

Analysis of Nitrate Nitrogen (NO_3^-) in Water by the EPA Approved Brucine Method

Gordon Bain, Ph.D., Michael W. Allen, Ph.D., Nicole Kreuziger Keppy, Thermo Fisher Scientific, Madison, WI, USA

Key Words

- Brucine Method
- Nitrate
- UV-Visible
- Water Analysis

Introduction

In this experiment quantitative analysis of nitrate nitrogen (NO_3^-) was performed by the Brucine method. Data was acquired and processed using the Thermo Scientific Evolution Array UV-Visible spectrophotometer and Thermo Scientific VISIONcollect software.

Background

The Brucine method for nitrate is approved by the United States EPA as Method No. 352.1.¹ When a water sample containing nitrate ion is treated with Brucine in sulfuric acid, a yellow solution results. The concentration of nitrate nitrogen may be calculated based upon the absorbance of the solution at 410 nm.

Experimental

Reagents and Apparatus

- Nitrate ion standard solution (0.001 mg NO_3^- N/mL)
- Sodium chloride solution (30% w/v)
 - Dissolve 30 g NaCl in 100 mL deionized water
- Sulfuric acid solution (4+1)
 - 4:1 = Sulfuric acid (H_2SO_4):deionized water
 - **CAUTION: Wear appropriate protective clothing and observe proper protocols for diluting sulfuric acid**
- Brucine-Sulfanilic acid solution
 - Dissolve 1 g brucine dihydrate ($\text{C}_{23}\text{H}_{26}\text{N}_2\text{O}_4 \cdot 2\text{H}_2\text{O}$) and 0.1 g sulfanilic acid ($\text{H}_2\text{NC}_6\text{H}_4\text{SO}_3\text{H} \cdot \text{H}_2\text{O}$) in 3 mL conc. HCl. Dilute to 100 mL with deionized water
- Evolution™ Array™ UV-Vis spectrophotometer with VISIONcollect™ software
- Cuvette (10 mm pathlength)

Standards and Measurement Procedure

1. Prepare a series of nitrate standards from the 0.001 mg/mL NO_3^- standard. Each of the standards should be an accurately measured volume between 1 mL and 10 mL of stock diluted to a total volume of 10.0 mL. Also prepare a 0 ppm reagent blank. Between four and six standards are generally considered to be sufficient to establish a robust calibration curve.
2. Treat each standard according to the procedure described in the sample work-up procedure below
3. In Quantification Standard mode
 - a. Measure the 0 ppm standard as the blank
 - b. Measure the absorbance of the remaining standards
4. In Quantification Sample mode, measure the absorbance of the unknown sample(s) and calculate the concentration

Sample Work-up Procedure

1. Measure a known, consistent volume (1–5 mL is a typical range) of sample into a calibrated tube or flask
2. Dilute with deionized water to a total volume of 10 mL
3. If samples are saline, add 2 mL sodium chloride solution
4. Add 10 mL sulfuric acid solution and mix well by shaking while cooling in flowing water.
Do not stopper the tube/flask.
CAUTION: This step is exothermic. The contents will heat up and pressure will result inside the container if it is stoppered.
5. Add 0.5 mL brucine in sulfanilic acid solution and mix by shaking
6. Heat 25 minutes at 100 °C in a water bath
7. Cool with flowing water, dilute to 25 mL with deionized water, stopper and invert several times to mix thoroughly
8. Pour ca. 3 mL of solution into the 10 mm pathlength cuvette and record the absorbance using the spectrophotometer

Instrument Parameters

Data collection was carried out with VISIONcollect software using the parameters displayed in Figure 1. The Curve Zero Offset and Curve Order settings specify a linear data fit that is not forced through the origin.

Experiment Type: Quantification Standard	
Experiment Setup	
Data Type	Absorbance
Sampling	Single Cell Holder
Mode	User Defined
Scan No.	30
Integration No.	1
Baseline Correction	
Quantification Standard	
Analysis Name	NO3_N
Concentration Unit	ppm
Use Wavelength (nm)	410
Curve Zero Offset	Yes
Curve Order	1
Standard Concentration...	

Figure 1: Experimental parameters for Nitrate assay

Results

Figure 2 shows the software screen for the standard solution data collection. Full spectra of all standards are displayed in the Standard Spectrum window (upper left). Details of the standards used and their measured absorbance values at the analytical wavelength (410 nm) are shown at the bottom of the screen and a calibration curve is plotted (upper right). The equation of the line of best fit is calculated by the least squares method and displayed below the Standard Spectrum.

