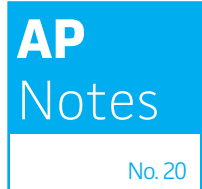


# DEVELOPMENT OF IMPROVED SYNTHETIC MOLECULAR STANDARDS FOR NOROVIRUS GENOGROUPS I AND II

Shamaila Ashraf, Ph.D., Prasanthi Bandi, M.S., Brian Chase, M.S., and Dev Mittar, Ph.D.  
ATCC, Manassas, VA



## ABSTRACT

ATCC has improved existing first-generation qualitative synthetic molecular standards for Norovirus genogroups I and II (NoV-GI and NoV-GII, respectively) through the development of two new quantitated synthetic molecular standards. These second-generation synthetic RNA standards are synthesized under ISO 13485:2003 guidance and comprise short conserved fragments representing the NoV genome. Further, they are compatible with numerous published quantitative reverse transcription polymerase chain reaction (qRT-PCR) assays, can be handled at biosafety level one (BSL-1), are stable, and can be shipped internationally.

## INTRODUCTION

NoVs are the most common cause of epidemic gastroenteritis, accounting for 95% of viral gastroenteritis outbreaks worldwide. These viruses are difficult to detect due to their genetic heterogeneity and unculturable nature. Currently, the principle detection method utilized by diagnostic laboratories is qRT-PCR. The accuracy of a qRT-PCR assay relies on the generation of a standard curve using a positive control with a known genome copy number. In addition, clinical laboratories require an independent control to monitor the accuracy of their assays. Since Norovirus cannot be grown in cell culture, the acquisition of such positive control material proves challenging. To circumvent these issues, ATCC has developed quantitative synthetic molecular standards that include conserved sequences from NoV-GI and NoV-GII for the detection and quantification of NoV from either clinical, food, or environmental samples (Table 1).

These novel synthetic molecular standards are an improvement over existing first-generation qualitative synthetic molecular standards for NoV-GI and NoV-GII in that they are synthesized under ISO 13485:2003 guidance, quantified by Droplet Digital™ PCR (ddPCR™, Bio-Rad®), and stabilized by the addition of RNastable® (Biomatrix®). Further, the improved design of these second-generation constructs enables compatibility with additional molecular assays for the detection and quantification of NoV. In this study, we used the second-generation synthetic molecular standards for NoV-GI and NoV-GII to generate standard curves using two different primer sets; one used by the laboratories in the CaliciNet<sup>1</sup>, and another recommended by the European Committee for Standards (CEN/TC/WG6/TAG4) working group<sup>2</sup>. Furthermore, we used the respective standard curves to quantify the working reagents for NoV-GI and NoV-GII from National Institute for Biological Standards and Control (NIBSC) using the above primer sets.

