

## EnzyChrom™ Ammonia/Ammonium Assay Kit (ENH3-100)

### Quantitative Colorimetric/Fluorimetric Determination of Ammonia

#### DESCRIPTION

AMMONIA (NH<sub>3</sub>) or its ion form ammonium (NH<sub>4</sub><sup>+</sup>) is an important source of nitrogen for living systems. It is synthesized through amino acid metabolism and is toxic when present at high concentrations. In the liver, ammonia is converted to urea through the urea cycle. Elevated levels of ammonia in the blood (hyperammonemia) have been found in liver dysfunction (cirrhosis), while hypoammonemia has been associated with defects in the urea cycle enzymes (e.g. ornithine transcarbamylase).

Simple, direct and automation-ready procedures for measuring NH<sub>3</sub> are popular in research and drug discovery. BioAssay Systems' ammonia assay is designed to directly measure NH<sub>3</sub> and NH<sub>4</sub><sup>+</sup>. In this assay, NADH is converted to NAD<sup>+</sup> in the presence of NH<sub>3</sub>, ketoglutarate and glutamate dehydrogenase. The decrease in optical density at 340 nm or fluorescence intensity at λ<sub>em/ex</sub> = 450/360 nm is directly proportionate to the NH<sub>3</sub> concentration in the sample.

#### KEY FEATURES

**High sensitivity and wide linear range.** Use 20 μL sample. Linear detection range 24 to 1000 μM ammonia.

**Homogeneous and simple procedure.** Simple "mix-and-measure" procedure allows reliable quantitation of NH<sub>3</sub> within 30 minutes.

#### APPLICATIONS

**Direct Assays:** NH<sub>3</sub> in biological samples (e.g. serum, plasma, urine, saliva, cell culture etc).

#### KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

**Assay Buffer:** 20 mL                      **Enzyme:** 120 μL  
**Ketoglutarate:** 120 μL                **Standard:** 400 μL  
**NADH Reagent:** Dried

**Storage conditions.** The kit is shipped on ice. Store all components at -20°C. Shelf life of six months after receipt, 3 weeks after reconstitution.

**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.

#### PROCEDURES

**Reagent Preparation.** Equilibrate all components to room temperature. Briefly centrifuge all tubes before opening. Reconstitute the NADH Reagent tube with 1000 μL dH<sub>2</sub>O (final 10 mM). Unused reconstituted NADH reagent is stable for three weeks when stored frozen at -20°C.

**Sample preparation:** solid samples can be extracted by homogenization in distilled water (dH<sub>2</sub>O) and filtered, centrifuged or, if necessary, deproteinized to remove any undissolved material. Samples should be clear and colorless with pH adjusted to 7 - 8.

Serum and plasma samples can be assayed directly. Cell culture media should be diluted 5-10 fold in dH<sub>2</sub>O prior to assay.

#### Colorimetric Procedure

1. **Standards and Samples.** Prepare a 1000 μM NH<sub>3</sub> Standard Premix by mixing 15 μL of the 20 mM Standard and 285 μL dH<sub>2</sub>O. Dilute Standard as follows.

No	Premix + dH <sub>2</sub> O	Vol (μL)	NH <sub>3</sub> (μM)
1	100 μL + 0 μL	100	1000
2	60 μL + 40 μL	100	600
3	30 μL + 70 μL	100	300
4	0 μL + 100 μL	100	0

Transfer 20 μL standards into separate wells of a clear, flat-bottom 96-well plate.

Transfer 20 μL of each sample into two separate wells, one serving as a sample blank well (R<sub>BLANK</sub>) and one as a sample well (R<sub>SAMPLE</sub>).

2. **Enzyme Reaction.** For each standard and sample well, prepare Working Reagent by mixing 180 μL Assay Buffer, 1 μL Enzyme, 8 μL reconstituted NADH Reagent and 1 μL Ketoglutarate. Add 180 μL Working Reagent to the *four Standards* and the *Sample Wells*.

Prepare blank control reagent by mixing 180 μL Assay Buffer, 8 μL reconstituted NADH Reagent and 1 μL Ketoglutarate (*No Enzyme*). Add 180 μL Blank control reagent only to the *Sample Blank Wells*.

Tap plate to mix. Incubate 30 min at room temperature.

3. Read OD<sub>340nm</sub>.

#### Fluorimetric Procedure

The fluorimetric procedure is the same as for the colorimetric assay, except that a black, flat-bottom 96-well plate is used. After incubation for 30 min at room temperature, read fluorescence intensity at λ<sub>ex</sub> = 350-360 nm and λ<sub>em</sub> = 450 nm.

#### CALCULATION

Subtract the standard values from the blank value (#4) and plot the ΔOD or ΔF against standard concentrations. Determine the slope and calculate the NH<sub>3</sub> concentration of Sample,

$$[\text{Ammonia}] = \frac{R_{\text{BLANK}} - R_{\text{SAMPLE}}}{\text{Slope } (\mu\text{M}^{-1})} \times n \quad (\mu\text{M})$$

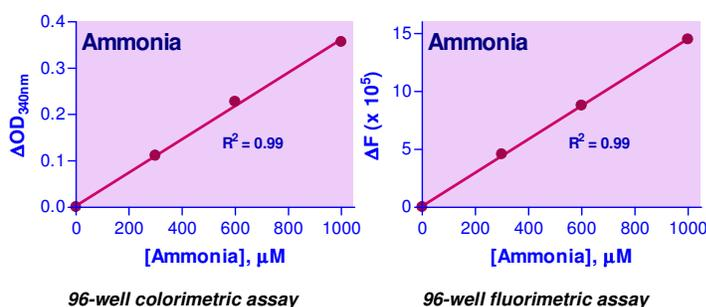
R<sub>SAMPLE</sub> and R<sub>BLANK</sub> are optical density or fluorescence intensity readings of the Sample and Sample Blank, respectively. *n* is the sample dilution factor.

*Note:* if the calculated NH<sub>3</sub> concentration is higher than 1000 μM, dilute sample in dH<sub>2</sub>O and repeat assay. Multiply result by the dilution factor *n*.

*Conversions:* 1000 μM NH<sub>3</sub> equals 1.7 mg/dL or 17 ppm.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipeting devices, and clear flat-bottom 96-well plates and optical density plate reader for colorimetric assays; black flat-bottom 96-well plate and fluorescence intensity plate reader for fluorimetric assays.



#### LITERATURE

- Bruce, A.W. et al. (1978). Two-point determination of plasma ammonia with the centrifugal analyzer. Clin Chem. 24:782-787.
- Mondzac, A. et al. (1965). An enzymatic determination of ammonia in biological fluids. J Lab Clin Med. 66:526-531.
- Seligson, D. and Hirahara, K. (1957). The measurement of ammonia in whole blood, erythrocytes, and plasma. J Lab Clin Med. 49:962-974.