

Alien Reference RNA QRT-PCR Detection Kit for Monitoring the Overall Performance of QRT-PCR Assays

Bahram Arezi, Melissa McCarthy & Holly Hogrefe
Agilent Technologies, Stratagene Products Division, La Jolla, CA 92037

Introduction

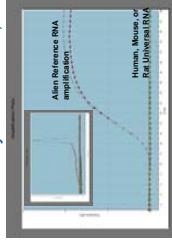
Exogenous RNA controls are increasingly used in QRT-PCR experiments to normalize variations that arise due to the presence of co-purified inhibitors (1,2). In addition, RNA controls allow assay standardization from experiment-to-experiment, across platforms, and between laboratories (3). We previously described the use of a highly sensitive exogenous Alien RNA transcript (Alien QRT-PCR Inhibitor Alert, cat. no. 300600) for monitoring the presence of common QRT-PCR inhibitors in RNA samples. The Inhibitor Alert kit also includes a sample RNA versus Alien RNA assay. This Inhibitor Alert kit employs SYBR Green chemistry, which provides high sensitivity, but does not allow multiplex amplification of Alien RNA alongside a transcript of interest.

We have recently expanded our line of exogenous controls to include TaqMan primers and probe (FAM- and VIC-labeled) sets for detection of exogenous Alien RNA transcript in 1-step and 2-step QRT-PCR assays. As we will demonstrate, the Alien probe and primer sets are highly sensitive and specific for detection of Alien RNA. In addition, the Alien probe is sensitive to a variety of QRT-PCR inhibitors, and allow multiplex detection of Alien and sample RNA in the same tube.

Results

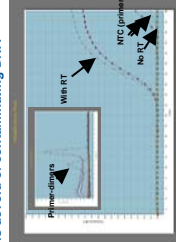
Alien Reference RNA template is a 500bp *in vitro* transcript made from a computer generated sequence that possesses 50% GC content and is non-homologous to sequences in GenBank.

Part 1: Alien Reference RNA does not cross hybridize to human, mouse, or rat sequences



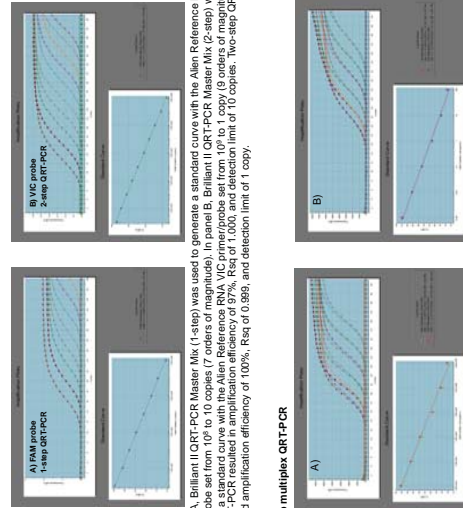
Amplification reactions contained 100nM of Alien RNA primers and either 10⁷ copies of Alien RNA target or 50ng of Human, Mouse, or Rat Universal Reference RNA (Stratagene) in separate tubes. Amplifications were performed with the Brilliant[®] SYBR[®] Green QRT-PCR Master Mix (1-step). No amplification products were detected using Human, Mouse, or Rat Universal Reference RNA.

Part 2: Alien Reference RNA is devoid of contaminating DNA



Amplification reactions were performed using Brilliant[®] SYBR[®] Green QRT-PCR Master Mix (1-step), 10⁷ copies of Alien Reference RNA, and 100nM of Alien RNA primers with and without the reverse transcriptase as indicated. No amplification is detected in the no-RT control.

Part 3: Alien Reference RNA primer/probe sets exhibit high amplification efficiency, sensitivity, and dynamic range



In panel A, Brilliant II QRT-PCR Master Mix (1-step) was used to generate a standard curve with the Alien Reference RNA FAM primer/probe set from 10⁷ to 10 copies (7 orders of magnitude). In panel B, Brilliant II QRT-PCR Master Mix (2-step) was used to generate a standard curve with the Alien Reference RNA VIC primer/probe set from 10⁷ to 1 copy (9 orders of magnitude). One-step QRT-PCR resulted in amplification efficiency of 97%, Rsq of 1.000, and detection limit of 10 copies. Two-step QRT-PCR generated amplification efficiency of 100%, Rsq of 0.999, and detection limit of 1 copy.

