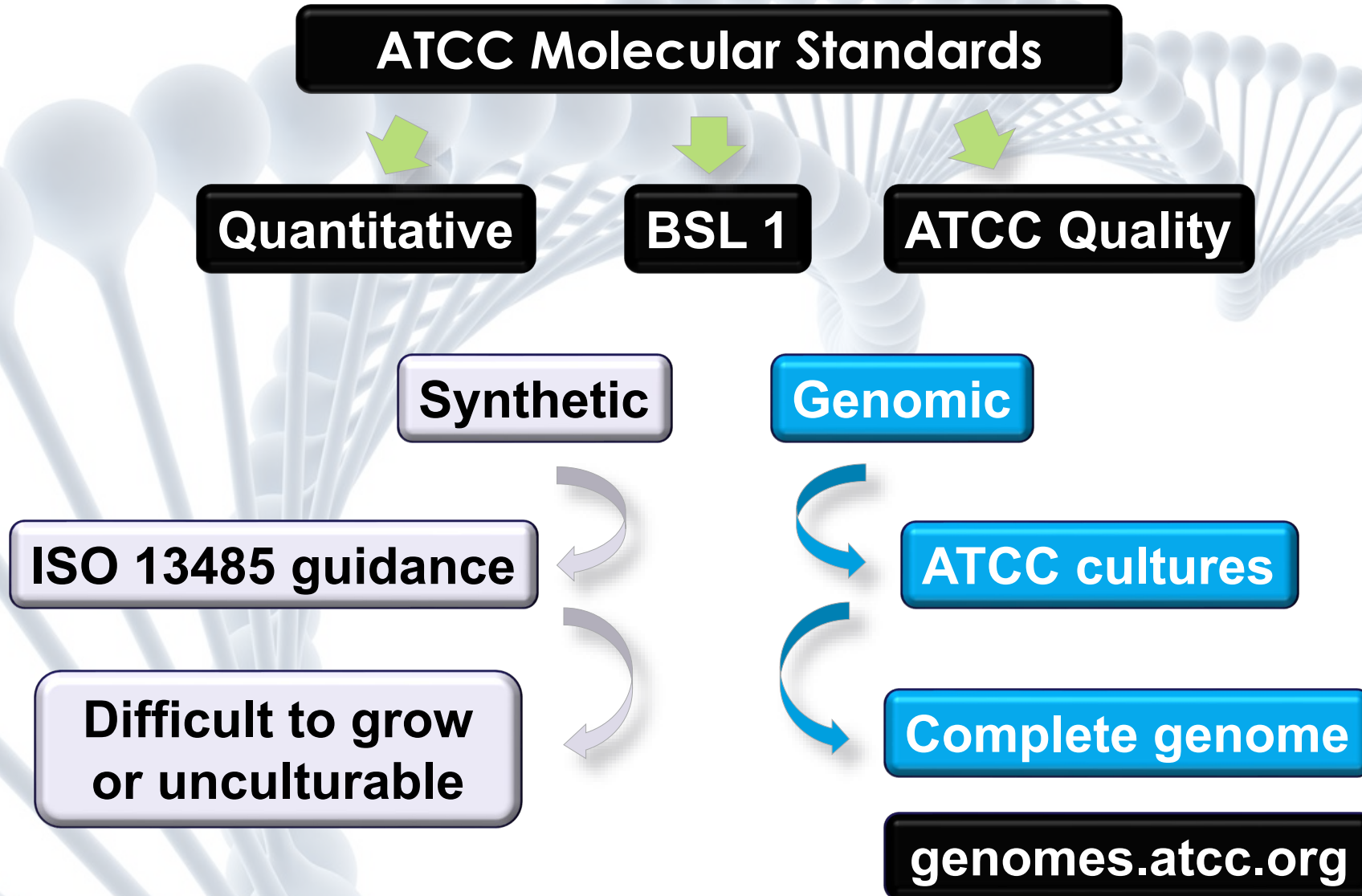
# About ATCC

- 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 industry-standard biological products and innovative solutions
- Growing portfolio of products and services
- Sales and distribution in 150 countries, 19 international distributors
- Talented team of 500+ employees, over one-third with advanced degrees



# Molecular Standards

# ATCC molecular standards



# Specifications

## Synthetic Standards

## Genomic Standards

<b>Authentication</b>	NGS to verify synthetic sequence
<b>Functionality &amp; Identity</b>	qPCR amplification, 3.32 cycles between Cq threshold
<b>Genome copy number by ddPCR™</b>	$1 \times 10^5$ to $1 \times 10^6$ construct copies/ $\mu\text{L}$
<b>Fill Volume</b>	100 $\mu\text{L}$ per vial
<b>Format</b>	Frozen

<b>Authentication</b>	Amplicon sequencing
<b>Integrity</b>	High molecular weight DNA by gel electrophoresis
<b>Genome copy number by ddPCR™</b>	$1 \times 10^5$ to $1 \times 10^6$ genome copies/ $\mu\text{L}$
<b>Fill Volume</b>	100 $\mu\text{L}$ per vial
<b>Format</b>	Frozen

# Assay development

## Molecular Diagnostics

Assay Sensitivity

Assay Specificity

Inclusivity

Limit of Detection

Exclusivity

# Assay Sensitivity - Inclusivity testing

Molecular Diagnostics

Assay Sensitivity

Assay Specificity

Inclusivity

Limit of Detection

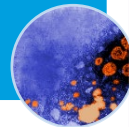
Exclusivity



# A menu for assay design

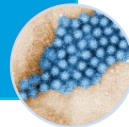
- BK virus
- Hepatitis B virus
- Hepatitis C virus
- Epstein-Barr virus
- Human immunodeficiency virus 1
- Human T-cell leukemia virus
- Human cytomegalovirus
- Varicella-zoster virus
- *Neisseria meningitidis*
- *Plasmodium malariae*
- *Human parechovirus 3*

## Blood-borne disease



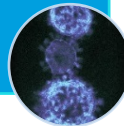
- Astrovirus
- *Cyclospora cayetanensis*
- Hepatitis A virus
- Hepatitis E virus
- Norovirus GI
- Norovirus GII
- Sapovirus
- *Mycobacterium avium* subsp. *paratuberculosis*
- *Clostridiodes difficile*
- *Salmonella enterica* subsp. *enterica* serovar Typhimurium
- *Cryptosporidium parvum*
- Human Enterovirus 71
- Rotavirus A
- *Dientamoeba fragilis*
- *Babesia canis*
- *Giardia lamblia*
- Murine norovirus
- *Legionella pneumophila* subsp. *pneumophila*
- Human enterovirus 71 strain H
- *Entamoeba histolytica*
- *E. coli*

## Gastro-Intestinal disease



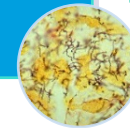
- SARS-CoV-2
- SARS-CoV
- MERS-CoV
- Human coronavirus OC43
- Human coronavirus HKU1
- Human coronavirus NL63
- Human coronavirus 229E
- Human metapneumovirus
- *Bordetella pertussis*
- *Mycobacterium bovis*
- *Mycobacterium talmoniae*
- *Mycobacterium microti*
- *Mycobacterium pinnipedii*
- *Mycobacterium tuberculosis*
- *Streptococcus pneumoniae*
- Human respiratory syncytial virus
- Influenza B virus (Victoria)
- Influenza B virus (Yamagata)
- Influenza A virus (H3N2)
- Influenza A virus (H1N1)
- Human bocavirus
- *Bordetella pertussis*
- *Haemophilus influenzae*
- Adenovirus
- Parainfluenza viruses
- Rhinoviruses
- *Chlamydia pneumoniae*
- *Legionella pneumophila*
- *Mycoplasma pneumoniae*

## Respiratory disease



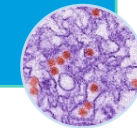
- *Neisseria gonorrhoeae*
- Human immunodeficiency virus 1
- Human papillomavirus 16
- Human papillomavirus 18
- Human papillomavirus 31
- Human T-cell leukemia virus 2
- *Treponema pallidum*
- *Chlamydia trachomatis* LGV I
- *Chlamydia trachomatis* LGV II
- *Chlamydia trachomatis* LGV III
- Human herpesvirus 1
- Human herpesvirus 2
- Hepatitis B virus
- Human herpesvirus 8
- Human herpesvirus 7
- Human herpesvirus 6
- *Mycoplasma genitalium*
- *Staphylococcus saprophyticus*
- *Hamophilus ducreyi*

## Sexually transmitted infections



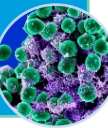
- Powassan virus
- Chikungunya virus
- Dengue virus types 1-4
- Eastern equine encephalitis virus
- *Plasmodium malariae*
- St. Louis encephalitis virus
- West Nile virus
- Yellow fever virus
- Zika virus
- *Borrelia burgdorferi*
- *Plasmodium falciparum*
- Yellow fever virus
- Rift Valley Fever virus

## Vector-borne disease



- *Staphylococcus aureus* subsp. *aureus*
- *Staphylococcus epidermidis*
- *Streptococcus pyogenes*
- *Candida albicans*
- *Pseudomonas aeruginosa*
- *Candida krusei*

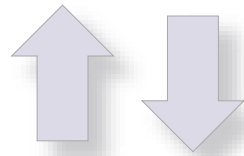
## Epidermal & Nosocomial disease



# Assay scope – strain selection for safety testing

## Water safety

- *Enterococcus faecalis*
- *Vibrio cholerae*
- *Cryptosporidium parvum*
- Rotavirus
- *Pseudomonas aeruginosa*
- *Escherichia coli* serotype O157:H7



## Food safety

- Norovirus
- Big Six *Escherichia coli*
- *Campylobacter jejuni*
- *Salmonella enterica*
- *Listeria monocytogenes*
- Sapovirus

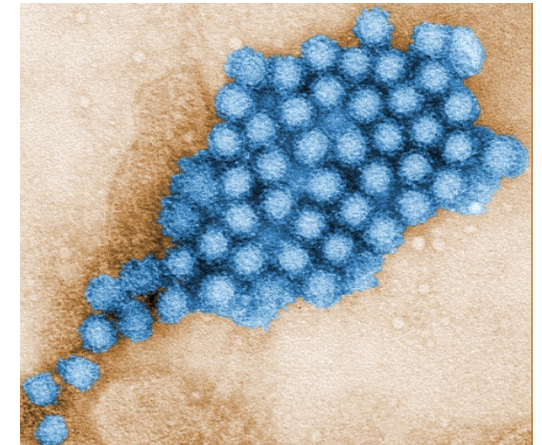


Photo credit: CDC, Dr. Charles D. Humphrey

# SARS-CoV-2 molecular standards



BSL 1

ATCC® No.	Description	Compatible assays
VR-3276SD™	Quantitative Synthetic SARS-CoV-2 RNA containing portions of ORF1ab, N, E, nsp12 (RdRp), and ORF1b-nsp14 genes	<ul style="list-style-type: none"> <li>•China CDC Primers and probes for detection 2019-nCoV (24 January 2020)</li> <li>•Diagnostic detection of Wuhan coronavirus 2019 by real-time RT-PCR – Charité, Berlin Germany (17 January 2020)</li> <li>•Detection of 2019 novel coronavirus (2019-nCoV) in suspected human cases by RT-PCR – Hong Kong University (23 January 2020)</li> <li>•PCR and sequencing protocol for 2019-nCoV - Department of Medical Sciences, Ministry of Public Health, Thailand (Updated 28 January 2020)</li> <li>•US CDC Real-Time RT-PCR Panel for Detection 2019-Novel Coronavirus (28 January 2020)</li> <li>•US CDC panel primer and probes– U.S. CDC, USA (28 January 2020)</li> </ul>
VR-3277SD™	Quantitative Synthetic SARS-CoV-2 RNA: containing a portion of Spike 5' end gene.	<ul style="list-style-type: none"> <li>•Detection of WN-Human1 sequence from clinical specimen. – National Institute of Infectious Diseases Japan (17 January 2020)</li> </ul>
VR-3278SD™	Quantitative Synthetic SARS-CoV-2 RNA: containing a portion of Spike 3' end gene.	<ul style="list-style-type: none"> <li>•PCR and sequencing protocols for 2019-nCoV- National Institute of Infectious Diseases Japan (24 January 2020)</li> </ul>
VR-3279SD™	Quantitative Synthetic SARS-CoV-2 RNA containing portions of the nsp9 and nsp12 (RdRp) genes	<ul style="list-style-type: none"> <li>•RT-PCR assays for the detection of SARS-CoV-2 with RdRp - Institut Pasteur, Paris (2 March 2020)</li> <li>•Diagnostic detection of Wuhan coronavirus 2019 by real-time RT-PCR – Charité, Berlin Germany (17 January 2020)</li> </ul>



