
Product Manual

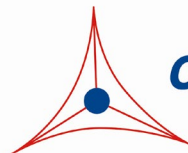
Urea Assay Kit (Fluorometric)

Catalog Number

MET-5180

200 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Urea, or carbamide, is the end product of protein nitrogen metabolism and is the primary vehicle for removing toxic ammonia from the body. Urea is synthesized in the liver from the ammonia produced from the catabolism of amino acids via the hepatic urea cycle. The conversion from ammonia to urea is regulated by N-acetylglutamate, which activates carbamoyl phosphate synthetase in the urea cycle. Urea is transported in the blood to the kidneys where it is excreted in the urine. In addition to its role as a carrier of waste nitrogen, urea also has a role in the countercurrent exchange system of the nephrons in which water and ions are re-absorbed from excreted urine. It is freely filtered by the glomeruli and partially passively resorbed as filtrate transverses the renal tubules. Urea reabsorption is inversely proportional to urine flow rate. Consequently, urea concentration depends upon protein intake, protein catabolism, and kidney function.

Urea quantitation is one of the most widely applied tests for kidney function evaluation. The analysis of urea in serum, plasma and urine is an important clinical test for renal disease and dysfunction. The test is frequently tested in conjunction with creatinine determination for diagnosis of pre-renal, renal, and post renal uremia. Toxic urea levels are associated with renal, liver, or other system dysfunction. Pre-renal uremia relates to water depletion, increased protein catabolism, infection, hypovolemia, or cardiac decomposition. Glomerulonephritis, tubular necrosis, nephrosclerosis, chronic nephritis, and polycystic kidney are examples of renal uremia, while post renal uremia is predominantly urinary tract obstructions or leakage. Increased urea levels can also be linked to other disease states such as liver disease, diabetes, and congestive heart failure. High plasma urea levels are known as Azotemia. Decreased urea levels are associated with acute hepatic insufficiency or excess parenteral fluid therapy.

Cell Biolabs' Urea Assay Kit (Fluorometric) is based on the enzymatic conversion of urea to ammonia and carbon dioxide, which further reacts with a compound to produce a fluorescent product that can be measured with a standard fluorometric plate. Each kit provides sufficient reagents to perform up to 200 assays, including blanks, urea standards and unknown samples and controls.

****Note: Each sample replicate requires 2 assays, one treated with Urease/Ammonia Reagent mixture and one with the Ammonia Reagent without Urease (-U). Urea is calculated from the difference in RFU readings from the 2 wells.***

Assay Principle

Cell Biolabs' Urea Assay Kit measures urea levels within urine, serum, plasma, cell lysates, or tissue homogenates. Samples are compared to a known concentration of urea standard within a 96-well microtiter plate format. Samples and standards are incubated for 10 minutes with the enzyme urease, which hydrolyzes urea to ammonia and CO₂. The ammonia reacts further with a compound to produce a fluorescent product. After 15-30 minutes, the plate is read with a standard black 96-well fluorometric microplate reader at an excitation wavelength of 360 nm and emission wavelength of 415 nm. Sample urea concentrations are determined by comparison with the known urea standards. The kit has a sensitivity of 0.1 mg/dL urea.

Related Products

1. STA-382: Urea Assay Kit (Colorimetric)

2. STA-375: Uric Acid/Uricase Assay Kit
3. STA-378: Creatinine Assay Kit

Kit Components

Box 1 (shipped at room temperature)

1. Urea Standard (Part No. 238201): One 250 μ L tube of a 1000 mg/dL solution.
2. Ammonia Reagent (Part No. 238202): One 20 mL amber bottle.
3. Fluorometric Reagent (Part No. 51801B): Two 20 mL amber bottles.
4. 10X Assay Buffer (Part No. 238205): One 10 mL bottle.

Box 2 (shipped on blue ice packs)

1. Urease (Part No. 238204): One 200 mg amber tube of powder.

Materials Not Supplied

1. Deionized water
2. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
3. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
4. Multichannel micropipette reservoirs
5. 37°C incubator
6. Fluorometric microplate reader capable of reading excitation 360 nm and emission 415 nm

Storage

Upon receipt, prepare aliquots and store the Urea Standard and Urease at -20°C. Store the remaining kit components at 4°C.

