automated electrophoresis

Myth Busted: A NanoDrop ND-1000 Spectrophotometric Reading Is Insufficient to Assess RNA Quality

Introduction

According to the recently published quantitative real-time PCR experiment publication guidelines, or MIQE (Bustin et al. 2009; http://www.rdml.org/miqe.php), providing RNA integrity data is essential when publishing real-time quantitative PCR (qPCR) and reverse transcription qPCR (RT-qPCR) data. RNA integrity is critical in qPCR and RT-qPCR experiments in order to obtain gene expression results that are reliable and reproducible, and therefore publishable.

The Experion[™] automated electrophoresis system (Bio-Rad Laboratories, Inc.) provides an automatic assessment of RNA integrity by providing the RNA quality indicator (RQI) in addition to the electropherogram, gel view, and 28S/18S ratio and concentration (Figure 1, Table 1). Here we illustrate that the NanoDrop (ND-1000) spectrophotometer (Thermo Fisher Scientific, Inc.) can assess concentration and also provide some purity data (via A_{260/280} nm and A_{260/230} nm readings) for the same mouse liver total RNA sample that has been degraded to varying degrees (Table 1). However, the ND-1000 spectrophotometer does not provide RNA integrity data.

When mouse liver samples (with varying degrees of integrity, as generated by heating samples at 90°C) were analyzed on both the Experion and ND-1000 systems, Experion system results show:

- Samples 1, 2, and 3 (RQI 9.8, 9.2, and 8.1, respectively) are highly intact and can be used in downstream applications with confidence
- Samples 4 and 5 (RQI 6.5 and 5.9, respectively) are somewhat degraded and may or may not be usable depending on the application
- Samples 6, 7, and 8 are highly degraded (RQI 2.2, 2.0, and 1.8, respectively) and should not be used for downstream applications (Figure 2)

ND-1000 spectrophometer results indicate that all samples are good quality.



Fig. 1. Experion system electropherogram overlay of mouse liver total RNA sample progressively degraded with heat. With longer heat exposure, there is an observed decrease in the 28S and 18S peaks with degradation peaks appearing between the lower marker and the 18S peak.

Table 1. Experion system and ND-1000 spectrophotometer assessment of the same mouse liver total RNA sample. Experion system results give fast (11–12 samples in 30 min) and the most complete information (electropherogram and virtual gel profile, RQI value, and 28S/18S ratio) to help determine RNA integrity. (
), very little or no degradation; (
), some degradation; (
), significant degradation.

		Experion Automated Electrophoresis System			NanoDrop ND-1000 Spectrophotometer		
Sample Number	Sample Name	Ratio 28S/18S	RQI	RQI Class	Conc, ng/µl	A _{260/280} *	A_260/230 *
1	Control — no heat	1.60	9.8		115	1.90	2.44
2	3 min @ 90°C	1.23	9.2		114	1.93	2.40
3	5 min @ 90°C	0.89	8.1		115	2.06	2.37
4	10 min @ 90°C	0.50	6.5		115	2.03	2.37
5	15 min @ 90°C	0.15	5.9		116	2.02	2.31
6	1.0 hr @ 90°C	0.46	2.2		109	1.99	2.18
7	2.0 hr @ 90°C	0.81	2.0) 117 (2.00	2.32
8	4.0 hr @ 90°C	0.00	1.8	•	118	1.89	2.23

* Note: Generally accepted ratios (A_{260/280} and A_{260/230}) for good quality RNA are >1.8; the ND-1000 spectrophotometric readings indicate that all samples are good quality.



Fig. 2. Experion system virtual gel clearly shows that samples 6, 7, and 8 are highly degraded.



Conclusions

NanoDrop ND-1000 spectrophotometric readings provide only part of the information needed for reliable RT-qPCR results. Conversely, the Experion system provides a complete evaluation of total RNA, with all results automatically generated, saved, and easily exported.

Reference

Bustin SA et al. (2009). The MIQE guidelines: minimum information for publication of quantitative real-time PCR experiments. Clin Chem 55:4, 611-622.

NanoDrop is a trademark of Thermo Fisher Scientific, Inc.



Bio-Rad Laboratories, Inc.

Life Science Group Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 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