**Table 1: ATCC synthetic NoV molecular standards**

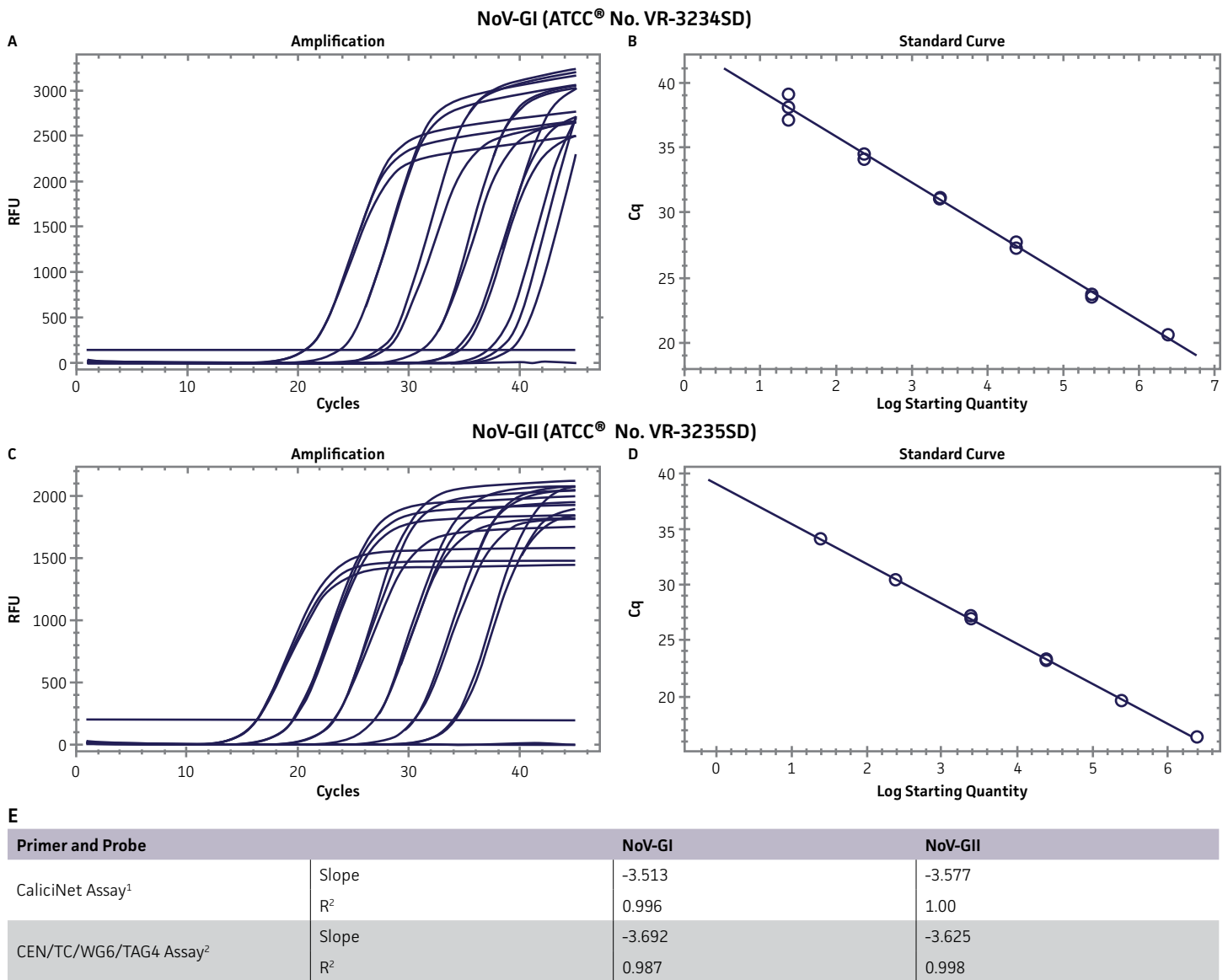
ATCC® No.	Designation
VR-3234SD	Synthetic NoV-GI RNA
VR-3235SD	Synthetic NoV-GII RNA

## DEVELOPMENT OF THE SYNTHETIC MOLECULAR STANDARDS

The quantitative synthetic molecular standards for NoV were manufactured under ISO 13485:2003 guidance and include respective fragments from the ORF1 and ORF2 junction for NoV-GI and NoV-GII, and an additional fragment from the VP2 region for NoV-GII. Both synthetic RNA standards were quantified by ddPCR in order to package precise copies of RNA. Given the inherent labile nature of RNA, a stabilization matrix (RNastable) was added to the quantitated RNA preparations.

## GENERATION OF STANDARD CURVES FROM SECOND-GENERATION NOV SYNTHETIC MOLECULAR STANDARDS

Standard curves were generated using serial ten-fold dilutions of the synthetic RNA standards, ranging from 20 copies to  $2 \times 10^6$  copies/reaction (Figure 1). Standards were tested in triplicate wells using the CFX96™ Real-Time PCR Detection System (Bio-Rad). The primer and probe sets from the Real-Time RT-PCR assay used by laboratories in CaliciNet<sup>1</sup> and an assay recommended CEN/TC/WG6/TAG4 group<sup>2</sup> were employed to examine each individual NoV genotype using the following conditions. Cycling conditions for all primer sets were 50°C for 15 min and 95°C for 2 min, followed by 45 cycles of 95°C for 15 sec and 60°C for 60 sec for both the NoV-GI and NoV-GII assays. RNA samples and standards were tested in triplicate. The relative fluorescence unit (RFU) baseline threshold was set and genome copy numbers calculated using CFX Manager™ 3.0 Software (Bio-Rad). From this analysis, it was determined that the second-generation synthetic NoV molecular standards are compatible with both primer sets over a broad linear dynamic range and exhibit minimal variability as evident from slope and R<sup>2</sup> values (Figure 1E).