Preparation of Reagents

- 1X Assay Buffer: Dilute the Assay Buffer 1:10 with deionized water. Mix to homogeneity. Store the 1X Assay Buffer at 4°C up to six months.
- Urease/Ammonia Reagent*: Immediately prior to use, reconstitute the Urease enzyme at 4 mg/mL in the Ammonia Reagent solution and mix thoroughly until dissolved (e.g. for a 10 mL solution or 100 assays, add 40 mg of Urease to 10 mL Ammonia Reagent). Prepare only enough for immediate use. Do not store the Urease/Ammonia Reagent solution.

**Note: High levels of ammonia may cause high background values, so we recommend two wells per sample for testing: one well treated with the Urease (+U) added to the Ammonia Reagent, and one well with Urease omitted from the Ammonia Reagent (-U). The urea concentration is calculated from the difference in OD readings between the two assay values.*

Preparation of Samples

These preparation protocols are intended as a guide for preparing unknown samples. The user may need to adjust the sample treatment accordingly. All samples should be assayed immediately or stored for up to 2 months at -80°C. A trial assay with a representative test sample should be performed to determine the sample compatibility with the dynamic range of the standard curve. A set of serial dilutions is recommended for samples to achieve optimal assay results and minimize possible interfering chromogens. High levels of interfering substances may cause variations in results. Samples may be diluted in 1X Assay Buffer or deionized water as necessary before testing. Run proper controls as necessary. Always run a standard curve with samples.

- Serum: Avoid hemolyzed and lipemic blood samples. Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C. Perform dilutions in deionized water.
- Plasma: Avoid hemolyzed and lipemic blood samples. Collect blood sample and add to a blood collection tube containing Heparin as the anticoagulant. Centrifuge at 3,000 rpm for 10-15 minutes at 4°C. Remove the upper yellow plasma supernatant layer without disturbing the white buffy coat (leukocytes). Samples should be tested immediately or frozen at -80°C. Perform dilutions in deionized water.
- Urine: Urine samples with visible particulates should be centrifuged or filtered prior to testing. A minimum 1:20 dilution of urine samples into deionized water is recommended to remove matrix interference and achieve optimal assay results. Diluted samples should be used within 2 hours upon preparation.
- Tissue or Lysates: Homogenize 20 mg of tissue or 2×10^6 cells in cold 1X Assay Buffer. Centrifuge at 14000 x g for 10 min to remove insoluble material. Collect the supernatant. Samples can be tested directly or diluted with 1X Assay Buffer. Test samples immediately or store at -80°C.

Notes:

- *Buffers containing MES, HEPES, CHES, EDTA, fluoride, 2-mercaptoethanol, acetohydroxamate, 1,4-benzoquinone, or phosphoramidate are not recommended because they can inhibit urease activity.*
- *Do not use ammonium or potassium salts or fluoride as anticoagulants. Citrate, sodium heparin or oxalate can be used. All samples must be free of heavy metals.*
- *Hemoglobin (>200 mg/dL), Bilirubin (>20 mg/dL), and Triglycerides (>800 mg/dL) may interfere with the assay. Use controls accordingly.*
- *Drug interferences are possible (see Young, D.S., et. al).*

Preparation of Urea Standard Curve

Prepare a dilution series of Urea standards in the concentration range of 0 mg/dL to 6.25 mg/dL by diluting the provided Urea Standard in deionized water (Table 1).

Standard Tubes	1000 mg/dL Urea Standard (μL)	Deionized Water (μL)	Resulting Urea Concentration (mg/dL)
1	4	636	6.25
2	250 of Tube #1	250	3.13
3	250 of Tube #2	250	1.56
4	250 of Tube #3	250	0.78
5	250 of Tube #4	250	0.39
6	250 of Tube #5	250	0.195
7	250 of Tube #6	250	0.098
8	0	500	0

Table 1. Preparation of Urea Standards.

Note: Do not store diluted urea standard solutions.

