

### Automated Purification of RNA from Strawberry

*Purify RNA from strawberry fruit using the Maxwell® RSC Instrument and the Maxwell® RSC Plant RNA Kit.*

<b>Kit:</b>	Maxwell® RSC Plant RNA Kit (Cat.# AS1500)
<b>Analyses:</b>	Dye-based quantitation and qRT-PCR
<b>Sample Type(s):</b>	Strawberry ( <i>Fragaria x ananassa</i> ) fruit
<b>Input:</b>	50-100mg of strawberry fruit
<b>Materials Required:</b>	<ul style="list-style-type: none"><li>▪ Maxwell® RSC Plant RNA Kit (Cat.# AS1500)</li><li>▪ CTAB buffer (Cat.# MC1411)</li><li>▪ Maxwell® RSC Instrument (Cat.# AS4500)</li><li>▪ Microcentrifuge</li><li>▪ Heat block capable of 65°C</li></ul>

This protocol was developed by Promega Applications Scientists and is intended for research use only.

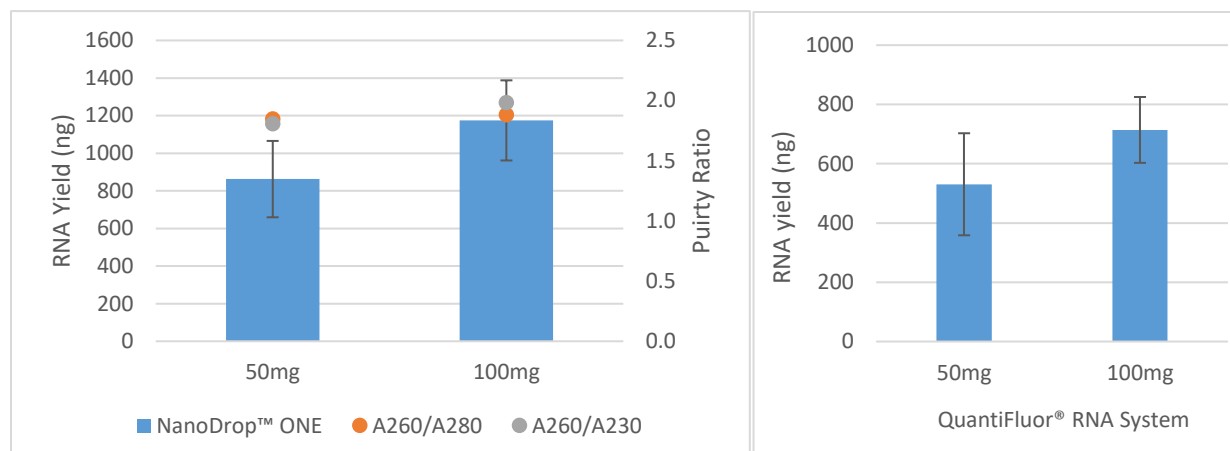
Users are responsible for determining suitability of the protocol for their application.

For further information, see Technical Manual TM459, available at:  
**[www.promega.com/protocols](http://www.promega.com/protocols)**  
or contact Technical Services at:  
**[techserv@promega.com](mailto:techserv@promega.com)**

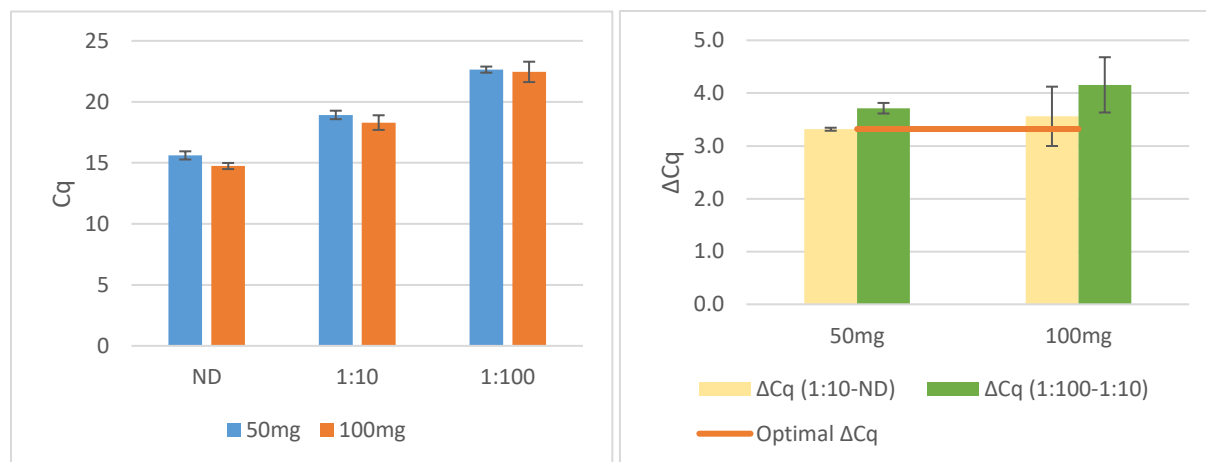
#### Protocol:

1. Weigh out 50-100mg of frozen strawberry fruit pieces and add 600µl of CTAB buffer supplemented with 2% 1-thioglycerol.
2. Vortex vigorously at maximum speed for 10 seconds to mix.
3. Place samples in a standard heat block at 65°C for 15 minutes.
4. Centrifuge for 5 minutes at 15,000 x *g* to separate any solid.
5. Transfer 400µl of the cleared sample supernatant to a clean 1.5 microcentrifuge tube.
6. Add 200µl of Lysis Buffer to the cleared supernatant and mix by vortexing at maximum speed for 10 seconds.
7. Incubate at room temperature for 10 minutes and transfer the lysate into well #1 of the Maxwell® RSC cartridge.
8. Add 5µl of DNase I solution to well #4.
9. Add plungers to well #8 of the Maxwell® RSC cartridge.
10. Place the supplied elution tubes into the sample rack and add 50µl of the Nuclease-Free Water.
11. Run the method *Maxwell® RSC Plant RNA Kit* on the Maxwell® RSC Instrument.

## Results:



**Figure 1. Quantification of RNA isolated from frozen strawberry fruit.** RNA was isolated from 50mg or 100mg of frozen strawberry fruit using the Maxwell® RSC Plant RNA Kit (Cat.# AS1500) modified with CTAB buffer (Cat.# MC1411) as the homogenization buffer. The RNA yield and purity ratio were quantified using both **(Left)** NanoDrop™ One and **(Right)** QuantiFluor® RNA System (Cat.# E3310). Average values  $\pm$  STD from three experiments are shown.



**Figure 2. qPCR amplification of RNA isolated from strawberry fruit.** Strawberry RNA was detected using GoTaq® 1-Step RT-qPCR System (Cat.# A6020). Cq value of different amount of input materials was shown. Undiluted RNA eluate prepared from both 50mg and 100mg of strawberry fruit was diluted at 1:10 and 1:100 and amplified with strawberry specific primers<sup>1</sup> to examine the amplification efficiency. No inhibitors were found to be co-purified with the RNA according to the  $\Delta$ Cq analysis. Average values from three experiments  $\pm$  STD are shown.

## References:

1. Amil-Rutz et al (2013) Identification and Validation of Reference Genes for Transcript Normalization in Strawberry (*Fragaria x ananassa*) Defense Responses. PLoS ONE (8): e70603.