

## DNA extraction from *Saccharomyces cerevisiae* using Maxwell® RSC instrument

*DNA extraction from Saccharomyces cerevisiae using Maxwell® RSC whole blood DNA kit*

<b>Kit:</b>	Maxwell® RSC Whole Blood DNA kit (Cat. #AS1520)
<b>Analyses:</b>	Nanovue® quantification and qPCR amplification
<b>Sample Type(s):</b>	<i>S. cerevisiae</i> pellet
<b>Input:</b>	up to $9 \times 10^7$ cells
<b>Materials Required:</b>	<ul style="list-style-type: none"><li>▪ Maxwell® RSC Whole Blood DNA kit (Cat. #AS1520)</li><li>▪ Maxwell® RSC Instrument (Cat. #AS4500)</li><li>▪ Lyticase (Sigma, Ref.L2524)</li><li>▪ TE Buffer, 1X, Molecular Biology Grade (Cat. #V6231)</li></ul>

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

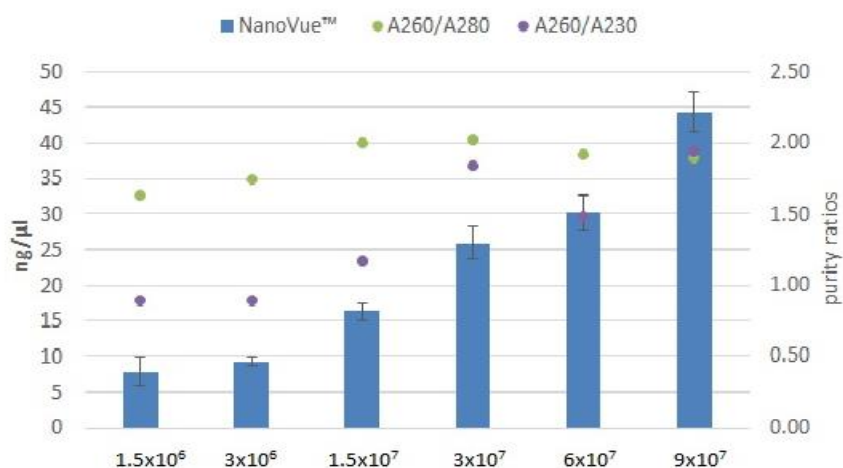
Users are responsible for determining suitability of the protocol for their application.

Further information can be found in Technical Manual #TM455, available at:  
[www.promega.com/protocols](http://www.promega.com/protocols)

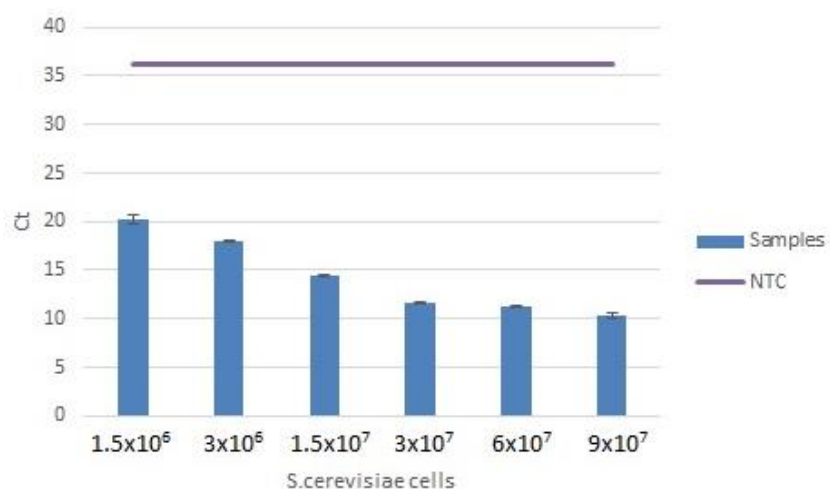
### Protocol:

1. Make pellets of a *S. cerevisiae* culture with a O.D~1 (approx.  $3 \times 10^7$  cells)
2. Add 12.5 µl of lyticase at 4 u/µl + 87.5 µl of TE Buffer 1X and resuspend the pellet
3. Incubate for 60 min at 30°C and 1000 rpm
4. Add the lysate on well #1 of the cartridge
5. Add a plunger on well #8 of the cartridge
6. Place an elution tube into the elution tube position and add 60µl of elution buffer
7. Run the whole Blood method on the Maxwell® RSC instrument

## Results:



**Figure 1. DNA concentration and purity ratios.** DNA extracted with Maxwell® RSC Instrument from 1.5x10<sup>6</sup>, 3x10<sup>6</sup>, 1.5x10<sup>7</sup>, 3x10<sup>7</sup>, 6x10<sup>7</sup> and 9x10<sup>7</sup> *S.cerevisiae* cells pellets were quantified and ratios were measured using NanoVue™ spectrophotometer. N=3



**Figure 2. qPCR amplification.** Ct values obtained for 2 μl of DNA extracted from *S.cerevisiae* using GoTaq® qPCR master mix and 18Sr primers<sup>1</sup>. NTC= 36.2. N=3±SD

## Reference:

<sup>1</sup> Cabib, E., Roh, D.H., Schmidt, M., Crotti, L.B. & Varma, A. The yeast cell wall and septum as paradigms of cell growth and morphogenesis. J Biol Chem 276, 19679-19682 (2001).