BSL 2

ATCC® No.	Genomic RNA from isolate
VR-1986D™	USA-WA1/2020 lineage A
VR-1991D™	Hong Kong/VM20001061/2020 lineage A
VR-1992D™	Italy/INMI1 lineage B
VR-1994D™	Germany/BavPat1/2020 lineage B (D614G mutation)
VR-3326D™	USA/CA_CDC_5574/2020 lineage B.1.1.7
VR-3327D™ *	USA/MD-HP01542/2021 lineage B.1.351
VR-3338D™ *	Japan/TY7-503/2021 lineage P.1

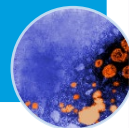
\* In development

# A menu for assay design

Inclusivity candidate

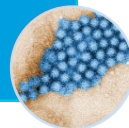
- BK virus
- Hepatitis B virus
- Hepatitis C virus
- Epstein-Barr virus
- Human immunodeficiency virus 1
- Human T-cell leukemia virus
- Human cytomegalovirus
- Varicella-zoster virus
- *Neisseria meningitidis*
- *Plasmodium malariae*
- *Human parechovirus 3*

## Blood-borne disease



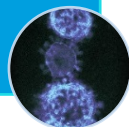
- Astrovirus
- *Cyclospora cayetanensis*
- Hepatitis A virus
- Hepatitis E virus
- Norovirus GI
- Norovirus GII
- Sapovirus
- *Mycobacterium avium* subsp. *paratuberculosis*
- *Clostridiodes difficile*
- *Salmonella enterica* subsp. *enterica* serovar Typhimurium
- *Cryptosporidium parvum*
- Human Enterovirus 71
- Rotavirus A
- *Dientamoeba fragilis*
- *Babesia canis*
- *Giardia lamblia*
- Murine norovirus
- *Legionella pneumophila* subsp. *pneumophila*
- Human enterovirus 71 strain H
- *Entamoeba histolytica*
- *E. coli*

## Gastro-Intestinal disease



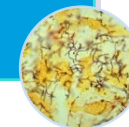
- SARS-CoV-2
- SARS-CoV
- MERS-CoV
- Human coronavirus OC43
- Human coronavirus HKU1
- Human coronavirus NL63
- Human coronavirus 229E
- Human metapneumovirus
- *Bordetella pertussis*
- *Mycobacterium bovis*
- *Mycobacterium talmoniae*
- *Mycobacterium microti*
- *Mycobacterium pinnipedii*
- *Mycobacterium tuberculosis*
- *Streptococcus pneumoniae*
- Human respiratory syncytial virus
- Influenza B virus (Victoria)
- Influenza B virus (Yamagata)
- Influenza A virus (H3N2)
- Influenza A virus (H1N1)
- Human bocavirus
- *Bordetella pertussis*
- *Haemophilus influenzae*
- Adenovirus
- Parainfluenza viruses
- Rhinoviruses
- *Chlamydia pneumoniae*
- *Legionella pneumophila*
- *Mycoplasma pneumoniae*

## Respiratory disease



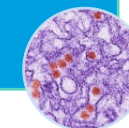
- *Neisseria gonorrhoeae*
- Human immunodeficiency virus 1
- Human papillomavirus 16
- Human papillomavirus 18
- Human papillomavirus 31
- Human T-cell leukemia virus 2
- *Treponema pallidum*
- *Chlamydia trachomatis* LGV I
- *Chlamydia trachomatis* LGV II
- *Chlamydia trachomatis* LGV III
- Human herpesvirus 1
- Human herpesvirus 2
- Hepatitis B virus
- Human herpesvirus 8
- Human herpesvirus 7
- Human herpesvirus 6
- *Mycoplasma genitalium*
- *Staphylococcus saprophyticus*
- *Hamophilus ducreyi*

## Sexually transmitted infections



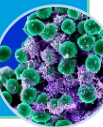
- Powassan virus
- Chikungunya virus
- Dengue virus types 1-4
- Eastern equine encephalitis virus
- *Plasmodium malariae*
- St. Louis encephalitis virus
- West Nile virus
- Yellow fever virus
- Zika virus
- *Borrelia burgdorferi*
- *Plasmodium falciparum*
- Yellow fever virus
- Rift Valley Fever virus

## Vector-borne disease



- *Staphylococcus aureus* subsp. *aureus*
- *Staphylococcus epidermidis*
- *Streptococcus pyogenes*
- *Candida albicans*
- *Pseudomonas aeruginosa*
- *Candida krusei*

## Epidermal & Nosocomial disease



# Assay sensitivity - limit of detection

Molecular Diagnostics

Assay Sensitivity

Assay Specificity

Inclusivity

Limit of Detection

Exclusivity

# Assay specificity - exclusivity testing

Molecular Diagnostics

```
graph TD; MD[Molecular Diagnostics] --> AS[Assay Sensitivity]; MD --> ASpec[Assay Specificity]; AS --> Inc[Inclusivity]; AS --> LOD[Limit of Detection]; ASpec --> Ex[Exclusivity];
```

The diagram illustrates the relationship between various assay performance metrics. It starts with 'Molecular Diagnostics' at the top, which branches into 'Assay Sensitivity' and 'Assay Specificity'. 'Assay Sensitivity' further branches into 'Inclusivity' and 'Limit of Detection'. 'Assay Specificity' leads to 'Exclusivity'. The 'Assay Specificity' and 'Exclusivity' boxes are highlighted in blue, while the others are light purple.

Assay Sensitivity

Assay Specificity

Inclusivity

Limit of Detection

Exclusivity

# Resources for SARS-CoV-2 inclusivity/exclusivity

List of recommended pathogens for assay design in the FDA's emergency-use-authorization letter.

Viral pathogen formats	Inactivated / Live	Quantitative Genomic	Synthetic
SARS-CoV-2	2 (heat-killed)	5	4
Human coronavirus 229E	1	1	
Human coronavirus OC43	1	1	
Human coronavirus HKU1			1
Human coronavirus NL63			1
SARS-CoV			1
MERS-CoV			1
Adenovirus	68	7	
Human metapneumovirus			1
Parainfluenza virus 1-4	4	3	
Influenza A & B	82	14	
Enterovirus	118	3	1
Respiratory syncytial virus	9	3	
Rhinovirus	132	6	

Non-viral pathogen formats	Inactivated / Live	Quantitative Genomic	Synthetic
<i>Chlamydia pneumoniae</i>	7	2	
<i>Haemophilus influenzae</i>	73	1	
<i>Legionella pneumophila</i>	35	1	
<i>Mycobacterium tuberculosis</i>	30	2	
<i>Streptococcus pneumoniae</i>	170	1	
<i>Streptococcus pyogenes</i>	155	2	
<i>Bordetella pertussis</i>	30	3	
<i>Mycoplasma pneumoniae</i>	16	2	
<i>Candida albicans</i>	225	2	
<i>Pseudomonas aeruginosa</i>	328	4	
<i>Pneumocystis jirovecii</i>			1*
<i>Staphylococcus epidermis</i>	33	1	
<i>Streptococcus salivarius</i>	11	1*	

\*In development and planning stages.



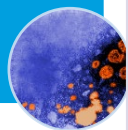
# A menu for assay design

Inclusivity candidate

Exclusivity candidate

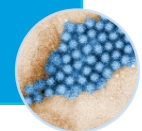
- BK virus
- Hepatitis B virus
- Hepatitis C virus
- Epstein-Barr virus
- Human immunodeficiency virus 1
- Human T-cell leukemia virus
- Human cytomegalovirus
- Varicella-zoster virus
- *Neisseria meningitidis*
- *Plasmodium malariae*
- *Human parechovirus 3*

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*
- Murine norovirus
- *Legionella pneumophila* subsp. *pneumophila*
- Human enterovirus 71 strain H
- *Entamoeba histolytica*
- *E. coli*

Gastro-Intestinal disease



- SARS-CoV-2
- SARS-CoV
- MERS-CoV
- Human coronavirus OC43
- Human coronavirus HKU1
- Human coronavirus NL63
- Human coronavirus 229E
- Human metapneumovirus
- *Bordetella pertussis*

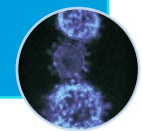
- *Mycobacterium bovis*
- *Mycobacterium talmoniae*
- *Mycobacterium microti*
- *Mycobacterium pinnipedii*

- *Mycobacterium tuberculosis*
- *Streptococcus pneumoniae*

- Human respiratory syncytial virus
- Influenza B virus (Victoria)
- Influenza B virus (Yamagata)
- Influenza A virus (H3N2)
- Influenza A virus (H1N1)
- Human bocavirus

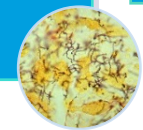
- *Bordetella pertussis*
- *Haemophilus influenzae*
- Adenovirus
- Parainfluenza viruses
- Rhinoviruses
- *Chlamydia pneumoniae*
- *Legionella pneumophila*
- *Mycoplasma pneumoniae*

Respiratory disease



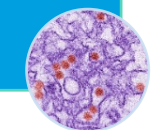
- *Neisseria gonorrhoeae*
- Human immunodeficiency virus 1
- Human papillomavirus 16
- Human papillomavirus 18
- Human papillomavirus 31
- Human T-cell leukemia virus 2
- *Treponema pallidum*
- *Chlamydia trachomatis* LGV I
- *Chlamydia trachomatis* LGV II
- *Chlamydia trachomatis* LGV III
- Human herpesvirus 1
- Human herpesvirus 2
- Hepatitis B virus
- Human herpesvirus 8
- Human herpesvirus 7
- Human herpesvirus 6
- *Mycoplasma genitalium*
- *Staphylococcus saprophyticus*
- *Hamophilus ducreyi*

Sexually transmitted infections



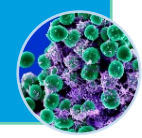
- Powassan virus
- Chikungunya virus
- Dengue virus types 1-4
- Eastern equine encephalitis virus
- *Plasmodium malariae*
- St. Louis encephalitis virus
- West Nile virus
- Yellow fever virus
- Zika virus
- *Borrelia burgdorferi*
- *Plasmodium falciparum*
- Yellow fever virus
- Rift Valley Fever virus

Vector-borne disease



- *Staphylococcus aureus* subsp. *aureus*
- *Staphylococcus epidermidis*
- *Streptococcus pyogenes*
- *Candida albicans*
- *Pseudomonas aeruginosa*
- *Candida krusei*

Epidermal & Nosocomial disease





# Get the materials

*How to find the strains you need*

The screenshot displays the ATCC website's navigation structure. At the top, there are links for 'Resources', 'Support', and 'United States'. On the right, there are links for 'Log in' and 'Create a Profile'. The main navigation menu includes 'CELL PRODUCTS', 'MICROBE PRODUCTS', 'SERVICES', 'FEDERAL SOLUTIONS', 'THE SCIENCE', and 'ABOUT US'. A search bar is located on the right side of the menu. A dropdown menu for 'Microbe Products' is open, listing 'Applications', 'Bacteriology and Archaea', 'Collections and Projects', 'Mycology', 'Protistology', and 'Virology'. A large banner on the right side of the page features the text 'Customize Your Research' and 'Don't find exactly what you need? Let us customize it for you!' with a 'DISCOVER MORE' button.

Resources Support United States

Log in | Create a Profile

ATCC

CELL PRODUCTS MICROBE PRODUCTS SERVICES FEDERAL SOLUTIONS THE SCIENCE ABOUT US

Search

Microbe Products

- Applications
- Bacteriology and Archaea
- Collections and Projects
- Mycology
- Protistology
- Virology

able leads to Incredible

Customize Your Research

Don't find exactly what you need? Let us customize it for you!

DISCOVER MORE

# Get the materials

How to find the strains you need

The screenshot shows the ATCC website interface. At the top, there is a navigation bar with links for Resources, Support, United States (with a dropdown arrow), Log in, and Create a Profile. Below this is a secondary navigation bar with the ATCC logo and links for CELL PRODUCTS, MICROBE PRODUCTS (which is underlined), SERVICES, FEDERAL SOLUTIONS, THE SCIENCE, and ABOUT US. A search bar is located to the right of these links, with a magnifying glass icon and a red arrow pointing to it. Below the navigation bar, there are two columns of menu items. The left column is titled 'Microbe Products' and includes Applications, Bacteriology and Archaea, Collections and Projects, Mycology, Protistology, and Virology (which is underlined). The right column is titled 'Virology' and includes Animal Viruses, Antibodies and Antisera, Bacteriophages, Nucleic Acids (with a red arrow pointing to it), Plant Viruses, Purified Viruses, and Viral Reference Materials. To the right of these menus is a large banner with the text 'Customize Your Research' and 'Don't find exactly what you need? Let us customize it for you!' with a 'DISCOVER MORE' button. The banner features a background image of a person in a lab coat looking at a computer screen. At the bottom of the banner, there are navigation arrows and a series of dots indicating the current slide position.