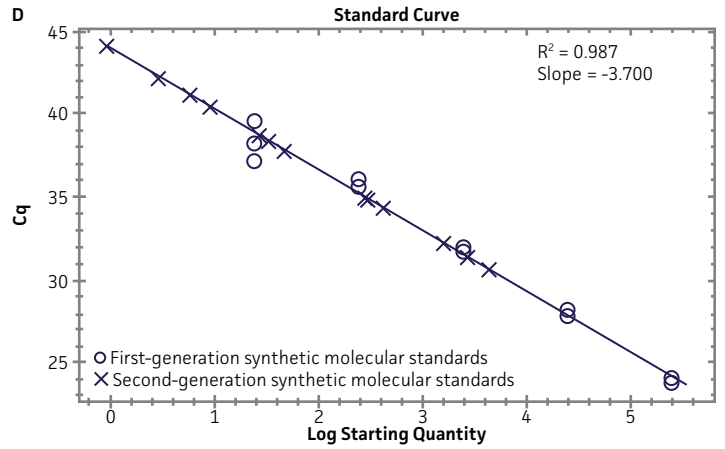
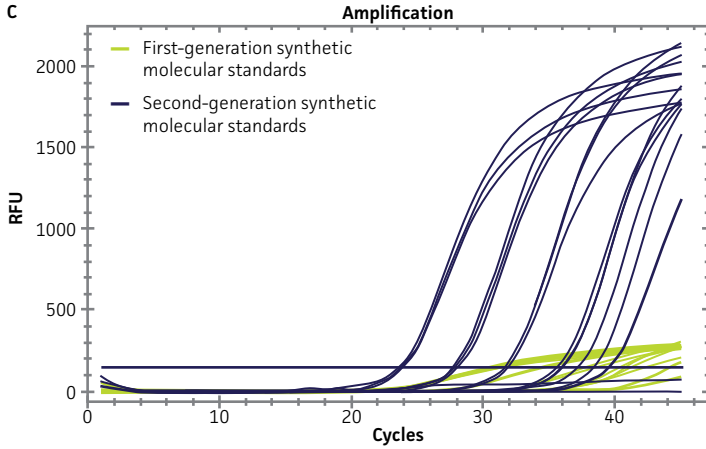
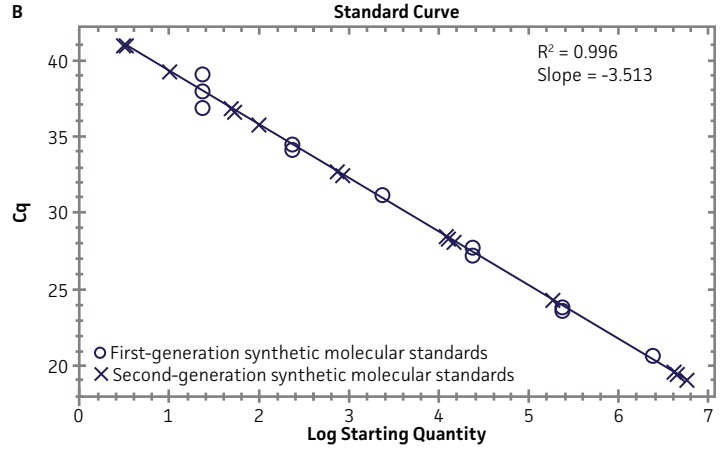
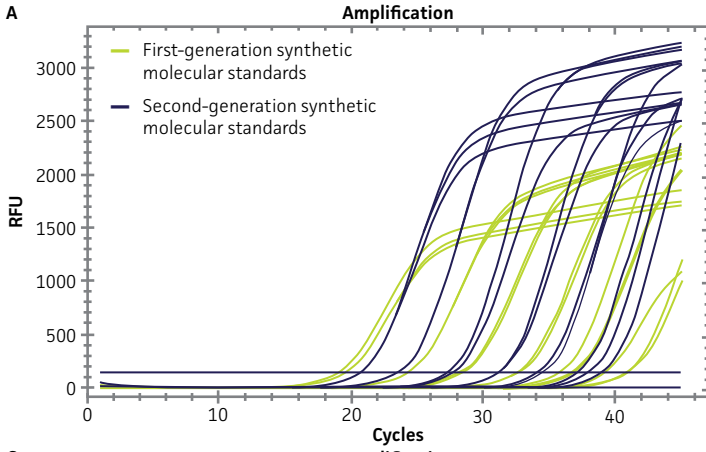
**Figure 1.** Generation of Standard Curves from Second-Generation NoV Synthetic Molecular Standards

A&C) Amplification plots and B&D) standard curves generated using the second-generation NoV-GI and NoV-GII synthetic molecular standards (ATCC® VR-3234SD and ATCC® VR-3235SD, respectively) with the CaliciNet primer and probe set<sup>1</sup>. E) The slope and R<sup>2</sup> values generated using the second-generation NoV-GI and NoV-GII synthetic molecular standards with the primer and probe sets from the CaliciNet<sup>1</sup> and CEN/TC/WG6/TAG4<sup>2</sup> real-time RT-PCR assays.

