

Product Information

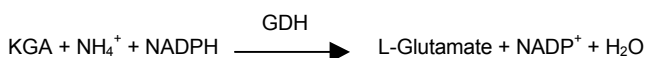
Ammonia Assay Kit

Catalog Number **AA0100**
Storage Temperature 2–8 °C
DO NOT FREEZE

TECHNICAL BULLETIN

Product Description

This kit is for the quantitative, enzymatic determination of ammonia in food and biological samples. Ammonia reacts with α -ketoglutaric acid (KGA) and reduced nicotinamide adenine dinucleotide phosphate (NADPH) in the presence of L-glutamate dehydrogenase (GDH) to form L-glutamate and oxidized nicotinamide adenine dinucleotide phosphate (NADP⁺) as follows:



The decrease in absorbance at 340 nm, due to the oxidation of NADPH, is proportional to the ammonia concentration. L-Glutamate dehydrogenase reacts specifically with ammonia. The Ammonia Assay Kit may be used to determine ammonia concentrations in the range of 0.2–15 $\mu\text{g/ml}$.

Components

Sufficient reagents are provided for 100 assays.

Ammonia Assay Reagent 10 vials
(Catalog Number A0853)

The dry reagent contains α -ketoglutaric acid, NADPH, buffers, stabilizers, and nonreactive fillers. The reagent should be stored at 2–8 °C and it should be discarded if the vials exhibit caking due to possible moisture penetration, if the vial contents do not dissolve completely upon reconstitution, or if the reconstituted solution appears turbid or yellow in color.

L-Glutamate Dehydrogenase, 1 vial
bovine liver
(Catalog Number G2294)

The enzyme is supplied as a solution in phosphate buffer, pH 7.4, with 50% glycerol. The enzyme solution should be stored at 2–8 °C. DO NOT FREEZE.

This reagent is stable for at least 1 year at 2–8 °C.

Activity: $\geq 1,200$ units/ml

Unit Definition: One unit will reduce 1.0 μmole of α -ketoglutaric acid to L-glutamate per minute at pH 7.3 and 25 °C in the presence of ammonium ions.

Ammonia Standard Solution 1 vial
(Catalog Number A0978)

The ammonia concentration of the standard solution is ~ 10 $\mu\text{g/ml}$ (~ 588 μM as ammonium sulfate). The solution should be stored at 2–8 °C and is stable for at least 1 year. To avoid evaporation or ammonia uptake, **the Ammonia Standard Solution should be kept capped when not being used.** The Ammonia Standard Solution is used to ensure assay reliability. It is not necessary to run the standard to determine the ammonia concentration of samples.

Equipment Required but Not Provided

- Spectrophotometer suitable for measuring absorbance at 340 nm
- Cuvettes
- Pipettes capable of accurately dispensing 10 μl to 1 ml

Precautions and Disclaimer

The Ammonia Assay Kit is R&D use only, not for *in vitro* diagnostic use, drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Ammonia Assay Reagent - Reconstitute each vial with 10 ml of water. After addition of the water, cap the vial and immediately mix several times by inversion. DO NOT SHAKE. The prepared Ammonia Assay Reagent should be kept capped to avoid possible ammonia uptake. The prepared Ammonia Assay Reagent contains ~ 3.4 mM α -ketoglutaric acid and ~ 0.23 mM NADPH. It is stable, in the absence of visible microbial growth, for at least 1 day at 18–26 °C and for 2 weeks at 2–8 °C. The reconstituted reagent may be aliquoted and stored at –20 °C for 6 months. Avoid freeze/thaw cycles.

The L-Glutamate Dehydrogenase solution and the Ammonia Standard Solution are supplied ready-to-use.

Sample preparation - Samples should be clear and colorless. Adjust the pH to 7–8.

Liquid samples should be diluted with water to an ammonia concentration of 0.20–150 $\mu\text{g/ml}$ (0.1 ml final sample volume). Filter or deproteinize sample solutions, if necessary, to clarify. Sample solutions that are strongly colored and have a low ammonia concentration should be decolorized. Carbonated or fermented products must be degassed.

Solid samples should be weighed to nearest 0.1 mg. Extract the solid sample with water. Dilute the extract with water to an ammonia concentration of 0.20–150 $\mu\text{g/ml}$ (0.1 ml final sample volume). Filter or deproteinize solution, if necessary, to clarify.

Procedure

Any compounds present in the sample that react with $\text{NADP}^+/\text{NADPH}$ under the conditions of this assay will result in inaccurate results for determination of the ammonia concentration. Samples containing protein or protein fragments that liberate ammonia from amino acids will result in falsely high values of ammonia. L-Glutamate dehydrogenase is inhibited by heavy metals such as copper, silver, mercury, iron, and zinc. High concentrations of tannins in the sample will slow the reaction due to inhibition of the L-glutamate dehydrogenase.