# Get the materials

How to find the strains you need

## Nucleic Acids

Show per page: 12

PRODUCTS | RESOURCES

97 Products  
Results 1-12 of 97

Search

Refine by

- Product category: Viruses (97)
- Product type: Nucleic acid (+156), **Molecular standard (97)**
- Product application: Assay development (88), Infectious disease re... (81), Next-generation seq... (76), Respiratory disease ... (38)

**Quantitative Genomic DNA Human herpesvirus 2 strain G**  
VR-734DQ BSL 1  
Product format: Frozen  
Derived from: Human herpesvirus 2 G (ATCC VR-734)  
Specification range:  $\geq 1 \times 10^5$  copies/ $\mu$ L  
Price: \$541.00 ea  
Quantity  **ADD TO CART**

**Quantitative Genomic RNA from Human coronavirus 229E**  
VR-740DQ BSL 1  
Product format: Frozen  
Derived from: Human coronavirus 229E 229E (ATCC VR-740)  
Specification range:  $\geq 1 \times 10^5$  copies/ $\mu$ L  
Price: \$541.00 ea  
Quantity  **ADD TO CART** **DOWNLOAD GENOME**

**Quantitative Genomic DNA from Human adenovirus 7 strain Gomen**  
VR-7DQ BSL 1  
Product format: Frozen  
Price: \$541.00 ea  
Quantity  **ADD TO CART**

### Organism

- Influenza A virus (H1... 5
- Human respiratory sy... 4
- Influenza B virus 3
- Human herpesvirus 1 2
- Human enterovirus 71 2
- Search

### Type strain

- No 7

### Nucleic acid type

- Genomic 50
- Synthetic 46

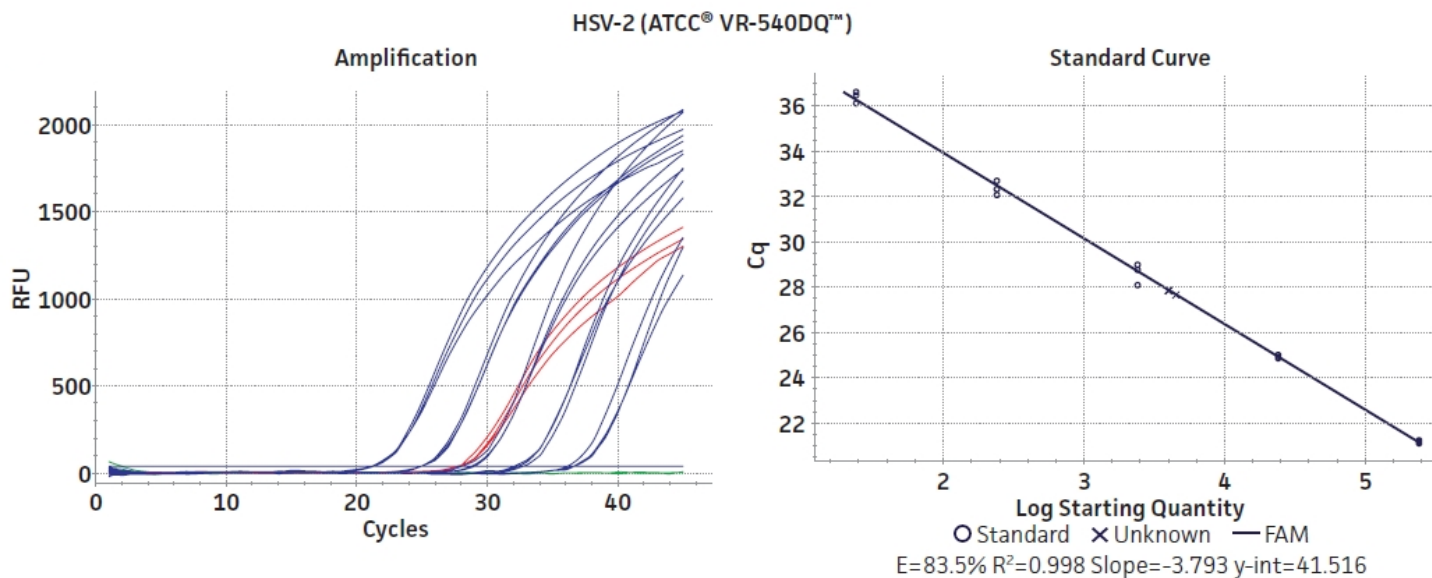
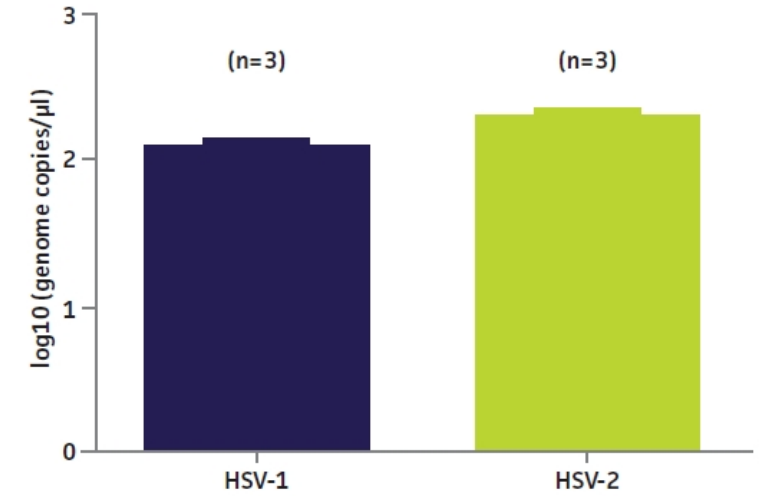
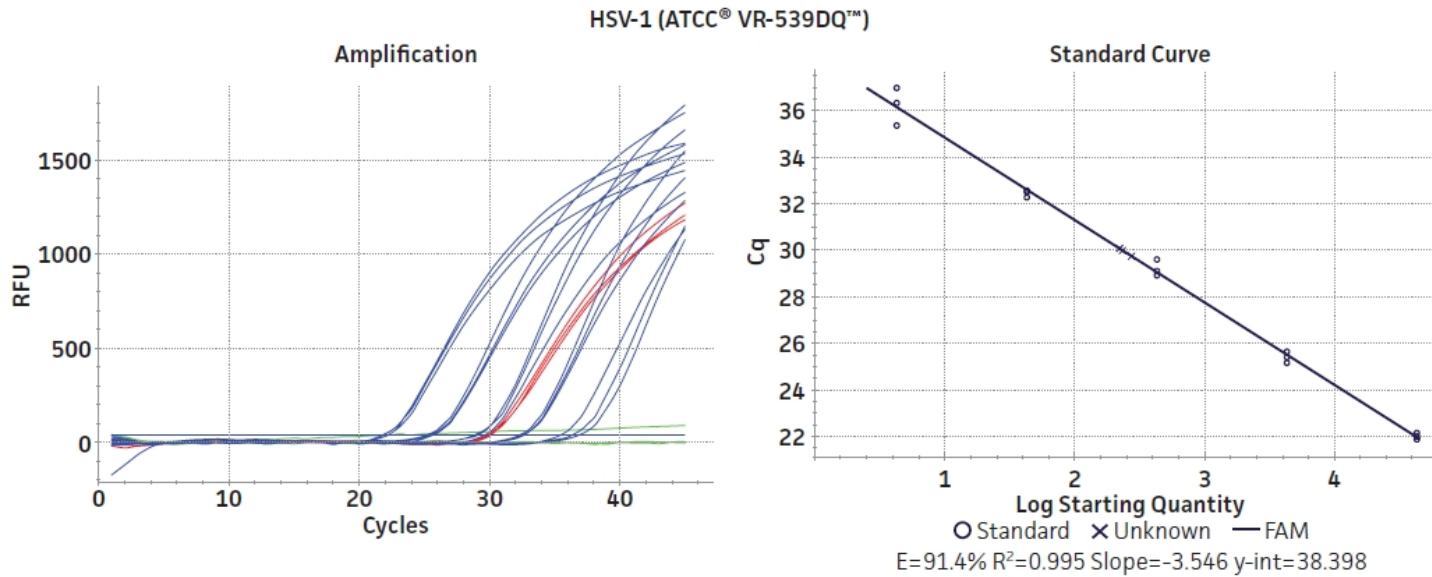
### DNA or RNA

- RNA 67
- DNA 30

# Ongoing and future efforts

- Quality control strains – pharmacopeia, CLSI, or other citations
- Panels for pathogens
  - Respiratory
  - STI
  - Enteric
  - Oncoviruses
  - Opportunistic
  - Anti-microbial resistant
- What do *you* need?
  - ATCC exists to be a resource for scientists.

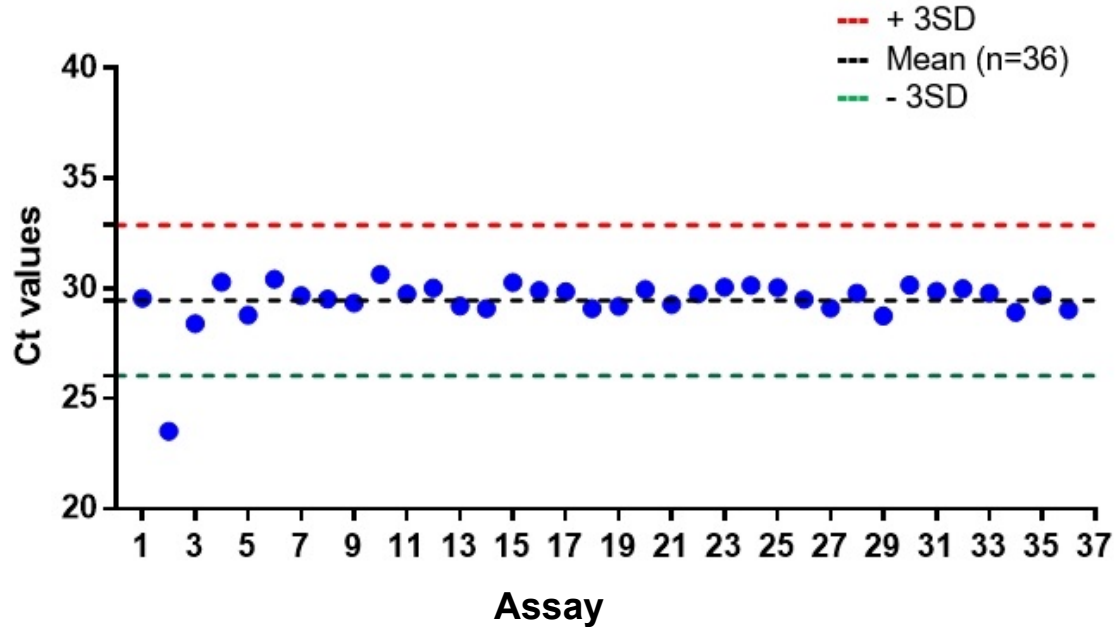
# Assay validation - Human herpes viruses



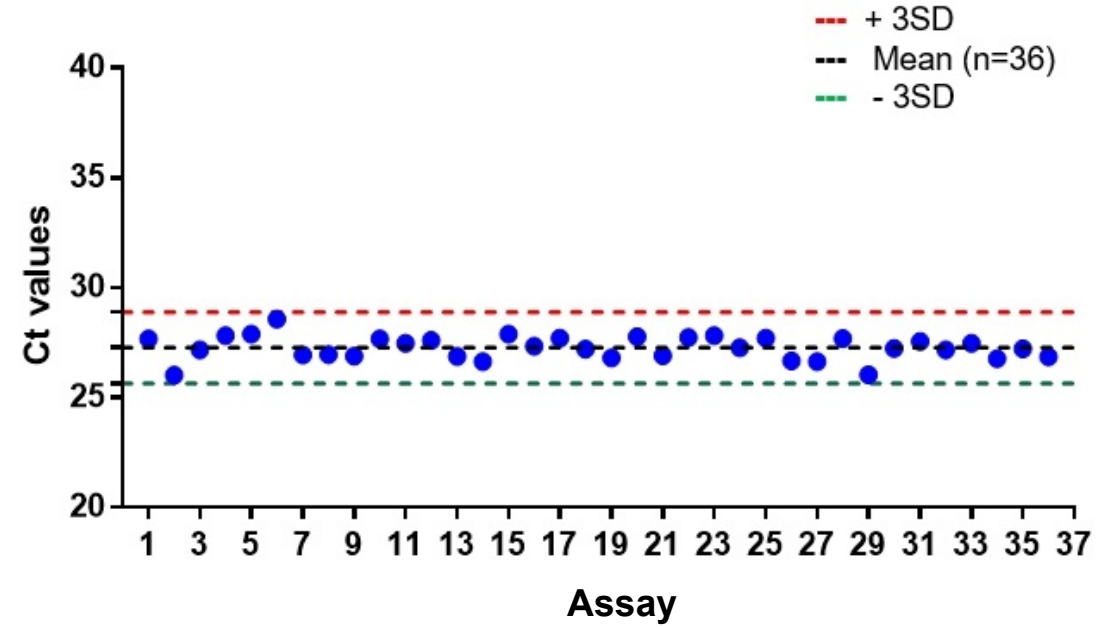
Ryncarz AJ, *et al.* Development of a high-throughput quantitative assay for detecting herpes simplex virus DNA in clinical samples. *J Clin Microbiol* 37(6): 1941-1947, 1999. Pubmed: 10325351

# Assay control - Human herpes viruses

## HSV-1 (ATCC® VR-539D™)



## HSV-2 (ATCC® VR-540D™)



	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%



# Synthetic Molecular Standard Design

# Norovirus standards – feedback and adjustment

Interest in the first 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.

