

# Quantify chlorophyll *a* and chlorophyll *b* with a custom method

## Using the NanoDrop One Spectrophotometer

### Abstract

Scientists can accurately quantify chlorophyll *a* and chlorophyll *b* on the Thermo Scientific™ NanoDrop™ One/One<sup>c</sup> Microvolume UV-Vis Spectrophotometer using a user-defined custom method.

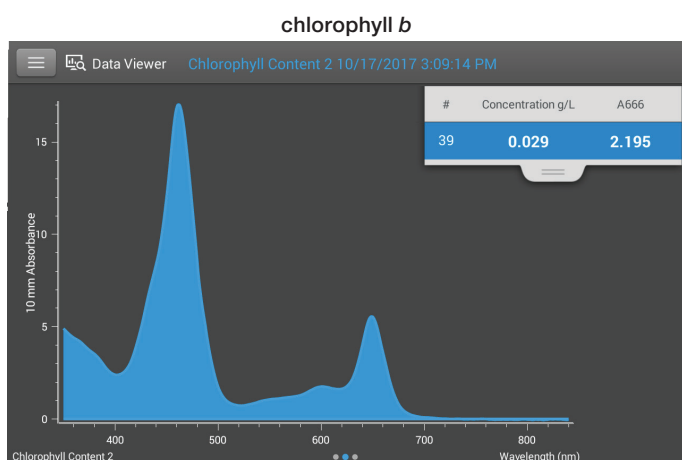
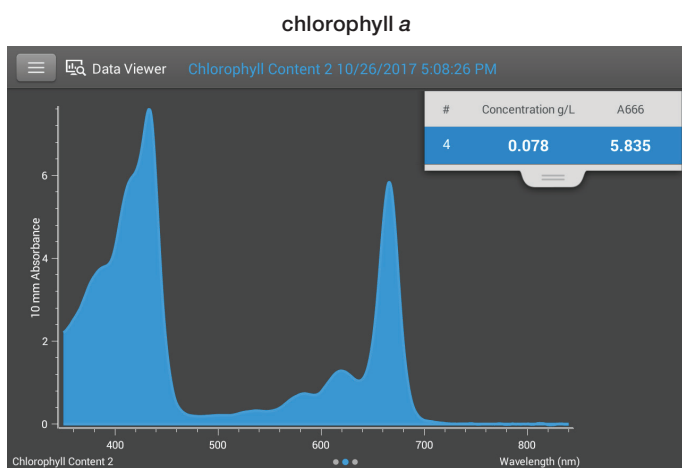
### Introduction

Chlorophyll *a* is the principal pigment that converts light energy to chemical energy, and chlorophyll *b* is the accessory photosynthetic pigment that transfers light it absorbs to chlorophyll *a*. Chlorophyll *a* is found in all plants, green algae, and cyanobacteria, and chlorophyll *b* is found in plants and green algae. Chlorophyll quantitation is valuable in a vast array of disciplines including but not limited to plant biology, environmental science, ecotoxicology, disease prevention, and medical drug discovery.

Spectrophotometry is a common method used to measure the absorbance of light by the chlorophyll molecules. The NanoDrop One/One<sup>c</sup> UV-Vis Spectrophotometer can be used to measure the absorbance of chlorophyll. Chlorophyll *a* and chlorophyll *b* absorb light at slightly different wavelengths. Chlorophyll *a* absorbs light at 433 nm and 666 nm and chlorophyll *b* absorbs light at 462 nm and 650 nm. The NanoDrop One/One<sup>c</sup> UV-Vis application can be used to observe the spectrum of each chlorophyll *a* and chlorophyll *b* and identify major absorbance



peaks (Figure 1). With this information, a user-defined custom method including user-defined formulas can be created to measure the absorbance and determine the concentration of chlorophyll.



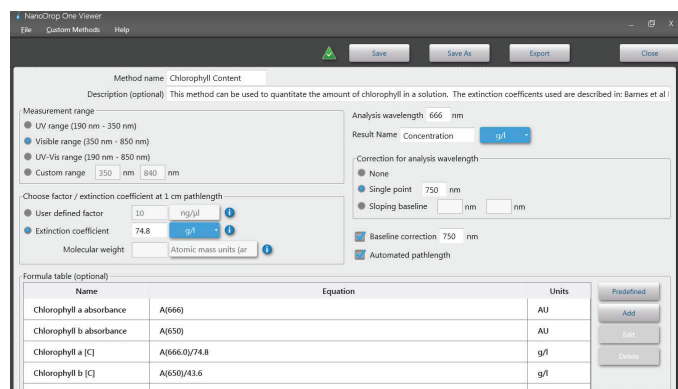
**Figure 1. Absorbance spectrum of chlorophyll a and chlorophyll b.** Chlorophyll a displayed major absorbance peaks at 433 nm and 666 nm (top). Chlorophyll b displayed major absorbance peaks at 462 nm and 650 nm (bottom).