Make sure the work area is free from ammonia sources.

1. Pipette the following solutions into appropriately marked cuvettes.

Cuvette	Ammonia Assay Reagent	Ammonia Standard Solution	Sample Volume	Water
Reagent Blank	1.0 ml			100 μl
Test	1.0 ml		100 μl	
Standard	1.0 ml	0.05 ml		

Sample volume may be varied from 10–200 μl as needed. If the sample volume is changed, make the appropriate adjustments in the calculations.

2. Set the spectrophotometer to 340 nm and the absorbance to zero using water as the reference.
3. Mix the contents in each cuvette and incubate for ~5 minutes at 18–35 °C. Measure the absorbance of each solution at 340 nm.
4. Add 10 μl of L-Glutamate Dehydrogenase solution (Catalog Number G2294) to each cuvette.
5. Mix the contents of each cuvette and incubate for ~5 minutes at 18–35 °C. Then measure the absorbance of each solution at 340 nm.
Note: The L-glutamate dehydrogenase reaction is typically complete in five minutes. If the absorbance at 340 nm continues to decrease after five minutes, a competitive reaction involving the oxidation of NADPH may be occurring. Continue measuring the absorbance at 1 minute intervals until the rate of decrease is constant for 2 minutes. From this constant rate of decrease, calculate the change in absorbance for 5 minutes due to the competitive reaction [$\Delta A_{340}(\text{CR})$]. When calculating the ammonia concentration, $\Delta A_{340}(\text{CR})$ must be subtracted from $\Delta A_{340}(\text{Test})$.

Calculation:

Determine ΔA_{340} for the Reagent Blank, Test, and Standard. For each: $\Delta A_{340} = A_{\text{Initial}} - A_{\text{Final}}$

$\Delta(\Delta A_{340})_{\text{Test or Standard}}$

$$= \Delta A_{340}(\text{Test or Standard}) - \Delta A_{340}(\text{Blank})$$

mg of NH_3 /ml of original sample

$$= \frac{(A) (TV) (MW \text{ of Ammonia}) (F)}{(\epsilon) (d) (SV) (\text{Conversion Factor for } \mu\text{g to mg})}$$

$$= \frac{(A) (TV) (17) (F)}{(6.22) (1) (SV) (1,000)}$$

$$= \frac{(A) (TV) (F)}{(SV)} \times 0.00273$$

$A = \Delta(\Delta A_{340})_{\text{Test or Standard}}$

TV = Total Assay Volume in ml

SV = Sample Volume in ml

MW of Ammonia = 17 g/mole or
equivalently 17 $\mu\text{g}/\mu\text{mole}$

F = Dilution Factor from Sample Preparation

ϵ = Millimolar Extinction Coefficient for NADPH at
340 nm [$\text{mM}^{-1} \text{cm}^{-1}$ or equivalently
(ml/ μmoles)(1/cm)]

d = Light path (cm) = 1 cm

Notes:

This kit may be used to determine ammonia concentrations as low as 0.2 $\mu\text{g}/\text{ml}$

$[\Delta(\Delta A_{340})_{\text{Test}} = 0.007]$ and as high as 15 $\mu\text{g}/\text{ml}$

$[\Delta(\Delta A_{340})_{\text{Test}} = 0.50]$.

The relative standard deviation is approximately 1–2%.

The Ammonia Standard Solution should be included with each set of assays. If the value obtained with the Ammonia Standard Solution is within 5% of the stated concentration, the test performance is acceptable.

References

1. Bergmeyer, H.U., and Beutler, H-O., in *Methods of Enzymatic Analysis*, 3rd Edition, vol. VIII, Bergmeyer, H. U., ed., Academic Press (New York, NY: 1985), pp 454-461.
2. Cheuk, W.L., and Finne, G., Enzymatic Determination of Urea and Ammonia in Refrigerated Seafood Products. *J. Agric. Food Chem.*, **32**, 14-18 (1984).
3. Mondzac, A., et al., An Enzymatic Determination of Ammonia in Biological Fluids. *J. Lab. & Clin. Med.*, **66**, 526-531 (1965).
4. Olson, J.A., and Anfinson, C.B., Kinetic and Equilibrium Studies on Crystalline L-glutamic Acid Dehydrogenase. *J. Biol. Chem.*, **202**, 841-856 (1953).
5. Van Anken, H.C., and Schiphorst, M.E., A Kinetic Determination of Ammonia in Plasma, *Clinica Chimica Acta*, **56**, 151-157 (1974).
6. Neeley, W.E., and Phillipson, J., Automated Enzymatic Method for Determining Ammonia in Plasma, with 14-day Reagent Stability. *Clin. Chem.*, **34**, 1868-1869 (1988).
7. Ratliff, C.R., and Hall, F.F., Blood Ammonia returns to the Laboratory. *Laboratory Management*, August, 16-23 (1979).

RG,CMH,MAM 11/08-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.