## **Stability**

Changed from dried to frozen

## **Stability**

Added RNA stabilizer

## **Quantification**

Added ddPCR™ to specifications

**ISO 13485**  
guidance for  
manufacturing

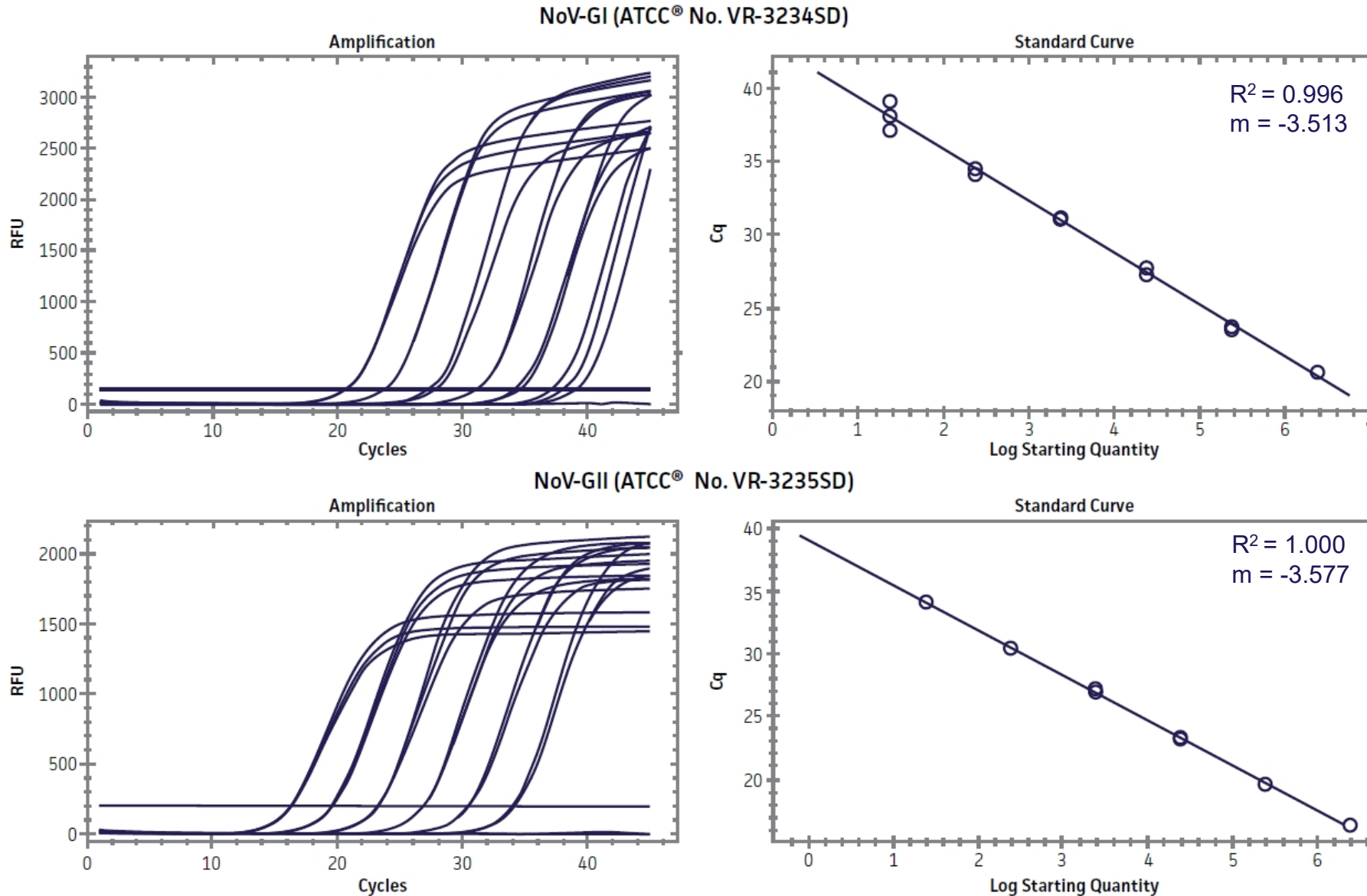
## **Versatility**

Added RdRp fragments to construct



# Validating the next generation of standards

www.atcc.org/2015posters



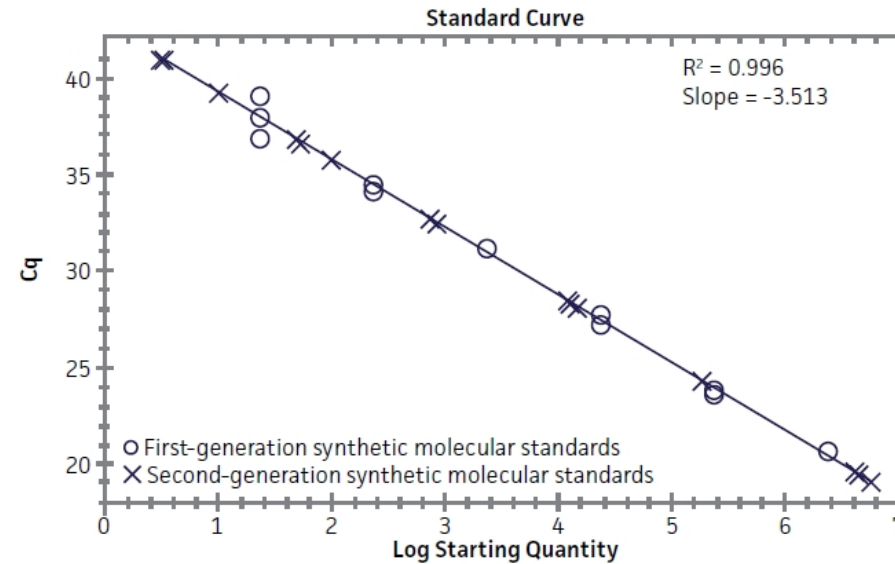
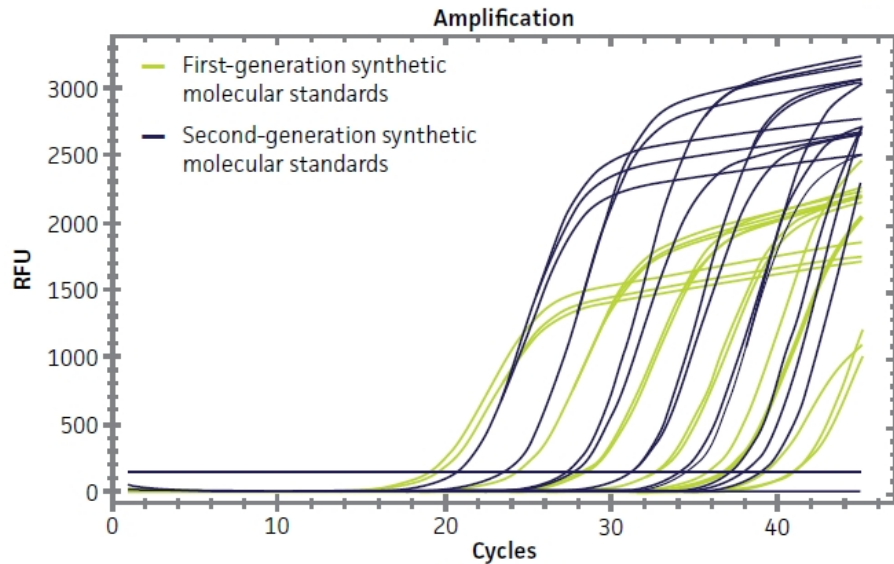
Pictured: Standard curves generated with CaliciNet primer and probe set.

Also tested: ECS working group primer and probe set.

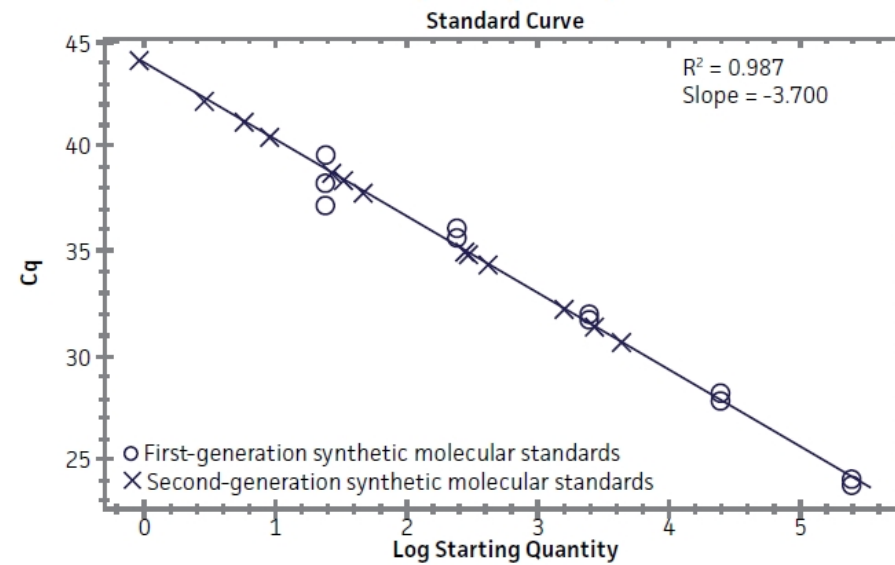
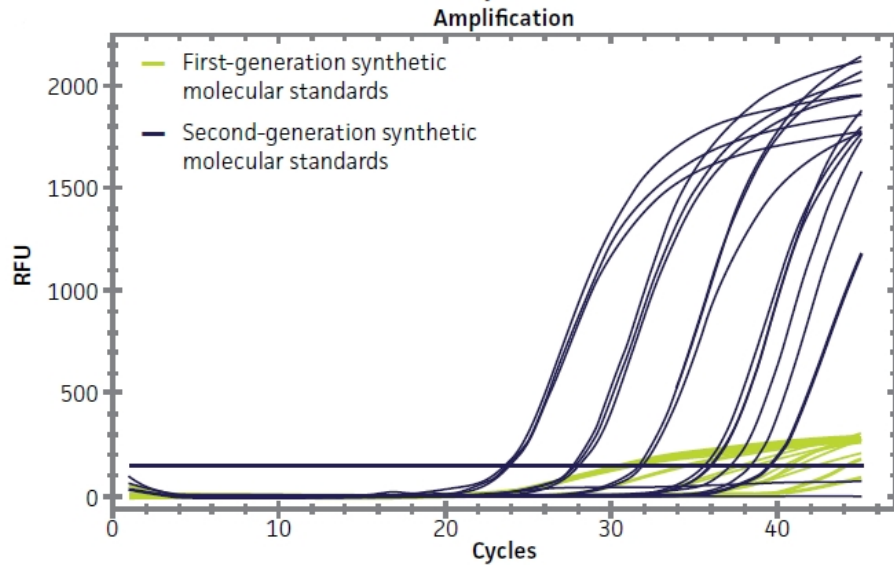
- VR-3234SD™
  - $R^2 = 0.987$
  - $m = -3.692$
- VR-3235SD™
  - $R^2 = 0.998$
  - $m = -3.625$