## Methods

### Sample Preparation

Various solvents can be used to extract chlorophyll a and chlorophyll b. Porra et al. (1989) discuss assays of chlorophyll a and chlorophyll b suspended in various solvents including dimethylformamide, methanol, and 80% aqueous acetone. Barnes et al. (1991) describe the use of dimethyl sulfoxide (DMSO) for the extraction and determination of chlorophyll a and chlorophyll b. Based on the findings of Barnes et al., and the low toxicity and low vapor pressure of DMSO, we determined DMSO was an appropriate solvent for chlorophyll a and chlorophyll b measurements on the NanoDrop One/One<sup>c</sup> pedestal.

Pure chlorophyll a from spinach (Sigma-Aldrich® Product # C5753) and pure chlorophyll b from spinach (Sigma-Aldrich Product # C5878) were each dissolved in 100% DMSO. Half-fold serial dilutions for each suspension were prepared using DMSO. The samples were stored in amber tubes at -20°C until measured.



**Figure 2. Chlorophyll Content custom method created to quantify chlorophyll a and chlorophyll b samples suspended in 100% DMSO.**

### Chlorophyll Quantification Custom Method

The NanoDrop One/One<sup>c</sup> PC Viewer Custom Method application allows a user to specify how to calculate and report results. The Custom Method application was used to create a user-defined custom method to quantify chlorophyll a and chlorophyll b.

For this study, the following custom method parameters were used:

<b>Wavelength range</b>	Visible (350–850 nm)
<b>Extinction coefficient</b>	74.8 g/L
<b>Analysis wavelength</b>	666 nm
<b>Correction for analysis wavelength</b>	750 nm
<b>Baseline correction</b>	750 nm
<b>Automated pathlength</b>	On

Custom formulas were entered in the Formula table to report the absorbance at the analysis wavelength for chlorophyll a and chlorophyll b, and to calculate the concentration of pure chlorophyll a and pure chlorophyll b. Additional formulas were added to report the concentration of chlorophyll a, chlorophyll b, and chlorophyll a+b in samples containing a mixture of chlorophyll a and chlorophyll b. The formulas were taken from Barnes et al. (1991) and the peak locations were determined in the UV-Vis application (Figure 1).

The custom method was loaded in the Custom Method application on the NanoDrop One/One<sup>c</sup> local control. The custom method was run to measure each chlorophyll a and chlorophyll b serial dilution in triplicate.

**Note:** For this study, the NanoDrop One/One<sup>c</sup> UV-Vis application was used to identify the major peaks of chlorophyll a and chlorophyll b to determine the analysis

wavelength. If you are using a solvent other than DMSO, we recommend using the UV-Vis module to identify the maximum peaks for your samples, and modifying the custom method accordingly.

### Custom Method Download

1. Navigate to [www.thermofisher.com/nanodrop](http://www.thermofisher.com/nanodrop)
2. On the left, select "NanoDrop Software Download"
3. Choose the "NanoDrop One/One<sup>c</sup>" tab
4. Select "Local Control Software Download Instructions"
5. Scroll to "How to add a NanoDrop One/One<sup>c</sup> Custom Method file" and click on "Chlorophyll Content Method"
6. Unzip the custom method file and copy the .method file to a USB device and then follow the online "Instructions for uploading a Custom Method to the instrument from a USB device".

### Performance Data

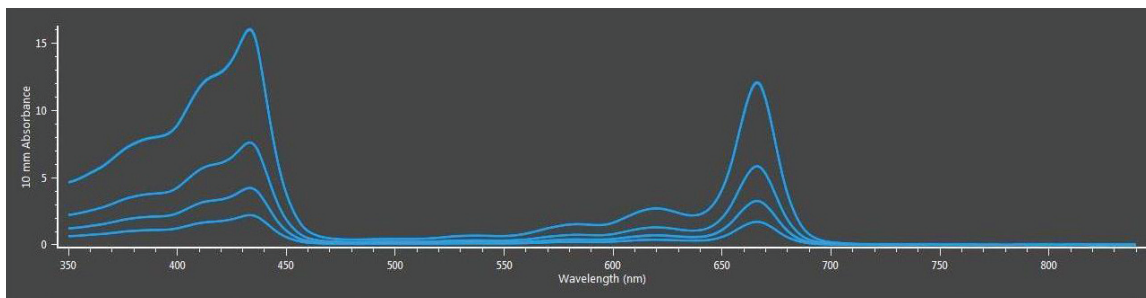
Each chlorophyll *a* and chlorophyll *b* dilution was measured using the user-defined custom method described above. Each sample was measured in triplicate to assess reproducibility (Tables 1 and 2). Standard deviation values in all cases were below 0.002 Abs (10 mm equivalent).

