

Fast Optimization of a Multiplex Influenza Identification Panel Using a Thermal Gradient

Caine Leong,¹ Kim Pern Toh,² Sue Yuen Wee,³ Kuan Kiat Chew,³ and Masafumi Inoue²

¹Bio-Rad Laboratories (Singapore) Pte Ltd, Singapore; ²Experimental Therapeutic Centre (ETC), A*Star, Singapore; and ³Tan Tock Seng Hospital, Singapore

Introduction

The year 2009 was marked by the emergence of a novel influenza A (H1N1) virus that infects humans. There is a need to identify the different strains of influenza virus for purposes of monitoring the H1N1 strain pandemic and for other epidemiological and scientific purposes. In preparation for an accurate and rapid identification of H1N1, a laboratory in Singapore established a multiplex influenza identification panel that detects H1N1 among influenza A subtype H1 (flu A), subtype H3 (H3N2), and influenza B (flu B) viruses. The panel uses a 5' nuclease (TaqMan) probe-based four-target multiplex real-time PCR assay.¹

In this tech report we illustrate the use of the thermal gradient feature on the CFX96™ real-time PCR detection system to optimize reverse transcription (RT) and fast PCR conditions for a four-target multiplex PCR assay.

Methods

Viral RNA

The following virus strains were used: influenza A (H1N1), influenza A subtype H1 (flu A), influenza A subtype H3 (H3N2), and influenza B (flu B). Input concentration ranged between 10 and 1,000 copies.

Optimization of RT

Optimization of RT was performed with the four-target multiplex RT-qPCR MDX kit (Experimental Therapeutic Centre). The TaqMan probes were tagged with the following fluorophores: H1N1, HEX; flu A, FAM; H3N2, Cy5; and flu B, Texas Red.

RT-qPCR reactions were set up as indicated in Table 1 using the iScript™ one-step RT-PCR kit for probes (Bio-Rad Laboratories, Inc.).² Reactions were incubated in the CFX96 system (Bio-Rad) for 30 min at 45, 45.6, 47, 48.9, 51.3, 53.3, 54.4, and 55°C (temperature gradient), followed by 2.5 min at 95°C, then subjected to 42 three-step cycles of 95°C for 17 sec, 55°C for 31 sec, and 68°C for 32 sec. Fluorescence data were collected at the 55°C annealing step. Positive control RNAs equivalent to approximately 1,000 viral copies were used for optimization.

Multiplex RT-qPCR

RT-qPCR reactions were set up as indicated in Table 1. Reactions were incubated at 50°C for 30 min followed by 2.5 min at 95°C, then subjected to 42 three-step cycles of 95°C for 17 sec, 55°C for 31 sec, and 68°C for 32 sec. Fluorescence data were collected at the 55°C annealing step.

Table 1. Reaction setup for ten reactions.

Components	Volume, μ l
2x RT-qPCR reaction buffer	125.0
H1N1 primer pair mix (7.5 μ M each)	5.0
H1N1 probe-HEX (10 μ M)	2.5
Flu A primer pair mix (7.5 μ M each)	5.0
Flu A probe-FAM (10 μ M)	2.5
H3N2 primer pair mix (7.5 μ M each)	5.0
H3N2 probe-Cy5 (10 μ M)	2.5
Flu B primer pair mix (7.5 μ M each)	5.0
Flu B probe-Texas Red (10 μ M)	2.5
RNA template (range from 0.01 pg–1 μ g)	Up to 90
Reverse transcriptase and Taq polymerase mix	5.0
RNase-free H ₂ O	Top up to 250

Optimization of Fast qPCR

RT-qPCR reactions were set up as indicated in Table 1, incubated for 10 min at 50°C followed by 2.5 min at 95°C, then subjected to 42 two-step cycles of 95°C for 17 sec and 31 sec at 55, 56, 58, 60.9, 64.5, 67.5, 69.2, and 70°C (temperature gradient). Fluorescence data were collected at the gradient range of the 55–70°C annealing-extension step. Positive control RNAs equivalent to about 1,000 viral copies were used for fast qPCR optimization.