# Old vs. new standards, Genogroup 1

NoV-GI



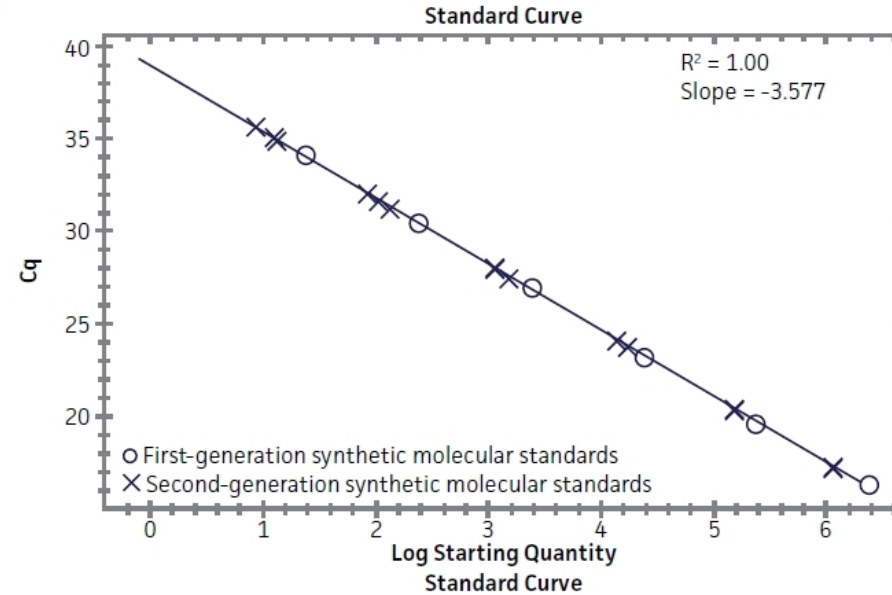
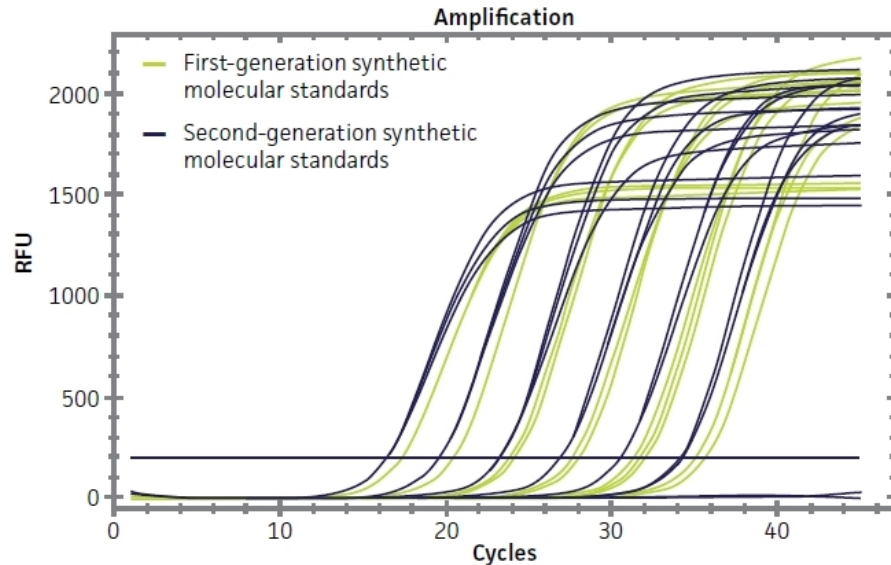
CaliciNet  
primers & probe



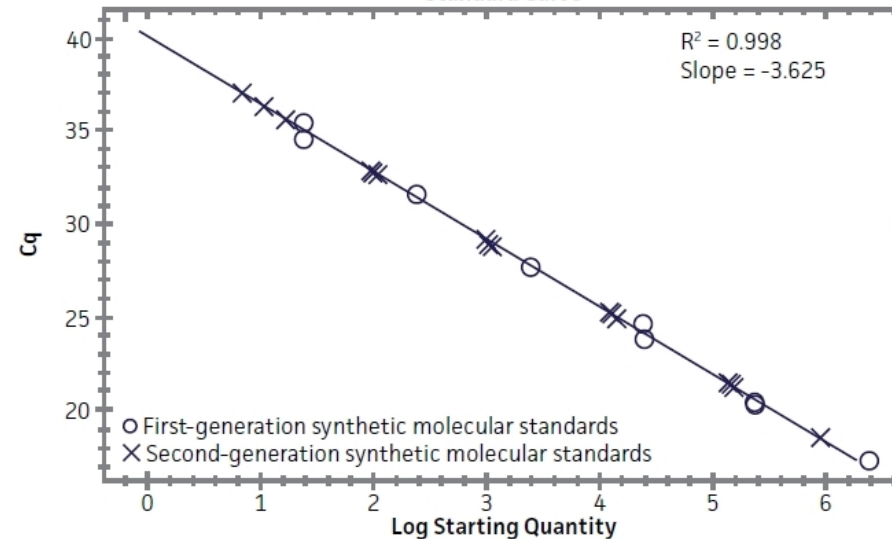
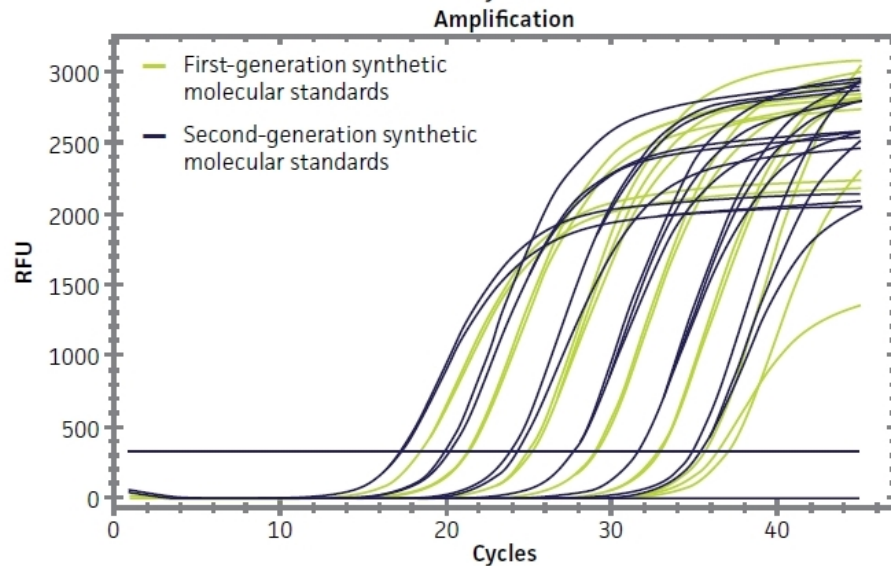
ECS working group  
primers & probe

# Old vs. new standards, Genogroup 2

NoV-GII



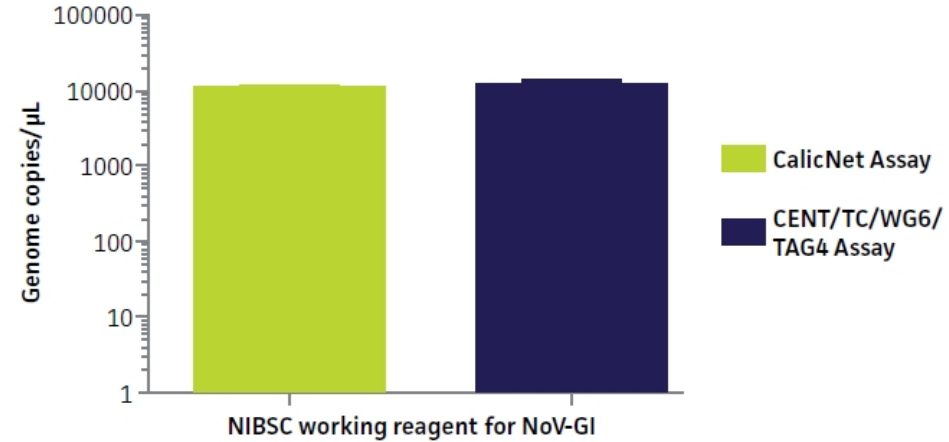
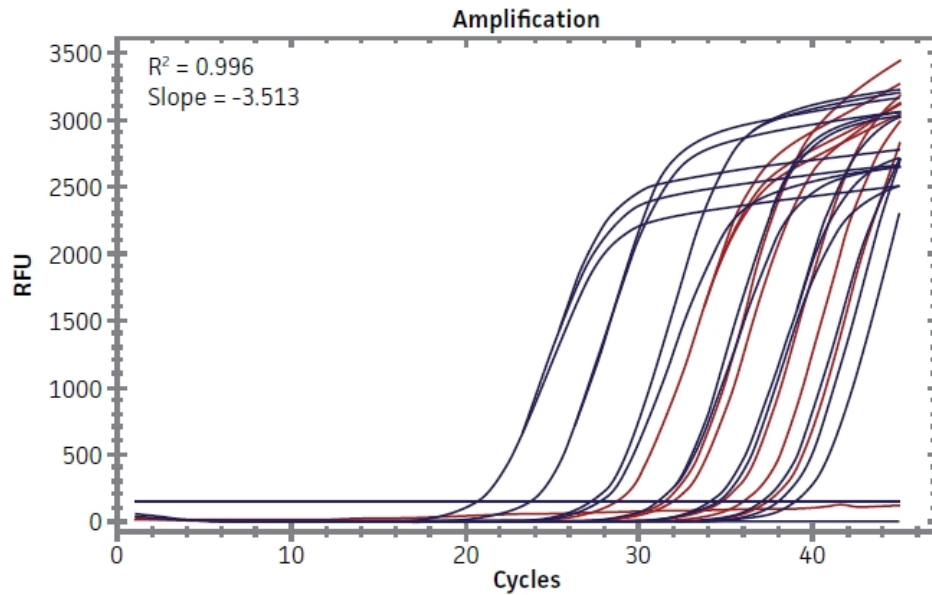
CaliciNet  
primers & probe



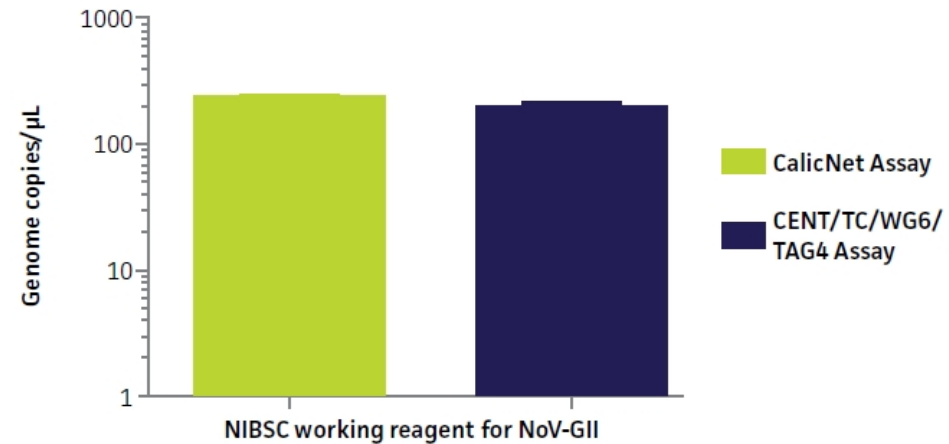
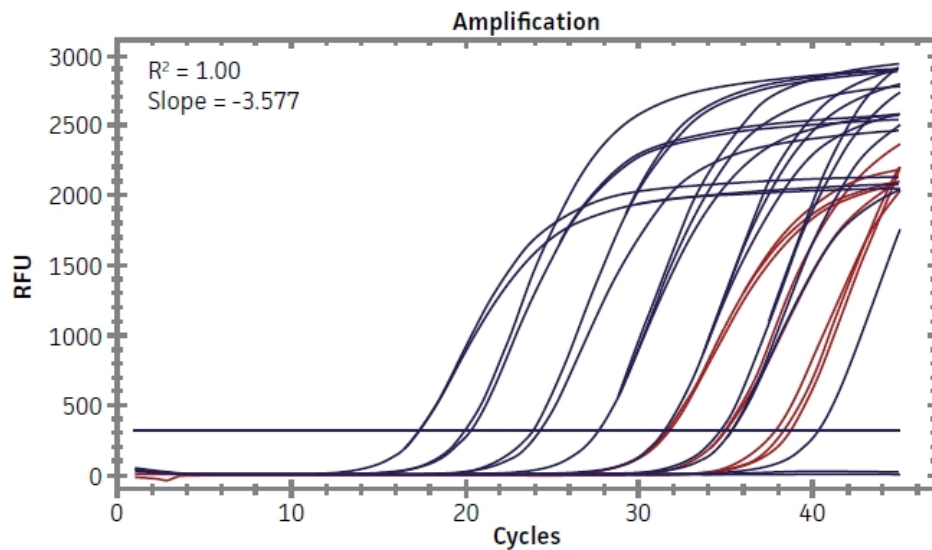
ECS  
primers & probe

# A test drive with NIBSC working reagents

NoV-GI



NoV-GII



# Design approach following design control

**Data collection**

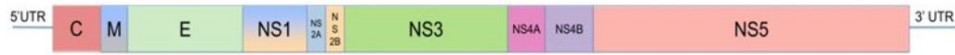
- Literature review
- Bioinformatics sequence data mining
- Technical review

**Construct design**

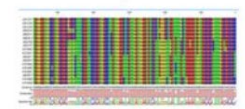
- Construct design
- Assay selection and optimization
- Technical review

