

Ammonium in sewage sludge

Photometric determination using the indophenol blue method

Introduction

Ammonium is a key parameter in water testing, as it can have adverse effects on marine ecosystems at high levels. One major issue is eutrophication, excessive plant and algal growth.¹ As a result, determining ammonium levels during the process of sewage treatment is critical to maintaining aquatic environments. In this application note, we describe quantification of ammonium in sewage sludge with the Spectroquant[®] photometry system and ammonium test kits.

Experimental

Method

Ammonium nitrogen (NH₄-N) occurs partly in the form of ammonium ions and partly as ammonia. A pH-dependent equilibrium exists between the two forms. In strongly alkaline solution ammonium nitrogen is present almost entirely as ammonia, which reacts with hypochlorite ions to form monochloramine. This in turn reacts with a substituted phenol to form a blue indophenol derivative that is determined photometrically. Due to the intrinsic yellow coloration of the reagent blank, the measurement solution is yellow-green to green in color.

The method is analogous to EPA 350.1, APHA 4500-NH₃ F, ISO 7150-1, and DIN 38406-5.

Reagents and Instruments

Cat. No.	Product Description
Test Kits	
1.14558	Ammonium Cell Test Method: photometric 0.20 - 8.00 mg/l NH ₄ -N: 0.26 - 10.30 mg/l NH ₄ ⁺ Spectroquant [®] or
1.14544*	Ammonium Cell Test Method: photometric 0.5 - 16.0 mg/l NH ₄ -N: 0.6 - 20.6 mg/l NH ₄ ⁺ Spectroquant [®] or
1.14559	Ammonium Cell Test Method: photometric 4.0 - 80.0 mg/l NH ₄ -N: 5.2 - 103.0 mg/l NH ₄ ⁺ Spectroquant [®] or
1.14752	Ammonium Test Method: photometric 0.010 - 3.00 mg/l NH ₄ -N: 0.013 - 3.86 mg/l NH ₄ ⁺ Spectroquant [®] or
1.00683	Ammonium Test Method: photometric 2.0 - 150 mg/l NH ₄ -N: 2.6 - 193 mg/l NH ₄ ⁺ Spectroquant [®]
Instruments	
1.73026	Spectroquant [®] VIS Spectrophotometer Prove 100 plus or
1.73027	Spectroquant [®] UV/VIS Spectrophotometer Prove 300 plus or
1.73028	Spectroquant [®] UV/VIS Spectrophotometer Prove 600 plus or
1.09748	Spectroquant [®] Photometer NOVA 30 or
1.09751	Spectroquant [®] Photometer NOVA 60 or
1.09752	Spectroquant [®] Photometer NOVA 60A or
1.73632	Spectroquant [®] Colorimeter Move 100
Materials	
1.14946	Rectangular cells 10 mm or
1.14947	Rectangular cells 20 mm or
1.14944	Rectangular cells 50 mm or
WHA10314745	Folded filter, Whatman [®] prepleated qualitative filter paper, Grade 1573 1/2
1.16754	Water for analysis EMSURE
	Charcoal activated for soil tests

*not compatible with Move 100
Also first generation Prove instruments are compatible and preprogrammed with this method.

Analytical Approach

Sample preparation

In a 1000-ml volumetric flask mix 50 g of the sample with water for analysis and fill up to volume. Let the insoluble parts settle and decant 250 ml of the supernatant solution. Then mix with 2.5 g of charcoal activated for soil tests and stir for 10 minutes, before filtering through a folded filter.

Analysis

Determine with the above-mentioned test kits.

Calculation

Ammonium content in mg/l NH_4 = analysis value in mg/l NH_4 x 20.

References

1. Camargo JA, Alonso A. Ecological and toxicological effects of inorganic nitrogen pollution in aquatic ecosystems: A global assessment. Environ Int. Aug 2006;32(6):831-49. doi:10.1016/j.envint.2006.05.002.

To place an order or receive technical assistance

In Europe, please call Customer Service:

France: 0825 045 645

Germany: 069 86798021

Italy: 848 845 645

Spain: 901 516 645 Option 1

Switzerland: 0848 645 645

United Kingdom: 0870 900 4645

For other countries across Europe, please call: +44 (0) 115 943 0840

Or visit: [MerckMillipore.com/offices](https://www.MerckMillipore.com/offices)

For Technical Service visit: [MerckMillipore.com/techservice](https://www.MerckMillipore.com/techservice)

[MerckMillipore.com](https://www.MerckMillipore.com)

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