Chlorophyll <i>a</i>	A(666) (n=3)	Conc. (g/L) (n=3)	Std. dev.
Sample 1	12.096	0.162	0.001
Sample 2	5.841	0.078	0.000
Sample 3	3.270	0.044	0.000
Sample 4	1.705	0.023	0.000

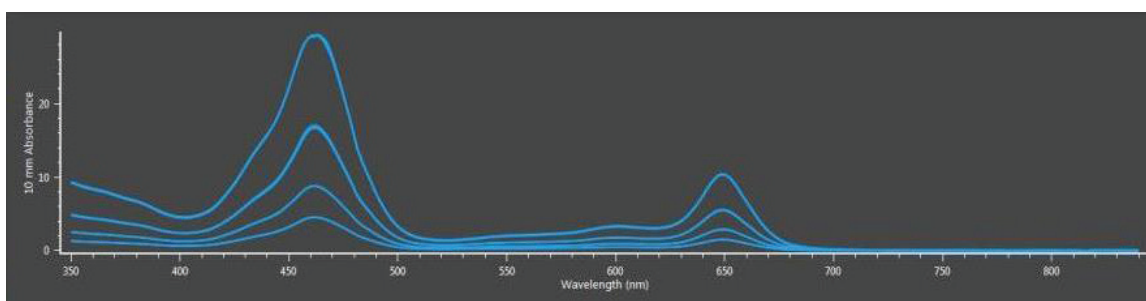
**Table 1. Chlorophyll *a* content.** The table displays calculated results for dilutions of chlorophyll *a* measured in triplicate.

Chlorophyll <i>b</i>	A(650) (n=3)	Conc. (g/L) (n=3)	Std. dev.
Sample 1	10.342	0.237	0.001
Sample 2	5.455	0.125	0.002
Sample 3	2.852	0.065	0.001
Sample 4	1.474	0.034	0.000

**Table 2. Chlorophyll *b* content.** The table displays calculated results for dilutions of chlorophyll *b* measured in triplicate.



**Figure 3. Custom method absorbance spectrum for each sample measurement of Chlorophyll *a* from spinach, Sigma-Aldrich Product # C5753, suspended in 100% DMSO.** Chlorophyll *a* displayed major absorbance peaks at 433 nm and 666 nm.



**Figure 4. Custom method absorbance spectrum for each sample measurement of Chlorophyll *b* from spinach, Sigma-Aldrich Product # C5878, suspended in 100% DMSO.** Chlorophyll *b* displayed major absorbance peaks at 462 nm and 650 nm.

## Conclusion

This study demonstrates the NanoDrop One/One<sup>c</sup> UV-Vis Microvolume Spectrophotometer can be used to accurately quantify chlorophyll *a* and chlorophyll *b* using a user-defined custom method. The low standard deviation indicates there was a high agreement between replicate measurements. This validates the NanoDrop One/One<sup>c</sup> Spectrophotometer can quantify samples accurately and reproducibly. The ability to measure chlorophyll *a* and *b* using a custom method serves as a valuable tool for research and the advancement of science.

## References

1. J.D. Barnes, L. Balaguer, E. Manrique, S. Elvira, and A.W. Davison (1991). A reappraisal of the use of DMSO for the extractions and determination of chlorophylls *a* and *b* in lichens and higher plants. *Environmental and Experimental Botany*. Vol. 32. No. 2, 85–100.
2. R.J. Porra, W.A. Thompson and P.E. Kriedemann. Determination of accurate extinction coefficients and simultaneous equations for assaying chlorophylls *a* and *b* extracted with four different solvents: verification of the concentration of chlorophyll standards by atomic absorption spectroscopy (1989). *Biochimica et Biophysica Acta*. 975, 384–394.

## Further Assistance and Technical Support

For further assistance, contact NanoDrop technical support at [nanodrop@thermofisher.com](mailto:nanodrop@thermofisher.com) or visit [thermofisher.com/nanodrop](http://thermofisher.com/nanodrop).

---

### In the United States:

For customer service, call 1-800-766-7000  
To fax an order, use 1-800-926-1166  
To order online: [thermofisher.com](http://thermofisher.com)

### In Canada:

For customer service, call 1-800-234-7437  
To fax an order, use 1-800-463-2996  
To order online: [thermofisher.ca](http://thermofisher.ca)

Find out more at [thermofisher.com/nanodrop](http://thermofisher.com/nanodrop)

**ThermoFisher**  
S C I E N T I F I C