**Validation**

- Verify construct specificity
- Technical review
- Quality control testing



Multiple sequence alignment



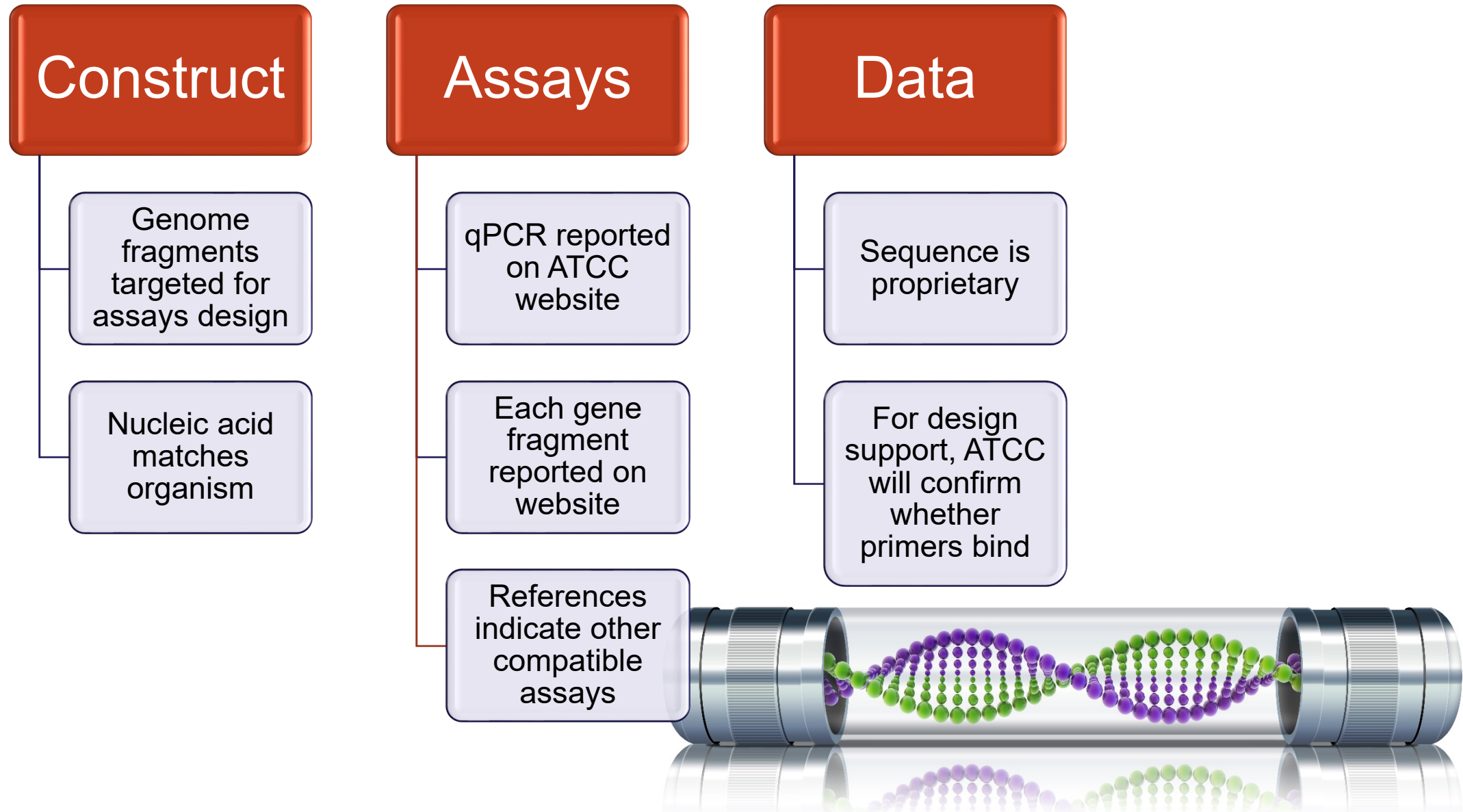
Artificial RNA synthesis



Stabilization



# Design approach – synthetic standards





## Validation of Molecular Standards

# Validation of synthetic standards for hepatitis viruses

## Hepatitis B virus

- ATCC catalog VR-3232SD™
- Hepadnaviridae, Orthohepadnavirus
- DNA construct
- Portions of precure, core, P, S, and X regions

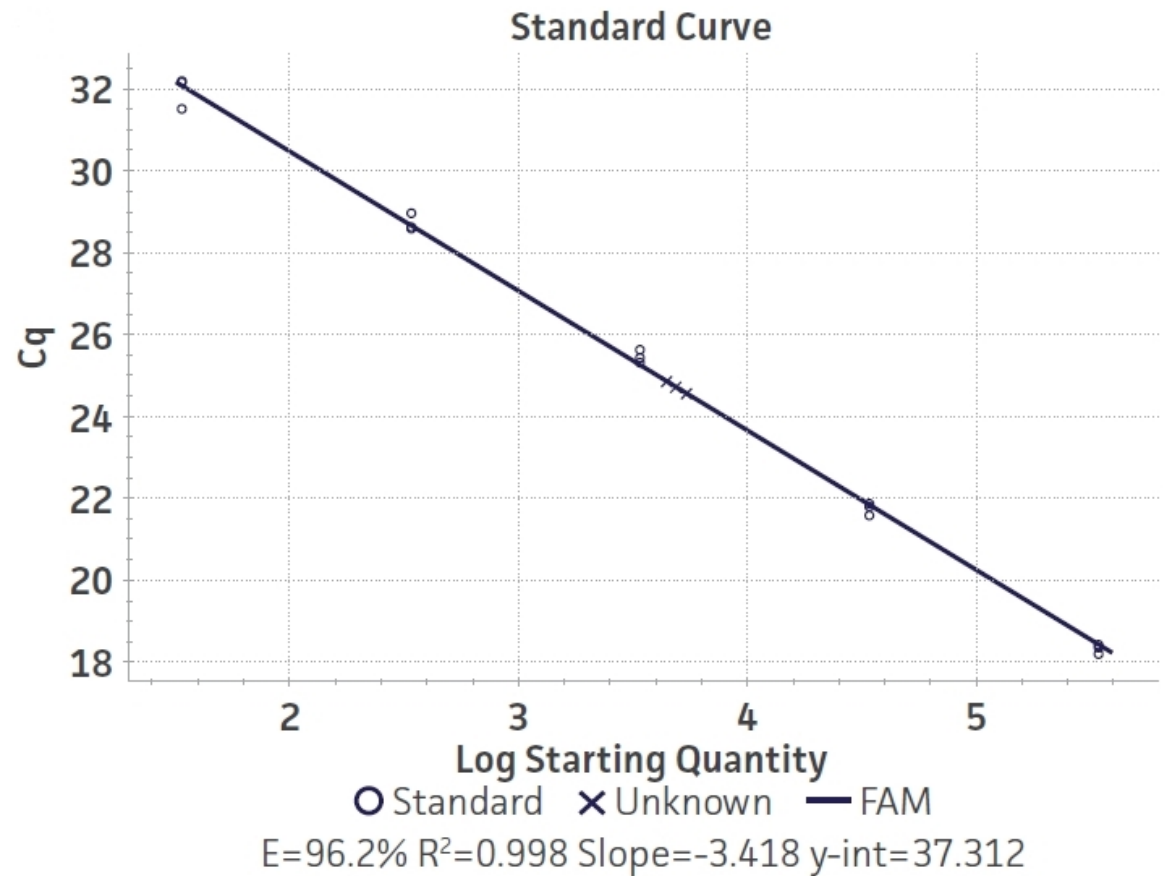
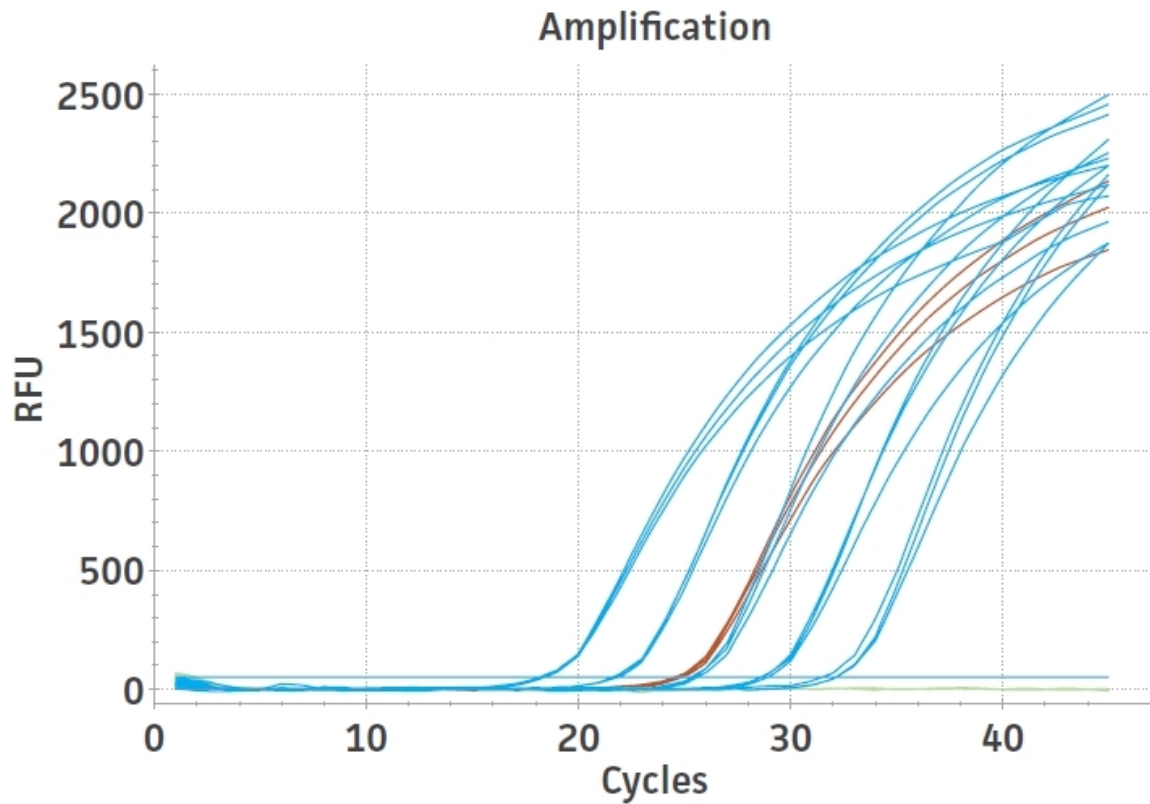
## Hepatitis C virus

- ATCC catalog VR-3233SD™
- Flaviviridae, Hepacivirus, Hepacivirus C
- RNA construct
- Portions of 5' UTR, and X-tail region (3' UTR)

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



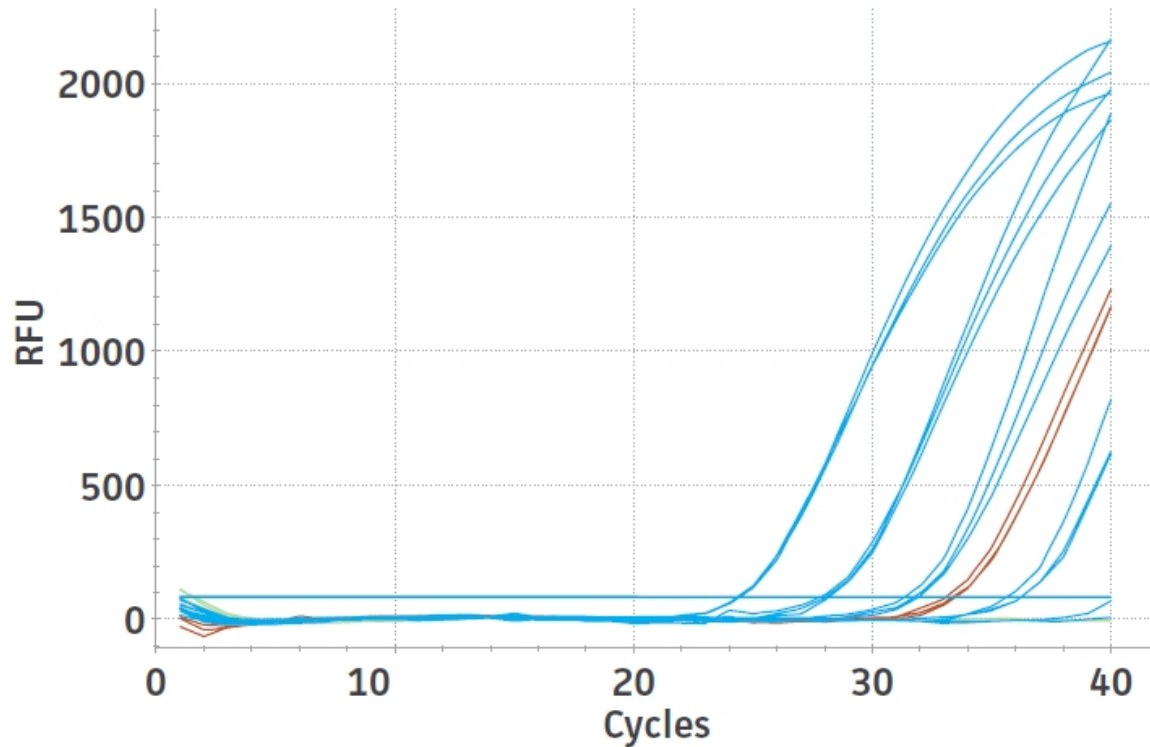
Blue = VR-3232SD™

Red = NIBSC code 10/264 (3<sup>rd</sup> WHO international working reagent for HBV)