Results and Discussion

RT Optimization Using the Thermal Gradient

The temperature gradient feature of the CFX96 real-time PCR system was used to determine the optimal RT temperature for the four-target multiplex RT-qPCR assay. The CFX96 system was programmed to perform RT at temperatures ranging from 45 to 55°C, followed by a three-step PCR protocol (denaturation-annealing-extension). Results from the RT gradient in Table 2 show that the optimal RT temperature range (in blue) was between 45.0 and 51.3°C for all primer-probe sets, with influenza A reactions working as high as 53.3°C. Selecting higher RT temperatures improves the specificity of the assay but compromises the assay sensitivity. In order to maintain both specificity and sensitivity, we have selected 50°C RT for the rest of the four-target multiplex assay validation.

Table 2. Optimization of RT reaction temperature for the four-target multiplex influenza RT-qPCR assay.

RT Reaction Temperature, °C	C_T^*			
	H1N1 (HEX)	Flu A (FAM)	H3N2 (Cy5)	Flu B (Texas Red)
55.0	32.68	33.70	32.41	31.71
54.4	32.34	33.70	32.46	30.67
53.3	32.09	32.50	31.81	30.57
51.3	31.67	32.48	31.70	29.26
48.9	31.45	32.66	31.04	28.89
47.0	31.18	32.02	31.09	29.45
45.6	31.16	32.24	31.15	29.47
45.0	30.93	32.97	30.98	29.43

* Threshold cycle (C_T) values in blue reflect optimal RT temperatures yielding the lowest $C_T \pm 0.5^\circ\text{C}$ for each primer-probe set, while C_T values in red indicate suboptimal temperatures.

Multiplex RT-qPCR Assay Sensitivity

To determine the multiplex RT-qPCR assay sensitivity, we tested different concentrations of the four viral RNAs ranging from 1,000 to 10 copies. Results shown in Figure 1 clearly show that the primer-probe sets were able to detect all the viral RNA controls with no cross talk between channels.

Optimization of Fast qPCR Conditions

Optimization of fast qPCR conditions was achieved by combining the annealing and extension steps. Combining these steps is essential for 5' nuclease (TaqMan) probe-based assays because the probes need to anneal to the target while the product is being extended.

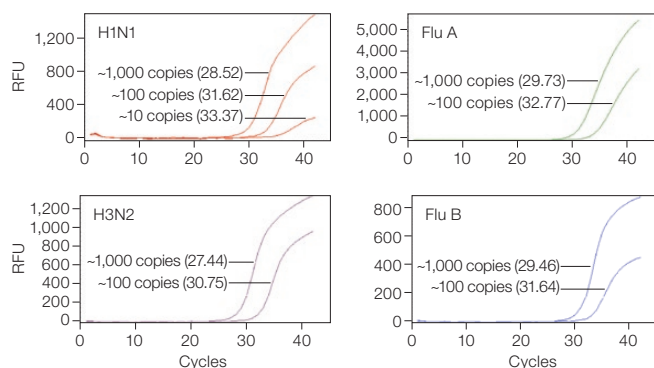


Fig. 1. Multiplexing of four-target influenza RT-qPCR. Amplification profiles of H1N1, flu A, flu B, and H3N2. Viral RNA isolates ranging from 1,000 to 10 copies were detected. RFU, relative fluorescence units.

Additional timesavings were achieved by reducing the RT time from 30 min to 10 min. C_T values in Table 3 and amplification plots in Figure 2 illustrate the optimal fast qPCR annealing-extension temperature range for H1N1 (58–60.9°C), flu A (60.9–64.5°C), flu B (55–58°C), and H3N2 (58–64.5°C). We selected 58°C for annealing-extension to perform the fast four-target multiplex RT-qPCR assay validation (total run time for this reaction was shortened from 128 to 73 min). The results generated using fast cycling conditions in Table 4 show that all viral RNA controls tested were detected with a 1–2 C_T delay compared to the standard 30 min RT using three-step qPCR cycling conditions.

