

THE ESSENTIALS OF LIFE SCIENCE RESEARCH **GLOBALLY DELIVERED**[™]

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 390 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.

Native DENV RNA can be used as a standard for these assays; however, the full-length dengue viral RNA is on the Commerce Control List and requires a permit from the US Department of Commerce for international shipment. To make DENV RNA standards more accessible, ATCC has developed four synthetic molecular standards that represent DENV serotypes 1-4. Each standard contains short fragments from the capsid, membrane, and envelope genes of the DENV genome, as well as target regions encompassing the primer sequences from numerous published RT-PCR assays, including the DENV-1-4 Real-Time RT-PCR Assay developed by the CDC¹. The synthetic RNA standards were quantified by Droplet Digital[™] PCR (ddPCR[™]) in order to package precise copies of RNA. Moreover, given the inherent labile nature of RNA, a stabilization matrix was added to the quantitated RNA preparation. As compared to native RNA, these synthetic standards are easier to use as controls for qRT-PCR assays, exhibit less variability, have a longer shelf life, eliminate the need to culture viruses and can be used under BSL-1 conditions. Further, this synthetic quantitative RNA approach can be extended to other pathogenic viruses which are unculturable or need to be grown in a high-containment facility.

In the following proof-of-concept study, we amplified the synthetic molecular standards with the published primers from the CDC assay¹ and Waggoner *et al*², and used the generated standard curves to quantify viral RNA extracted from various DENV strains.

Materials and Methods

Viruses:

The following DENV strains, representing DENV serotypes 1, 2 and 4 were quantified using the standard curves generated from the DENV 1, 2 and 4 synthetic molecular standards, respectively:

Dengue virus 1 = TH-S-man (ATCC[®] VR-1586^m) Dengue virus 2 = New Guinea C (NR-84, BEI Resources) Dengue virus 4 = H241 (ATCC[®] VR-1257^m)

Viral RNA was extracted using the QIAamp[®] Viral RNA Mini Kit (QIAGEN[®]). Viral RNA was diluted 1:100, 1:1000, and 1:10,000 for the qRT-PCR assay. Uninfected Vero (ATCC[®] CCL-81[™]) and LLC-MK2 Derivative (ATCC[®] CCL-7.1[™]) cell lines were used as negative controls in the qRT-PCR assays.

qRT-PCR assay:

qRT-PCR assays were performed according to instructions provided¹, with slight modifications, using the CFX96[™] Real-Time PCR Detection System (Bio-Rad[®]). Two different primer sets were employed to examine each individual Dengue serotype using the same conditions for both: the commercial primer set included in the CDC Real-Time RT-PCR assay¹ and the primer set published by Waggoner et al.². 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 30 sec. The exception was for the DENV-4 published primer set; the annealing temperature was 55°C for 30 sec. Standard curves were generated using serial ten-fold dilutions of the synthetic RNA standards, ranging from 1 copy to 1 x 10⁶ copies/µL. RNA samples and standards were tested in triplicate. The relative fluorescence unit (RFU) baseline threshold was calculated using CFX Manager[™] 3.0 Software (Bio-Rad).

> ATCC 10801 University Boulevard, Manassas, Virginia 20110-2209

Development of Synthetic Molecular Standards for Dengue Virus

and Andrew G. Cawthon

ATCC Design Strategy



Advantages		
 ATCC fully authenticated & characterized Genuine Nucleics Consistent and accurate results BSL-1 ready-to-use control No shipping restrictions Quantitative format Stabilized RNA 	•	Gene quant Positi Indep and v Monit Iot va

DENV-4

Shamaila Ashraf, Melisa Wilson, Afshin Sohrabi, Stephen King, Brian Chase, Dev Mittar, Kurt Langenbach

Generation of Standard Curves from DENV Synthetic Molecular Standards



Primer and Probe		DENV-1	DENV-2	DENV-4
CDC Assay ¹	Slope	-3.244	-3.277	-3.642
	R ²	0.99	0.996	0.996
Published Assay ²	Slope	-3.536	-3.535	-3.775
	R ²	0.991	0.997	0.996

Figure 1. An example of A) amplification plot and B) standard curve generated using the DENV-4 molecular standard (ATCC[®] VR-3231SD^m) in conjunction with the CDC primer and probe set¹. C) The slope and R² values generated using DENV 1, 2 and 4 molecular standards with the primer and probe sets from the CDC DENV-1-4 Real-Time RT-PCR Assay¹ and Waggoner et al. assay².



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

VR-1586

References

- dengue viruses. PLoS Negl Trop Dis 7(4): e2116, 2013.

Disclaimers

Bio-Rad[®], CFX96[™], Droplet Digital[™]PCR, Real-Time PCR Detection System, and CFX Manager[™] 3.0 Software are registered trademarks or trademarks of Bio-Rad Laboratories, Inc. RNAstable[®] is registered trademark of Biomātrica, Inc. QIAamp[®] Viral RNA Mini Kit and QIAGEN[®] are registered trademarks of the QIAGEN Inc. The following reagents were obtained through BEI Resources, NIAID, NIH: Dengue Virus Type 2, New Guinea C, NR-

ATCC[®] VR-3231SD[™]

Applications

- eration of a standard curve for ititative RT-PCR (qRT-PCR) tive control for RT-PCR assays pendent standard for validation verification studies tor assay-to-assay and lot-to-
- riation



NR-84

VR-1257

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

1. CDC. DENV-1-4 Real-Time RT-PCR Assay for Detection and Serotype Identification of Dengue virus. Instructions for Use Package Insert, http://www.cdc.gov/dengue/resources/rt_pcr/CDCPackageInsert.pdf. 2. Waggoner JJ, et al. Single-reaction, multiplex, real-time RT-PCR for the detection, quantitation, and serotyping of