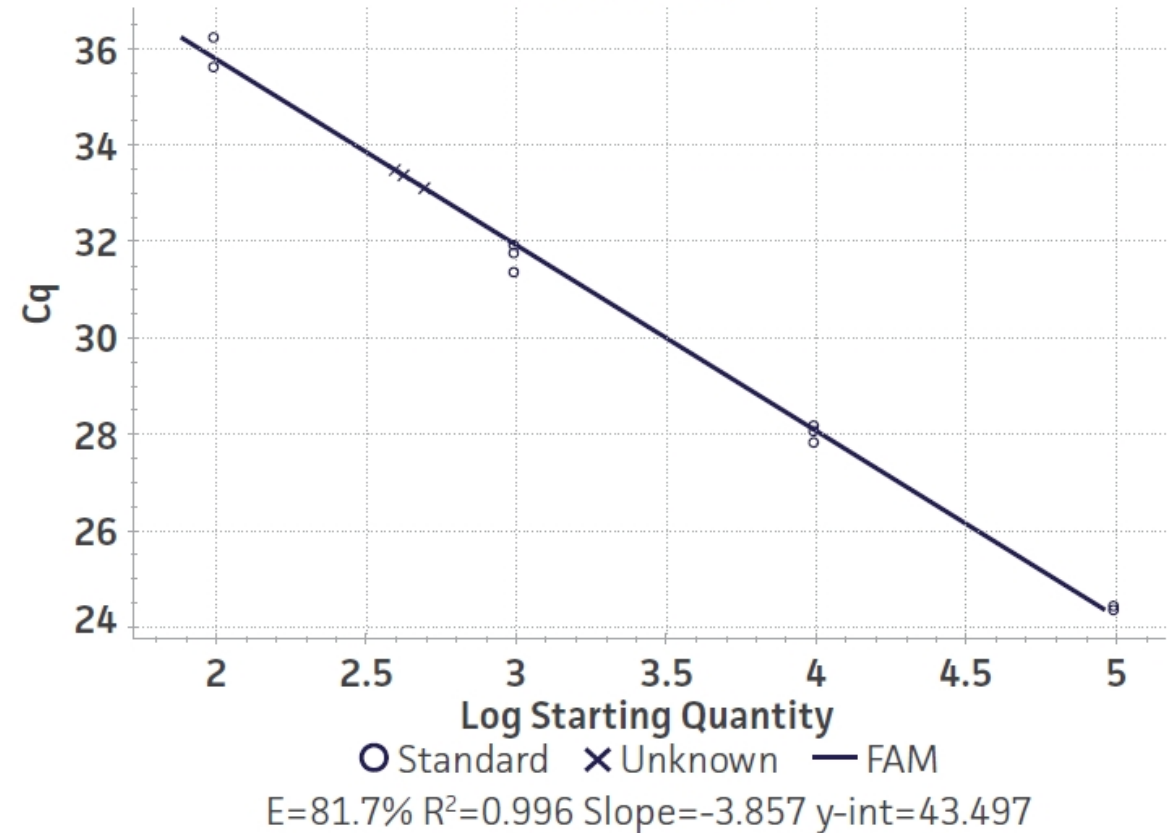
Sun S, *et al.* Development of a new duplex real-time polymerase chain reaction assay for hepatitis B viral DNA detection. *Viol. J.* 8: 227, 2011. PubMed: 21569595

# Hepatitis C virus

Amplification



Standard Curve



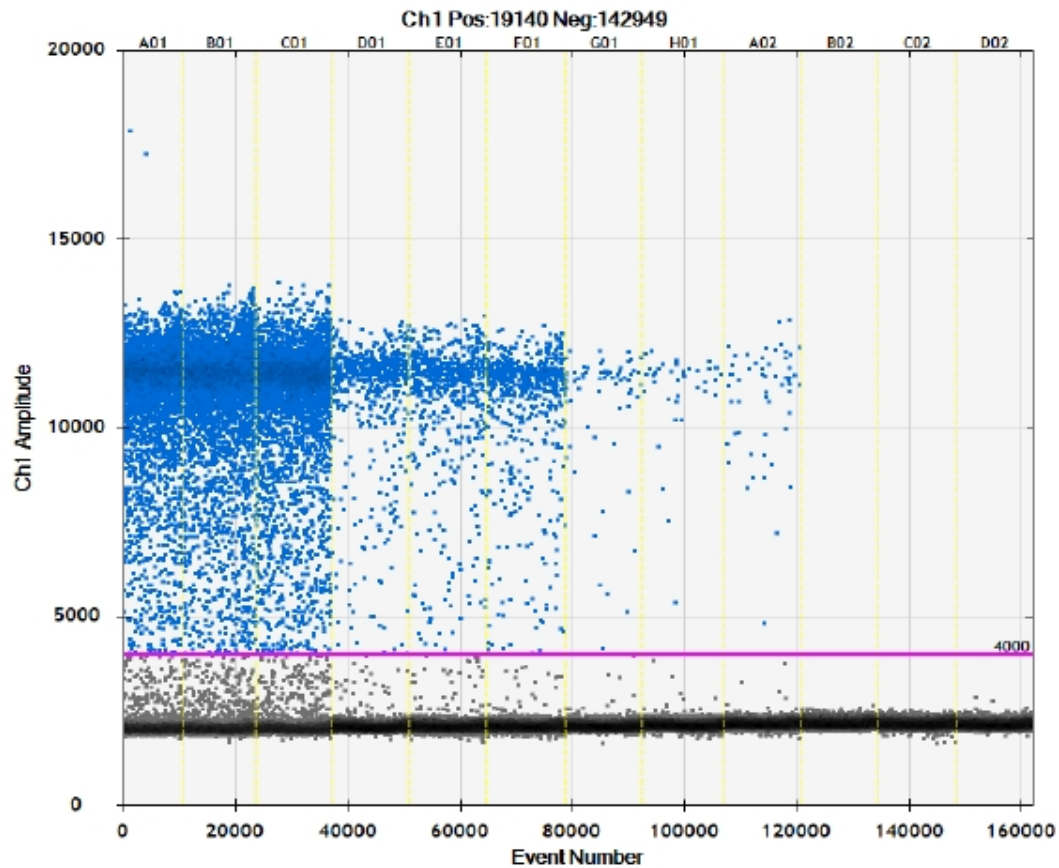
Blue = VR-3233SD™

Red = NIBSC code 06/102 (4<sup>th</sup> WHO international standard for HCV)

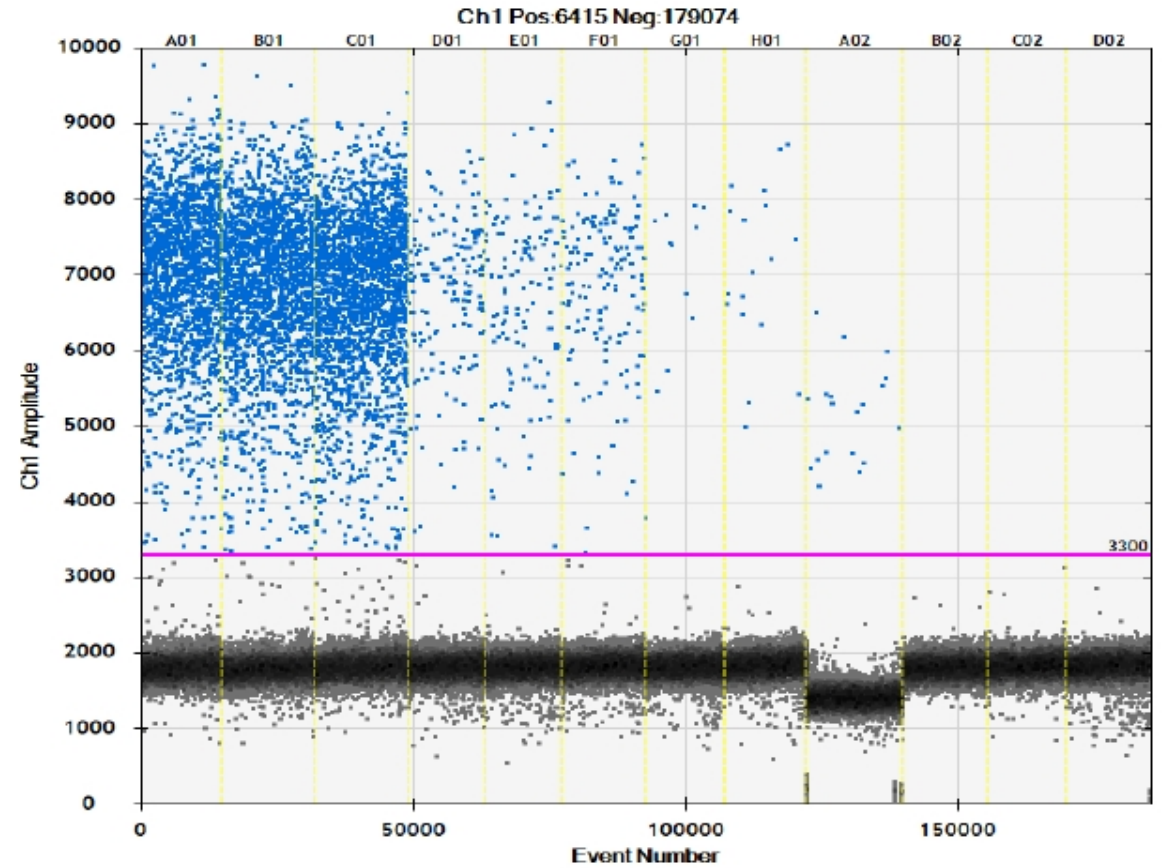
Lee SC, *et al.* Improved version 2.0 qualitative and quantitative AMPLICOR reverse transcription-PCR tests for hepatitis C virus RNA: calibration to international units, enhanced genotype reactivity, and performance characteristics. *J. Clin. Microbiol.* 38(11): 4171-4179, 2000. PubMed: 11060086

# Hepatitis viruses

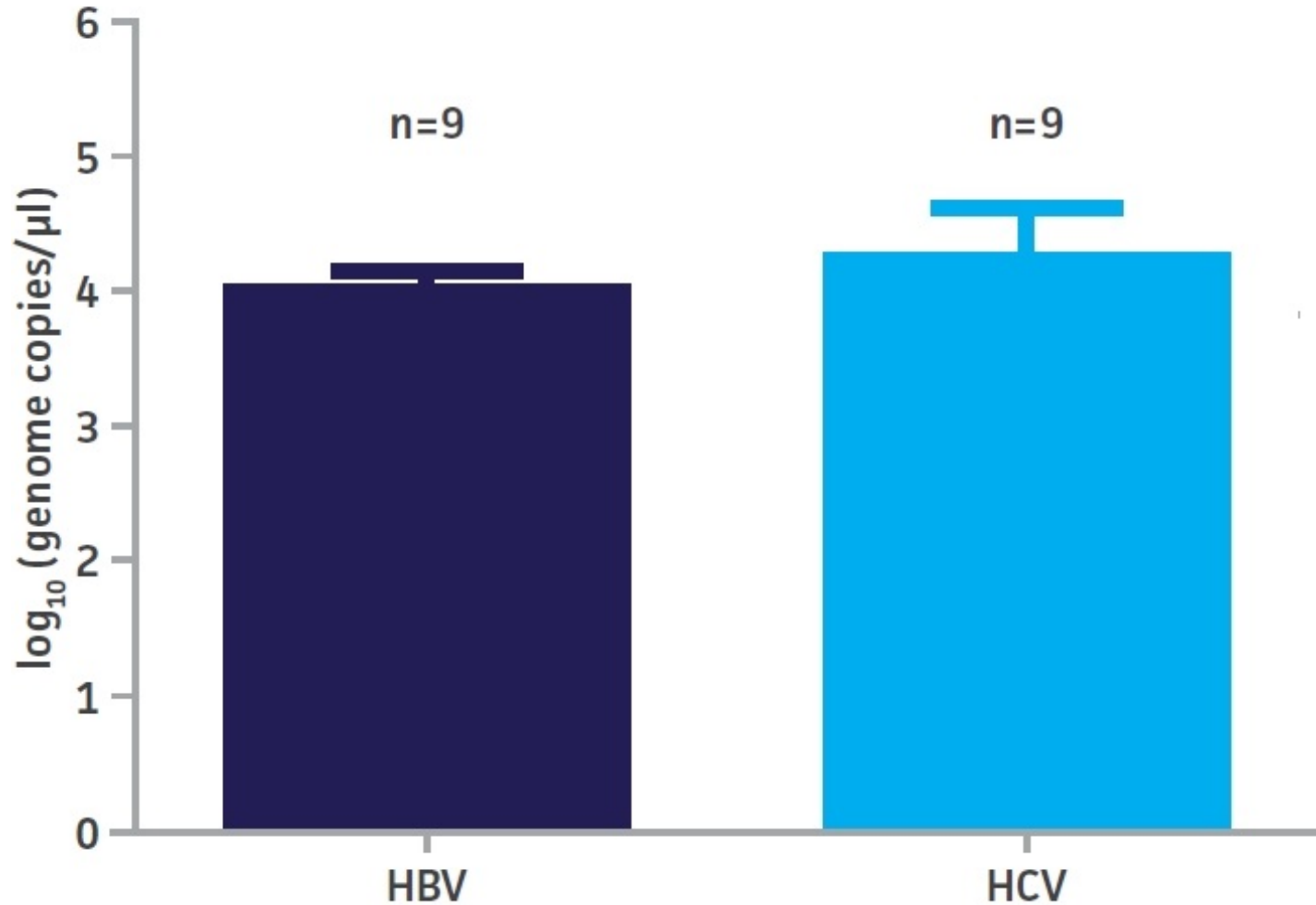
VR-3232SD™ (HBV synthetic standard)