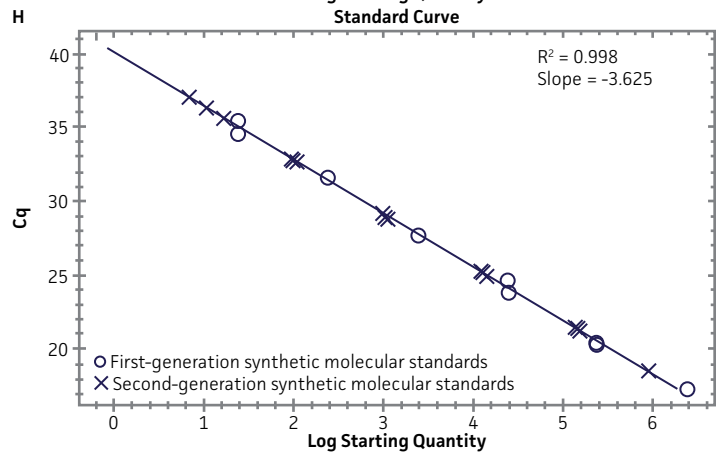
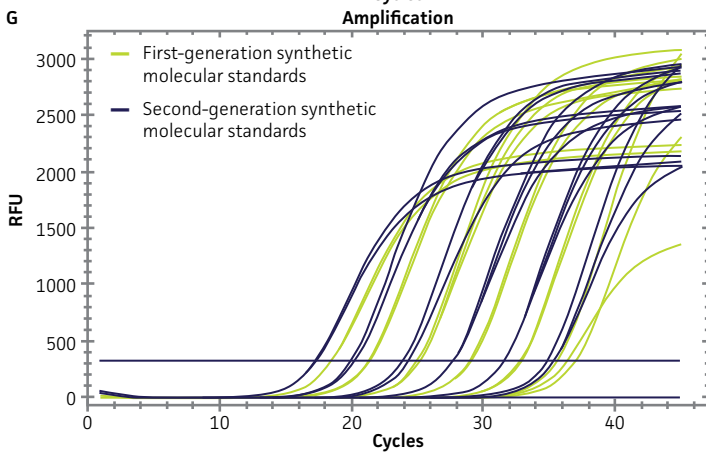
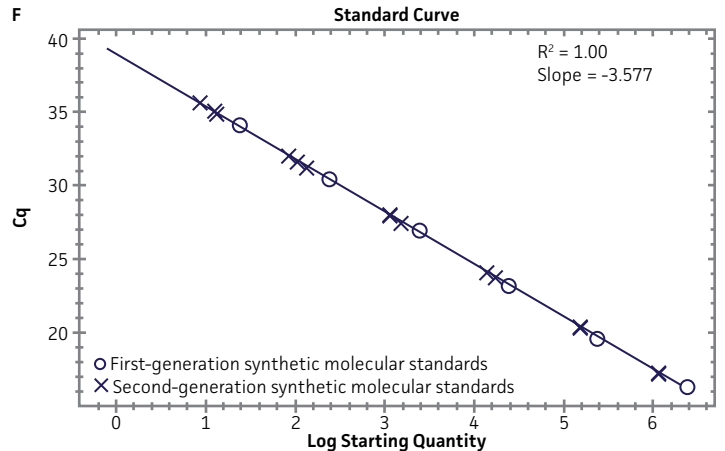
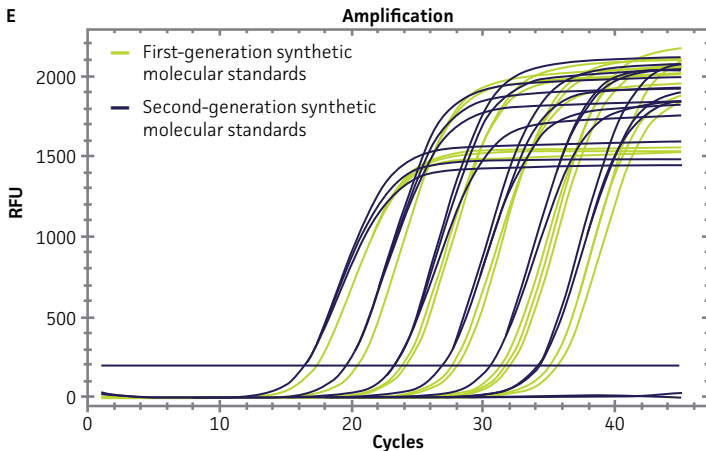
### COMPARISON OF FIRST-GENERATION AND SECOND-GENERATION SYNTHETIC MOLECULAR STANDARDS

