

Real-Time PCR Data Analysis Tool Utilizing a Novel Cloud-Based Software for Easy Interpretation of Animal Pathogen Detection



Denisse Meza, Rohan Shah, Robert Tebbs, Quoc Hoang, Rick Conrad – Thermo Fisher Scientific, Austin, TX, USA
 Anne Quijada, Emeline Ripoche, Sandrine Moine – Thermo Fisher Scientific, Lissieu, France

INTRODUCTION

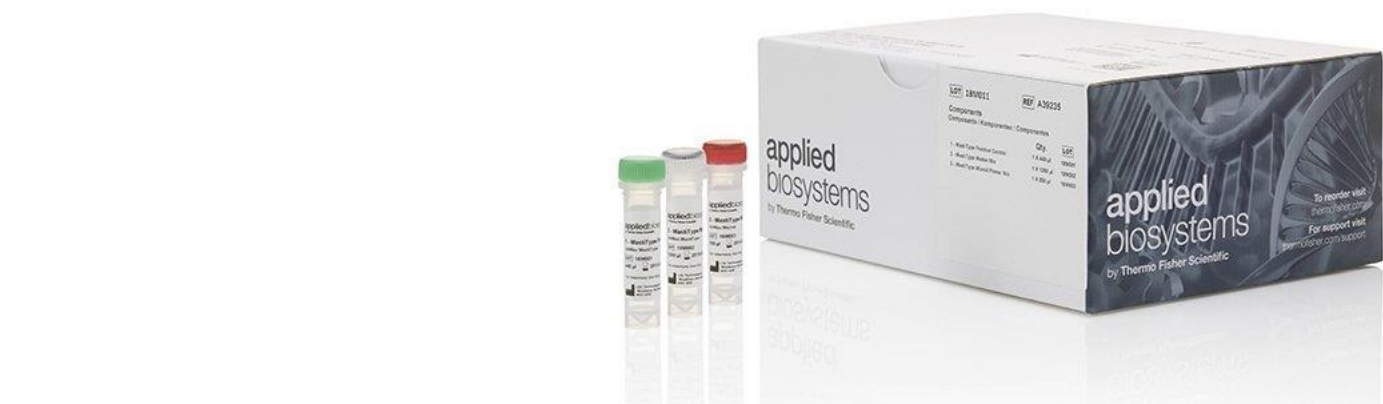
Current software versions on Real-Time PCR instruments were not specifically designed for the detection of pathogens, making data analysis and interpretation of multiplexed Animal Health assays difficult for users. Typically, qPCR data is analyzed manually which can be labor intensive and time-consuming. A complete package of data analysis software provides a faster and much more convenient alternative. The Animal Health group under the Applied Biosystems™ brand at Thermo Fisher Scientific now offers VeriVet Software, a user-friendly solution for the analysis of qPCR data for detecting animal pathogens.

MATERIALS AND METHODS

The solution is a new cloud-based software with a user-friendly graphical interface. The Animal Health VeriVet Software can be used to analyze data from runs performed on either the QuantStudio 5 or 7500 instrument series of Applied Biosystems Real-Time PCR Systems. Additionally, it has the capability to set up and remotely monitor a Real-Time PCR run in the QuantStudio 5 instrument.

1. Select assays for plate set up

The Animal Health VeriVet Software is compatible with all VetMAX kits and with Applied Biosystems 7500/7500 Fast and QuantStudio 5 Real-Time PCR Systems.



QuantStudio™ 5 Real-Time PCR System

7500 Fast & 7500 Real-Time PCR System

Administrator Settings

Assay List					
Add to Plate Setup	Assay Description	Assay Abbreviation	No of Targets	Assay Settings (Targets Cutoffs Thermal Protocol)	Assay Trends
<input type="checkbox"/>	VetMAX MastType Multi 4	Multi-4	5	View/Edit	View Trends
<input checked="" type="checkbox"/>	VetMAX MastType Micro4	Micro4	5	View/Edit	View Trends
<input checked="" type="checkbox"/>	VetMAX MastType Micro4 7500	Micro4 7500	5	View/Edit	View Trends
<input checked="" type="checkbox"/>	VetMAX MastType Multi 1	Multi-1	5	View/Edit	View Trends
<input type="checkbox"/>	VetMAX MastType Multi 1 7500	Multi-1 7500	5	View/Edit	View Trends
<input checked="" type="checkbox"/>	VetMAX MastType Multi 2	Multi-2	5	View/Edit	View Trends
<input type="checkbox"/>	VetMAX MastType Multi 2 7500	Multi-2 7500	5	View/Edit	View Trends
<input checked="" type="checkbox"/>	VetMAX MastType Multi 3	Multi-3	5	View/Edit	View Trends
<input type="checkbox"/>	VetMAX MastType Multi 3 7500	Multi-3 7500	5	View/Edit	View Trends
<input type="checkbox"/>	VetMAX MastType Multi 4 7500	Multi-4 7500	5	View/Edit	View Trends

Most VetMAX assays are pre-populated in the Assay List and new assays can be created easily.

2. Set up the plate layout, then run the Real-Time PCR

The plate layout is saved as a template file (EDT) in the Thermo Fisher Cloud.

- If using a cloud-connected instrument, the instrument run can be monitored from the app.
- If using an offline instrument, transfer the EDT file to the instrument, and then Import Results.

3. Review Results

Results are automatically analyzed and color-coded to produce calls on molecular testing determinations.

Well	Result	Sample	Assay	Target	Ct	Target Result	Custom Call	Task	Flag
A1	POSITIVE	Sample 9	VetMAX MastType Multi 1	M. bovis	23.53	POSITIVE	**	UNKNOWN	
				HC	26.52	POSITIVE	**	UNKNOWN	
				E. rhusiopathiae	22.23	POSITIVE	***	UNKNOWN	
				Enterococcus spp.	22.17	POSITIVE	***	UNKNOWN	
				C. bovis	23.13	POSITIVE	**	UNKNOWN	

RESULTS

The Animal Health VeriVet Software has been verified on a broad range of single and multiplexed assays containing 4 to 16 targets per run and provides equivalent results as manual analysis. The data analysis takes less than 5 minutes to set up, applies the recommended instrument settings, and provides a report with qualitative and quantitative results for each target.

Individual cells can be selected in either format to provide amplification curves.

CONCLUSIONS

Multiplexed Real-Time PCR assays yield a vast amount of data that can be cumbersome to analyze and interpret. The Thermo Fisher Animal Health VeriVet Software is designed to analyze these data rapidly with minimum user direction.

ACKNOWLEDGEMENTS

- Venkat Reddy
- Kalyana Rao Dulu
- Rohit Chaturvedi
- Somashree Patra
- Sarath Budam
- Shaohua Ma
- Deepali Gosain