b) 2-step multiplex QRT-PCR

Results-continued

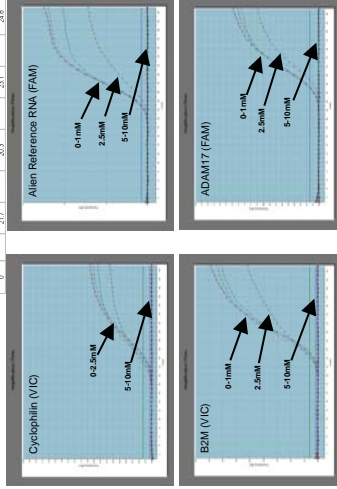
Brilliant Multiplex QPCR Master Mix was used to generate standard curves for cyclophilin (VIC) and Alien Reference RNA (FAM) individually or in multiplex. Amplifications were performed using Human QPCR Reference RNA ranging from 100 ng to 0.001 ng, and Alien Reference RNA ranging from 10⁷ to 10² copies. Panel A shows cyclophilin standard curves in the absence (SAF = Single reaction or presence (SAF = Duplex reaction) of Alien Reference RNA amplification. Panel B shows Alien Reference RNA standard curves in the absence (SAF = Single reaction) or presence (SAF = Duplex reaction) of cyclophilin amplification.

Part 4: Alien Reference RNA amplification is sensitive to the presence of common QRT-PCR inhibitors

In order to detect inhibition, amplification of Alien Reference RNA alone (singleplex) is compared to the amplification of Alien Reference plus sample RNA (in singleplex or multiplex format). A delay in Ct of Alien Reference plus sample RNA suggests that the RNA sample contains inhibitors. To mimic this situation, we added two common RT-PCR inhibitors (EDTA and ethanol) to Alien Reference plus sample RNA amplification reactions. In order to determine the Ct for 'no inhibitors' in sample RNA amplification reactions, we determined the Ct for 'no inhibitors' in sample RNA amplification reactions using standard multiplex probe formats, and hence should be able to detect a wide range of 'unknown' inhibitors in RNA samples.

a) EDTA in one-step singleplex QRT-PCR:

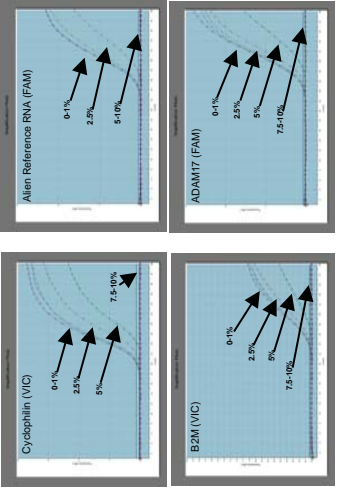
EDTA (nM)	Alien		Cyclophilin		B2M		ADAM17	
	Nt	Nt	Nt	Nt	Nt	Nt	Nt	Nt
100	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
5	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
0.5	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
0.005	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7



Amplifications were performed using Brilliant II one-step QRT-PCR kit, 10⁷ copies of Alien Reference RNA and 10ng Human Reference RNA were added to the reactions. EDTA was titrated into the reactions at the concentrations indicated above. Inhibition of amplification of Alien Reference RNA and ADAM17 (low abundance target) was detected at EDTA concentrations of 2.5nM with Ct delays of 1.7 (Alien Reference RNA) and 2.5 (ADAM17) cycles. No inhibition was observed at 0.1nM EDTA. At 5nM EDTA, amplification of B2M showed a slight reduction in final fluorescence but no Ct delay at 2.5nM EDTA. Concentrations of 2.5nM EDTA inhibited all target amplifications. Therefore, Alien Reference RNA and ADAM17 are more sensitive to the inhibitory effects of EDTA compared to B2M and cyclophilin.

b) Ethanol in one-step singleplex QRT-PCR:

Ethanol (%)	Alien		Cyclophilin		B2M		ADAM17	
	Nt	Nt	Nt	Nt	Nt	Nt	Nt	Nt
10	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
5	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
2.5	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
1	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
0.5	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
0.25	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7
0.1	5.5	21.7	5.5	21.7	5.5	21.7	5.5	21.7