First-generation and second-generation synthetic molecular standards were compared using the primer and probe sets from both the CaliciNet<sup>1</sup> and CEN/TC/WG6/TAG4<sup>2</sup> assays (Figure 2). Standard curves were generated using serial ten-fold dilutions of the synthetic RNA standards, and were tested in triplicate wells using the CFX96 Real-Time PCR Detection System as described above. From this analysis, it was determined that the first-generation molecular standards were not compatible with the primer and probe set used in the CEN/TC/WG6/TAG4 assay<sup>2</sup> for NoV-GI detection (Fig 2C), whereas the improvement in design of the newer second-generation constructs enabled consonance with both assays.

### NoV-GI



### NoV-GII

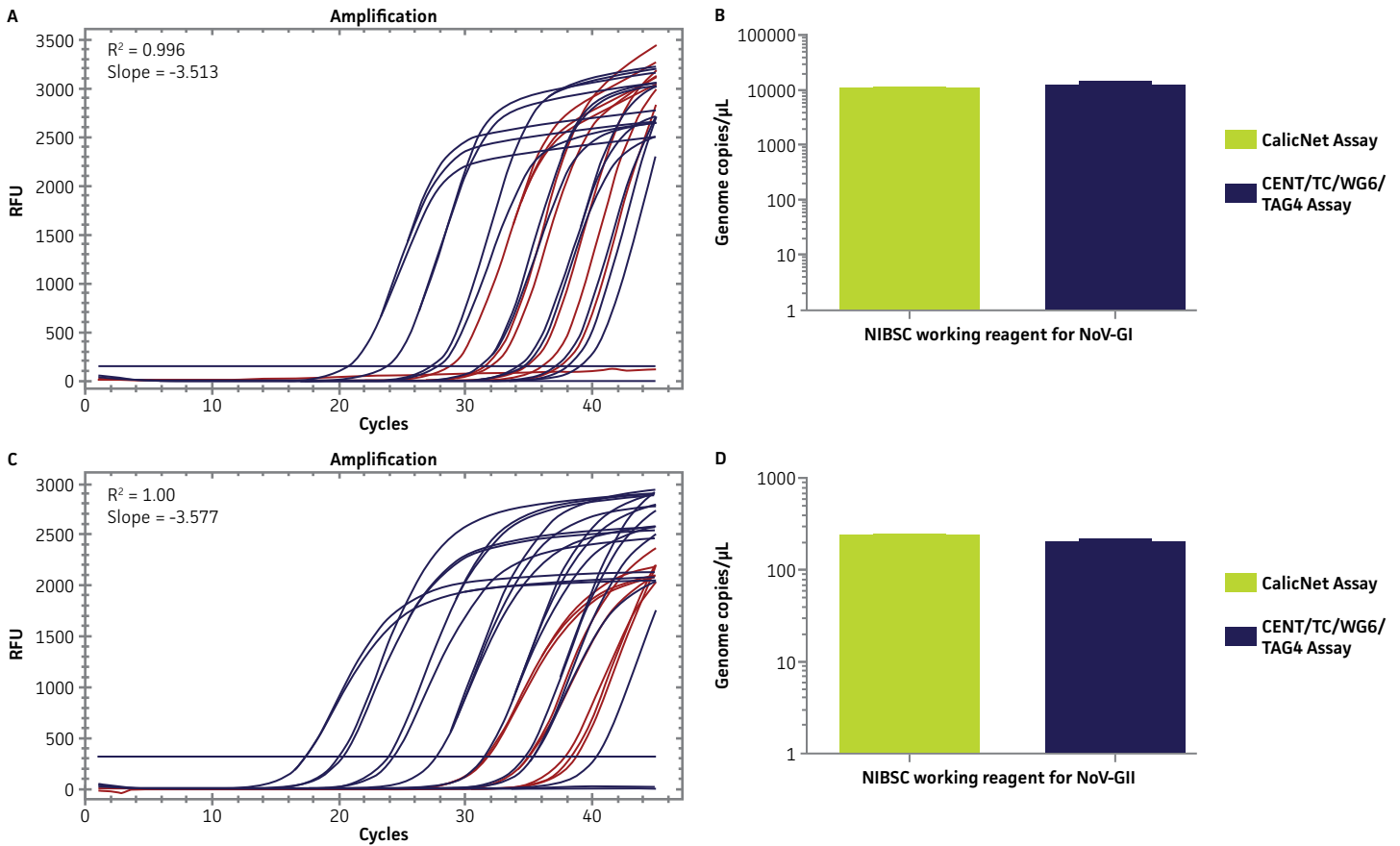


**Figure 2.** Comparison of First-Generation and Second-Generation Synthetic Molecular Standards

A&C) Amplification plots and B&D) standard curves generated using the first- and second-generation NoV-GI molecular standard ATCC® VR-3199SD and ATCC® VR-3234SD, respectively, versus the E&G) amplification plots and F&H) standard curves generated using the first- and second-generation NoV-GII synthetic molecular standards ATCC® VR-3200SD and ATCC® VR-3235SD, respectively. Amplification plots A&E) and standard curves B&F) were generated using the primer and probe set used in the CaliciNet assay<sup>1</sup>. Amplification plots C&G) and standard curves D&H) were generated using the primer and probe set used in the CEN/TC/WG6/TAG4 assay<sup>2</sup>.

### QUANTIFICATION OF THE NIBSC WORKING REAGENTS FOR NOV-GI AND NOV-GII USING ATCC SECOND-GENERATION SYNTHETIC MOLECULAR STANDARDS

As a proof of concept, we used the respective standard curves generated using the second-generation synthetic molecular standards paired with either the CaliciNet<sup>1</sup> or CEN/TC/WG6/TAG4 assay<sup>2</sup> primer and probe sets to quantify the NIBSC working reagents for NoV-GI and NoV-GII (Figure 3). Viral RNA from the NoV-GI and NoV-GII working reagents was extracted using the QIAamp® Viral RNA Mini Kit (QIAGEN®). The NIBSC working reagents were assayed in triplicate wells as undiluted samples and diluted 1:10 and 1:100 for the qRT-PCR assay. The average from the triplicate wells from undiluted samples was used to calculate the quantities of the NoV genome as  $2.3 \times 10^3$  copies / $\mu$ L for NoV-GI and 46.8 copies / $\mu$ L for NoV-GII.