Table 3. Optimization of the four-target multiplex influenza RT-qPCR assay for fast PCR.

Annealing Temperature, °C	C_T^*			
	H1N1 (HEX)	Flu A (FAM)	H3N2 (Cy5)	Flu B (Texas Red)
70.0	—	—	—	—
69.2	—	41.98	—	—
67.5	—	31.11	—	—
64.5	—	29.47	27.67	—
60.9	30.81	29.79	27.55	31.68
58.0	31.26	30.01	27.78	30.54
56.0	31.83	30.15	28.18	30.54
55.0	31.53	30.39	28.44	30.06

* C_T values in blue reflect optimal annealing-extension temperatures yielding the lowest $C_T \pm 0.5^\circ\text{C}$ for each primer-probe set, while C_T values in red indicate suboptimal temperatures.

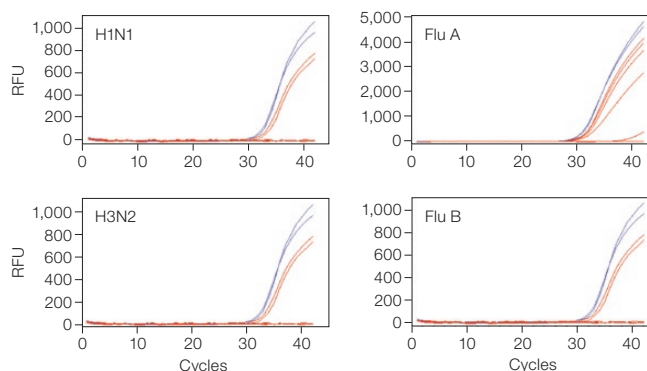


Fig. 2. Optimization of fast qPCR for the four-target multiplex influenza RT-qPCR. Amplification profiles of PCR reactions listed in Table 3. Blue and red traces indicate PCR reactions performed at optimal and suboptimal annealing temperatures, respectively. RFU, relative fluorescence units.

Table 4. Comparison of standard PCR and fast PCR efficiency.

Viral RNA	Copy Number	C_T		ΔC_T
		Standard Run (128 min)	Fast Run (73 min)	
H1N1	1,000	28.52	27.85	-0.67
	100	31.62	32.32	0.70
	10	33.37	34.42	1.05
Flu A	1,000	29.73	30.99	1.26
	100	32.77	34.32	1.55
	10	33.37	34.42	1.05
H3N2	1,000	27.44	28.72	1.28
	100	30.75	32.38	1.63
	10	33.37	34.42	1.05
Flu B	1,000	29.46	31.45	1.99
	100	31.64	33.54	1.90
	10	33.37	34.42	1.05

Conclusions

We have demonstrated that the gradient feature of the CFX96 system dramatically accelerates the optimization of the RT and PCR steps for multiplexing RT-qPCR assays. The thermal gradient also improved the ease of converting cycling conditions from a three-step to a two-step PCR assay, thus significantly reducing PCR run times while maintaining assay specificity and sensitivity.

¹The test protocol described in this paper has not been cleared or approved by the FDA or any other regulatory agency for human diagnostic or other clinical use.

²This product may be used for scientific research purposes only. Please refer to the product insert for additional information.

Cy is a trademark of GE Healthcare group companies. FAM is a trademark of Applied Biosystems Corporation. TaqMan is a trademark of Roche Molecular Systems, Inc. Texas Red is a trademark of Invitrogen Corporation.

Purchase of iScript one-step RT-PCR kit for probes includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied Biosystems' United States Patent Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

Information in this tech note was current as of the date of writing (2009) and not necessarily the date this version (Rev A, 2010) was published.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000