VR-3233SD™ (HCV synthetic standard)



# Quantitation of NIBSC Hepatitis standards



As determined by the WHO:

- HBV standard =  $8.5 \times 10^5$  IU/mL
- HCV standard =  $2.6 \times 10^5$  IU/mL

qRT-PCR and qPCR quantitation at ATCC:

- HBV:  $9.7 \times 10^6$  genome copies/mL
- HCV:  $1.6 \times 10^7$  genome copies/mL

Conversion ratio as quantified at ATCC:

- HBV: 1 IU/mL = 11.4 genome copies
- HCV: 1 IU/mL = 61.5 genome copies

# Other application data and posters

## Computational Design of a Synthetic Molecular Standard for Human Parechovirus 3

Bitray Tang, B.S., Michael Geimer, M.S., Maria Mayda, Ph.D., and Dev Mittar, Ph.D.  
ATCC, Manassas, VA 20110

### Synthetic Human Parechovirus 3 RNA (ATCC® VR-3260SD™)

Advantages	Applications
<ul style="list-style-type: none"> <li>Manufactured and authenticated with ISO13485:2010 compliance</li> <li>BSL-1 ready-to-use control</li> <li>No shipping restrictions</li> <li>Quantitative format</li> <li>Stabilized RNA</li> </ul>	<ul style="list-style-type: none"> <li>Generation of a standard curve for quantitative RT-PCR</li> <li>Positive control for RT-PCR assays</li> <li>Assay verification and validation studies</li> <li>Monitor assay-to-assay and lot-to-lot variation</li> <li>Molecular diagnostics assay development</li> </ul>

### Background & Introduction

Human parechovirus 3 (HPeV3) has been increasingly identified in cases of aseptic meningitis among neonates and young infants less than 1 year of age, and is associated with paralysis, viral-like sepsis, central nervous system (CNS) infection, and sudden death. Because these clinical manifestations are similar to those associated with enterovirus infections, HPeV3 infections are often misdiagnosed, which in turn results in poor patient outcome. Therefore, molecular detection assays that provide a rapid and accurate diagnosis of HPeV3 are critical for ensuring prompt and appropriate treatment. Due to its sensitivity and quick turnaround time, the preferred method for the detection of RNA viruses is quantitative reverse-transcription PCR (qRT-PCR), which relies on the generation of a standard curve that is prepared using a quantitative viral RNA standard. To this end, we have designed, developed, and quantified a synthetic molecular standard for HPeV3 that is compliant with ISO 13485. This preparation is supported by stringent quality control analysis to ensure product identity, stability, and functionality with molecular applications, making it an ideal control for assessing assay performance and ensuring accurate and reproducible results. In the following proof-of-concept study, the HPeV3 synthetic molecular standard was quantified using Droplet Digital™ PCR (ddPCR™, Bio-Rad) and validated via qRT-PCR using published primers.

### Computational Design Strategy

Reference Sequence: HPeV3 Strain BJ-C317 Genome

- Genome sequence data collection
- Multiple sequence alignment
- Phylogenetic analysis
- Sequence editing and assembly
- Artificial RNA synthesis
- Sequence verification and authentication via next-generation sequencing
- Stabilization of RNA
- Absolute quantification via ddPCR™ assay
- Validation using qRT-PCR assay

### Synthetic Design

Primer and Probe Sequences!

Forward Primer Sequence (5' to 3'): GAC AAC ATC TTT GGT AGA GGT TGG T  
Reverse Primer Sequence (3' to 5'): TTT TGC CTC CAG GTA TCC CAG T  
Probe Sequence (5' to 3'): TGT GAT CTC TGA TGA TTT T

ATCC 10801 University Boulevard, Manassas, Virginia 20110 Phone: 800.638.6597 Email: SalesRep@atcc.com Web: www.atcc.org

## Development and Evaluation of Quantitative Synthetic and Genomic Molecular Standards for Zika

Helen Christina, M.S., Sujatha Rashid, Ph.D., Dev Mittar, Ph.D.  
ATCC, Manassas, VA

### Background and Introduction

Early detection of the Zika virus (ZIKV), an emerging mosquito-borne pathogen, in infected people is of paramount importance for patient management and for curbing viral spread. Currently, one of the most reliable molecular-based methods used for viral detection during acute infection is quantitative reverse transcription polymerase chain reaction (qRT-PCR). To aid in the development and quality control of these assays, reliable positive control materials are essential for determining analytical specificity and sensitivity. To that end, we have developed quantitative synthetic and genomic ZIKV RNA standards that can be used for assay development and validation (Table 1).

### Advantages and Applications of the ZIKV RNA Standards

Advantages	Applications
<ul style="list-style-type: none"> <li>Fully authenticated &amp; characterized</li> <li>Quantified by ddPCR™</li> <li>Consistent and accurate results</li> <li>BSL-1 ready-to-use control</li> <li>No shipping restrictions</li> </ul>	<ul style="list-style-type: none"> <li>Generation of a standard curve for qRT-PCR</li> <li>Positive control for qRT-PCR assays</li> <li>Independent standard for validation and verification studies</li> <li>Monitor assay-to-assay and lot-to-lot variation</li> </ul>

### Results

#### Standard Curve Generation Using the Genomic and Synthetic ZIKV RNA Standards

Figure 1. Generation of Standard Curves from Genomic and Synthetic ZIKV RNA Standards. (A) qRT-PCR amplification plot showing the synthetic ZIKV RNA standard (ATCC® VR-3252SD™, Blue), the ZIKV working reagents (color), and a negative control (Black). (B) A qRT-PCR amplification plot showing the genomic ZIKV RNA standard (ATCC® VR-1830™, Blue), the ZIKV working reagents, and a negative control (Black).

#### Quantification of Native ZIKV using Synthetic and Genomic RNA Standards

Figure 2. Quantification of ATCC ZIKV Working Reagents using ATCC Quantitative Genomic and Synthetic ZIKV RNA Standards. (A) qRT-PCR amplification plot showing the synthetic ZIKV RNA standard (ATCC® VR-3252SD™, Blue), the ZIKV working reagents (color), and a negative control (Black). (B) A qRT-PCR amplification plot showing the genomic ZIKV RNA standard (ATCC® VR-1830™, Blue), the ZIKV working reagents, and a negative control (Black).

#### Genomic ZIKV

Figure 3. Generation of Standard Curves from Genomic ZIKV RNA Standards. (A) qRT-PCR amplification plot showing the genomic ZIKV RNA standard (ATCC® VR-1830™, Blue), the ZIKV working reagents (color), and a negative control (Black). (B) A qRT-PCR amplification plot showing the synthetic ZIKV RNA standard (ATCC® VR-3252SD™, Blue), the ZIKV working reagents, and a negative control (Black).

### Materials and Methods

#### Quantitative Synthetic ZIKV RNA

The synthetic ZIKV RNA standard was developed using an artificial RNA synthesis method and contains fragments from the membrane glycoprotein precursor M, envelope NS1, NS2B, NS3, NS4B, and NS5 regions of the ZIKV genome. The RNA was quantified using Droplet Digital™ PCR (ddPCR™, Bio-Rad) to determine genome copy number and was stabilized with RNAsin® (Qiagen).

#### Quantitative Genomic ZIKV RNA

Genomic RNA was isolated from a preparation of cell lysate and supernatant from Vero cells (ATCC® CCL-81™) infected with ZIKV strain MR 766 (ATCC® VR-1830™). The RNA was quantified using ddPCR™ to determine genome copy number.

#### ZIKV Working Reagents

Catalog Number	Product Description
ATCC® VR-1830™	IBH 30656 (Human/1993/Nigeria)
ATCC® VR-1843™	PH/ABC29 (Human/2015/Puerto Rico)
ATCC® VR-1844™	FLR (Human/2015/Columbia)
NR-50119 (BEI Resources)	FLR (Human/2015/Columbia)
NR-50219 (BEI Resources)	HPAN/2015/GDC-296359 (Human/2015/Panama)

### Disclaimers

© 2017 American Type Culture Collection. All rights reserved. This product is for research use only. It is not for clinical use. The information contained herein is for informational purposes only. It is not intended to be used for diagnosis or treatment of any disease. The information contained herein is not intended to be used for diagnosis or treatment of any disease. The information contained herein is not intended to be used for diagnosis or treatment of any disease.

## Development of Synthetic Molecular Standards for Dengue Virus

Shamaila Ashraf, Melissa Wilson, Afshin Sohrabi, Stephen King, Brian Chase, Dev Mittar, Kurt Langenbach and Andrew G. Cavthorn

### Background & Introduction

Dengue fever is an acute illness caused by any one of four serotypes (1-4) of genetically related dengue viruses (DENV), with an estimated 300 million cases reported annually. Currently, quantitative RT-PCR (qRT-PCR) is the preferred method for the detection and quantification of DENV in clinical diagnostics and epidemiological surveillance. The accuracy of a qRT-PCR assay relies on the generation of a standard curve using a positive control with a known viral genome concentration.

### ATCC Design Strategy

Figure 1. An example of a qRT-PCR amplification plot and standard curve generated using the DENV-4 molecular standard (ATCC® VR-3212SD™) in comparison with the qRT-PCR primer and probe set. (C) The slope and R<sup>2</sup> values generated using DENV-1, 2 and 4 molecular standards with the primer and probe sets for the qRT-PCR for the detection, quantification, and serotyping of dengue viruses. *PLoS Negl Trop Dis* 7(4): e2110, 2013.

### Generation of Standard Curves from DENV Synthetic Molecular Standards

Figure 2. An example of qRT-PCR amplification plot showing DENV-4 synthetic molecular standards (blue) and the native sample concentration (pink). Unknown samples were serially diluted 10-fold and run in triplicate wells. (B) Copy numbers of DENV-1, 2 and 4 samples as determined by the qRT-PCR standards curves generated by using DENV-1, 2 and 4 molecular standards respectively.

### Quantification of Native DENV-1, 2 and 4 Using Synthetic Molecular Standards

Figure 3. An example of qRT-PCR amplification plot showing DENV-4 synthetic molecular standards (blue) and the native sample concentration (pink). Unknown samples were serially diluted 10-fold and run in triplicate wells. (B) Copy numbers of DENV-1, 2 and 4 samples as determined by the qRT-PCR standards curves generated by using DENV-1, 2 and 4 molecular standards respectively.

### ATCC 10801 University Boulevard, Manassas, Virginia 20110-2209 phone: 800.638.6597 email: sales@atcc.org www.atcc.org



# Summary

- ATCC Molecular Standards are a fast, easy, reliable control for assay development and validation & control.
- Genomic standards eliminate the costs of growth, extraction, and quantitation.
- Synthetic standards provide controls for organisms that are difficult to culture or extract.
- Over 230 standards currently in the portfolio.
  - Over 170 genomic standards
  - Over 60 synthetic standards
  - Standards for pathogens, microbiome, & food safety
  - And more to come!

# 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, Ph.D.***



Questions?

Credible Leads to Incredible™