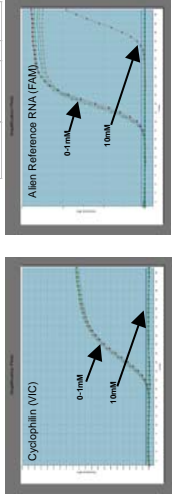


Amplifications were performed using Brilliant II one-step QRT-PCR kit, 10⁷ copies of Alien Reference RNA and 10ng Human Reference RNA were added to the reactions. Ethanol was titrated into the reactions at the concentrations indicated above. When 2.5% ethanol was added into the reaction, amplifications of Alien Reference RNA, cyclophilin, B2M, and ADAM17 targets were delayed by 2.5, 1.3, 0.7, and 0.2 Cts, respectively. At 5% ethanol, Alien Reference RNA failed to amplify whereas all other targets produced delayed Cts, suggesting that Alien Reference RNA is more sensitive to the inhibitory effects of ethanol compared to all the targets tested here.

Results-continued

c) EDTA in two-step singleplex QRT-PCR:

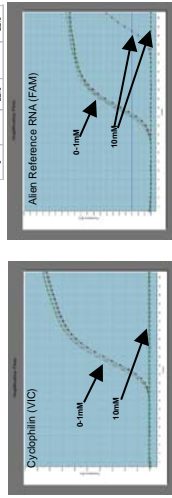
EDTA (nM)	Cyclophilin		Alien	
	Nt	Nt	Nt	Nt
100	17.37	17.37	17.37	17.37
10	17.37	17.37	17.37	17.37
1	17.37	17.37	17.37	17.37
0.1	17.37	17.37	17.37	17.37



Amplifications were performed using Brilliant II two-step QRT-PCR kit, 10⁷ copies of Alien Reference RNA and 10ng Human Reference RNA were added to the cDNA synthesis reactions. EDTA was titrated into the reactions at the concentrations indicated above. Inhibition of amplification of both Alien Reference RNA and cyclophilin was observed at 2.10nM EDTA.

d) EDTA in two-step multiplex QRT-PCR:

EDTA (nM)	Cyclophilin		Alien	
	Nt	Nt	Nt	Nt
1	22.1	22.1	22.1	22.1
0.5	22.1	22.1	22.1	22.1
0.005	22.1	22.1	22.1	22.1
0	22.1	22.1	22.1	22.1



Amplifications were performed using Brilliant Multiplex QPCR kit, 10⁷ copies of Alien Reference RNA and 10ng Human Reference RNA were added to the reactions. EDTA was titrated into the reactions at the concentrations indicated above. Inhibition of both Alien Reference RNA and cyclophilin amplifications are completely inhibited at 10nM EDTA similar to the results obtained above (part 4, section c).

Conclusions

- Alien Reference RNA QRT-PCR detection kit is used for validating the quality of experimental RNA samples, interpreting the quality of QRT-PCR data, and monitoring the overall performance of QRT-PCR assay reagents and instrumentation.
- The inclusion of Alien Reference RNA into long-term experimental studies, ensures the availability of a consistent quality of RNA for all experiments.
- The inclusion of Alien Reference RNA into multiplex assays ensures the availability of a consistent quality of RNA for all experiments.
- Multiplexing Alien Reference RNA with a gene-of-interest (GOI) general multiplexing guidelines should be followed to minimize the inhibition of the GOI. This includes the use of a multiplex QPCR kit (e.g., Brilliant Multiplex kit), and limiting primer/probe concentration for higher abundance targets.
- If the presence of an inhibitor in RNA samples is suspected, researchers are encouraged to re-purify or dilute the RNA sample (if target abundance allows it) and repeat the amplification. Furthermore, a two-step QRT-PCR format may be more appropriate to use due to increased tolerance of QRT-PCR inhibitors.
- Alien Reference RNA QRT-PCR detection kit will be commercially available on 11/2008.

References

1. Smith, R.D. et al. *Exogenous Reference RNA for Normalization of Real-time Quantitative PCR*. Biotechniques, 2003, 34:p 88-91.
2. Arezi, B. et al. *SPUD-A quantitative PCR assay for the detection of inhibitors in nucleic acid preparations*, 2006, 35:1-p 309-10.
3. Hartman, L.J. et al. *Development of a novel internal positive control for TaqMan based assays*. Molecular and Cellular Probes, 2005, 19: p 54-59.

For Research Use Only. Not for use in diagnostic procedures.