**Figure 3.** Quantification of the NIBSC Working Reagents for NoV-GI and NoV-GII Using ATCC Second-Generation Synthetic Molecular Standards. An example of a qRT-PCR amplification plot showing the second-generation A) NoV-GI and C) NoV-GII synthetic molecular standards (Purple), and the working reagent from NIBSC (Red). B&D) Quantitation of the NIBSC working reagent for NoV-GI and NoV-GII samples, respectively, using the CaliciNet<sup>1</sup> and CEN/TC/WG6/TAG4 assays<sup>2</sup>.

### CONCLUSION

The second-generation NoV synthetic molecular standards have been significantly improved over the first-generation synthetic molecular standards. The quantitative nature of these standards directly allows for the generation of standard curves for determining NoV genome copies from unknown samples. Moreover, the improved design of the synthetic construct provides the ability to work with additional molecular assays for the detection and quantification of NoV.

### REFERENCES

1. Vega *et al.* Novel surveillance network for Norovirus gastroenteritis outbreaks, United States. *Emerg Infect Dis* 17(8):1389-1395, 2011.
2. Le Guyader *et al.* Detection and quantification of noroviruses in shellfish. *Appl Environ Microbiol* 75(3): 618-624, 2009.





10801 University Blvd.  
Manassas, VA 20110

AP-0316-02

© 2016 American Type Culture Collection. The ATCC trademark and trade name, any and all ATCC catalog numbers and any other trademarks listed in this publication are trademarks of the American Type Culture Collection unless indicated otherwise. Bio-Rad®, CFX96™, Droplet Digital™, and CFX Manager™ are registered trademarks or trademarks of Bio-Rad Laboratories, Inc. RNASTable® is registered trademark of Biomātrica, Inc. Stabilized by Biomātrica the biostability company. QIAamp® and QIAGEN® are registered trademarks of QIAGEN Inc.

These products are for laboratory use only. Not for human or diagnostic use. ATCC products may not be resold, modified for resale, used to provide commercial services or to manufacture commercial products without prior ATCC written approval.

**PHONE**

800.638.6597  
703.365.2700

**EMAIL**

[SalesRep@atcc.org](mailto:SalesRep@atcc.org)

**WEB**

[www.atcc.org](http://www.atcc.org)