A sample containing an unknown amount of nitrate was prepared according to the procedure and measured in duplicate. The results were displayed as shown in Figure 3. The concentration of the sample was found to be 0.23 ppm.

Conclusion

In this experiment quantitative analysis of nitrate nitrogen (NO_3^-) in water was successfully performed with the Evolution Array spectrophotometer and VISIONcollect software. Rapid acquisition of full-range spectra and excellent instrument sensitivity made the analysis quick, accurate and reproducible.

References

1. Electronic: http://www.epa.gov/waterscience/methods/method/files/352_1.pdf. Date acquired 12-12-2009.

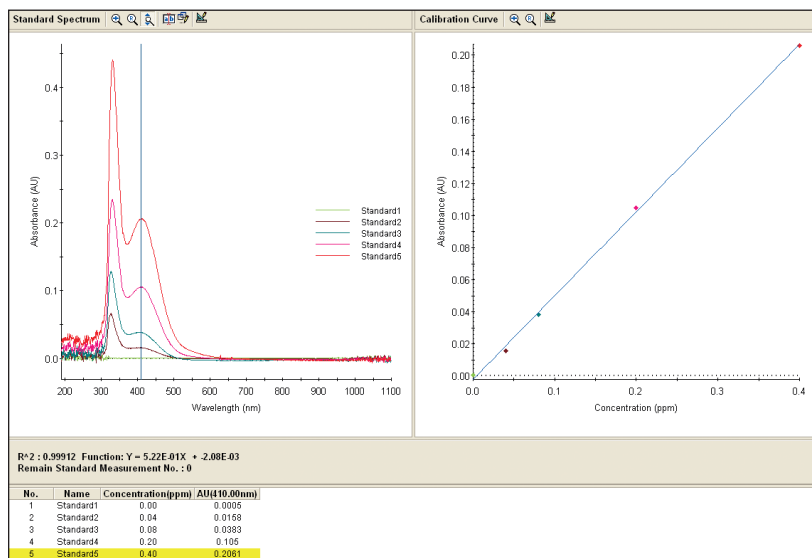


Figure 2: Standard solution data

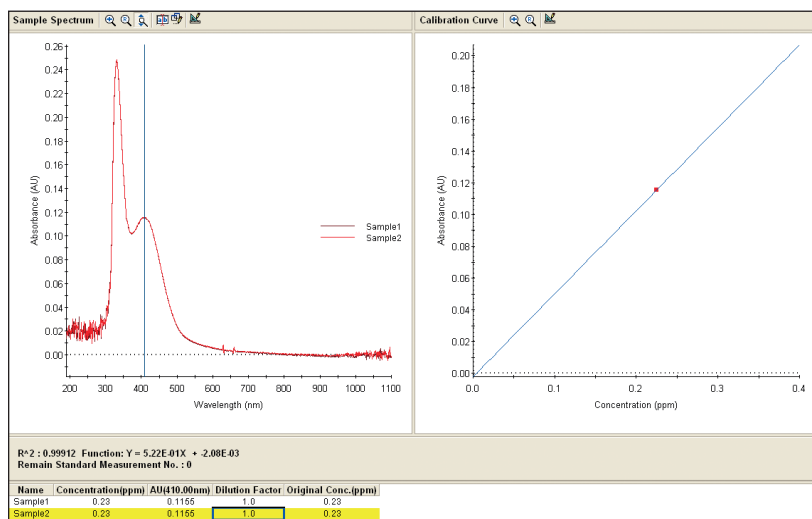


Figure 3: Sample solution data

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

Africa-Other
+27 11 570 1840

Australia
+61 2 8844 9500

Austria
+43 1 333 50 34 0

Belgium
+32 2 482 30 30

Canada
+1 800 530 8447

China
+86 10 8419 3588

Denmark
+45 70 23 62 60

Europe-Other
+43 1 333 50 34 0

Finland/Norway/Sweden
+46 8 556 468 00

France
+33 1 60 92 48 00

Germany
+49 6103 408 1014

India
+91 22 6742 9434

Italy
+39 02 950 591

Japan
+81 45 453 9100

Latin America
+1 608 276 5659

Middle East
+43 1 333 50 34 0

Netherlands
+31 76 579 55 55

South Africa
+27 11 570 1840

Spain
+34 914 845 965

Switzerland
+41 61 716 77 00

UK
+44 1442 233555

USA
+1 800 532 4752

www.thermo.com



Thermo Electron Scientific Instruments LLC, Madison, WI USA is ISO Certified.

AN51862_E 12/09M