

QuantiFluo™ Ammonia Assay Kit II (D2NH3-100)

Quantitative Fluorimetric Determination of Ammonia/Ammonium Concentration

DESCRIPTION

AMMONIA (NH₃) or its ion form ammonium (NH₄⁺) is an important source of nitrogen for living systems and is ubiquitously present in the nature. Simple, direct and automation-ready procedures for measuring NH₃ are very desirable. BioAssay Systems' updated ammonia/ammonium assay is based on an improved *o*-phthalaldehyde method. This reagent reacts with ammonia/ ammonium and forms a fluorescent product. The fluorescence intensity ($\lambda_{\text{ex/em}} = 415/475\text{nm}$) is proportional to the ammonia concentration in the sample.

KEY FEATURES

Fast and sensitive. Linear detection range of 0.005 - 2mM ammonia.

Convenient and high-throughput. A single component reagent. Homogeneous "mix-incubate-measure" type assay. Can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

APPLICATIONS

Ammonia/ammonium determination in biological (e.g. urine, serum, plasma) and environmental samples.

KIT CONTENTS

Detection Reagent: 10 mL **Standard:** 100 μL NH₄Cl

Storage conditions: This product is shipped at room temperature. Store kit at -20°C. Shelf life of 6 months after receipt.

Precautions: Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

ASSAY PROCEDURE

Use black flat-bottom 96-well plates. Prior to assay, bring all reagents to room temperature.

Note: (1). This assay is compatible with most detergents, chelators and buffer components. Protein concentration up to 17 $\mu\text{g/mL}$ was tolerated and primary amine-containing buffers (e.g. Tris, glycine) should be kept below 10 mM, if possible. For best results, include the same concentration of the sample buffer in the standards and blank. (2). Samples should be clear and not contain any particles or precipitates. Particles or precipitates can be removed by centrifugation for 5 min at 14,000 rpm or by filtration. (3). Urine samples should be diluted 50-fold in water prior to assay. Serum and plasma samples should be diluted 10-fold in water prior to assay.

1. **Standards.** Prepare 200 μL 1 mM Standard Premix by mixing 10 μL 20 mM NH₄Cl Standard and 190 μL H₂O. Dilute standards as follows.

No	Premix + H ₂ O	Standard mM)
1	100 μL + 0 μL	1.00
2	50 μL + 50 μL	0.50
3	25 μL + 75 μL	0.25
4	0 μL + 100 μL	0

Transfer 10 μL standards into separate wells of the plate.

Transfer 10 μL of each sample in separate wells of the plate.

2. **Assay.** Add 90 μL Detection Reagent to each well. Immediately tap plate to mix. Incubate for 20 min in the dark at room temperature. Measure fluorescence intensity at $\lambda_{\text{ex/em}} 415/475\text{nm}$ on a plate reader.

CALCULATION

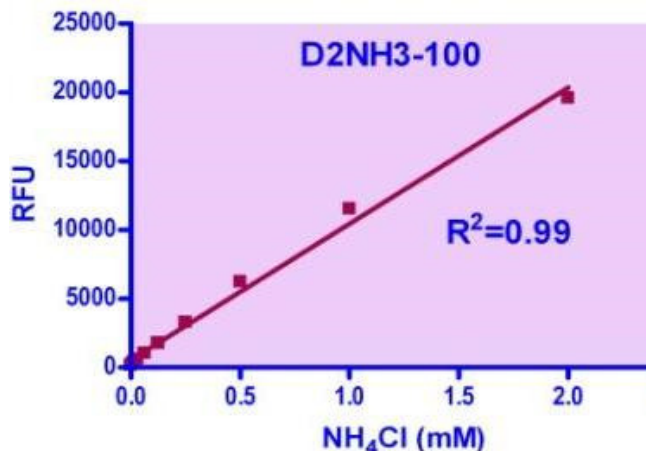
Plot the ammonia standard curve and determine its Slope. The ammonia concentration of a Sample is calculated as

$$[\text{NH}_3] = \frac{F_{\text{SAMPLE}} - F_{\text{BLANK}}}{\text{Slope}} \text{ (mM)}$$

where F_{SAMPLE} and F_{BLANK} are the fluorescence intensity values of the Sample and the blank (i.e. #4 H₂O), respectively. If ammonia concentration is higher than 2 mM, dilute Sample in water and repeat assay. Multiply the results by the dilution factor.

MATERIAL REQUIRED BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, black flat bottom 96-well plates and plate reader.



Standard curve performed on a 96-well plate reader (Spectramax M2);

LITERATURE

1. Gips CH, et al (1970). Preservation of urine for ammonia determination with a direct method. Clin Chim Acta. 29(3):501-5.
2. Beecher GR, Whitten BK (1970). Ammonia determination: reagent modification and interfering compounds. Anal Biochem. 36(1):243-6.
3. Rodger MR, Jenkins P (1984). Enzymic fluorometric assay of plasma ammonia with a centrifugal analyzer. Clin Chem. 30(10):1670-2.
4. Sugawara K, Oyama F (1981). Fluorogenic reaction and specific microdetermination of ammonia. J. Biochem. 89:771-774.