Assay Protocol

Each urea standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

Note: Samples require two paired wells, one to be treated with the Urease/Ammonia Reagent mixture (+U), and one with the Ammonia Reagent without Urease (-U). This controlled pairing will allow for the determination of background ammonia levels within the sample.

1. Add 10 μL of the diluted urea standards or samples to the 96-well black plate wells. Remember to provide two sample paired wells if testing for ammonia background.
2. Add 100 μL of the Urease/Ammonia Reagent (+U) mixture (see Preparation of Reagents section) to each standard and to one half of the paired sample wells using either a multichannel pipette. Mix thoroughly and carefully to avoid foaming in the well.
3. Add 100 μL of the Ammonia Reagent (-U) to the other half of the paired sample wells using either a multichannel pipette. Mix thoroughly and carefully to avoid foaming in the well.
4. Incubate 10 minutes at 37°C.
5. Add 200 μL of the Fluorometric Reagent to each well using either a multichannel pipette. Mix the solution thoroughly and carefully to avoid foaming in the well.
6. Incubate 15-30 minutes at 37°C.
7. Read the plate at 360 nm excitation wavelength 415 nm emission wavelength and record data.

Example of Results

The following figures demonstrate typical Urea Assay results. One should use the data below for reference only. This data should not be used to interpret actual sample results.

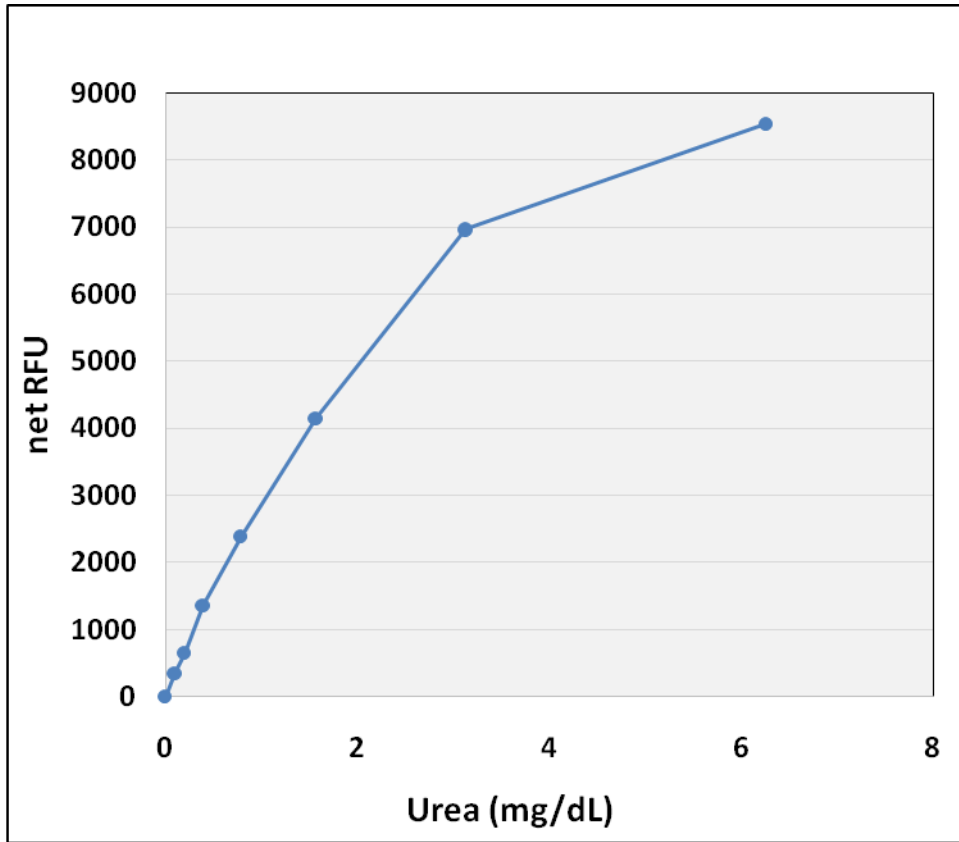


Figure 2: Urea Assay Standard Curve.

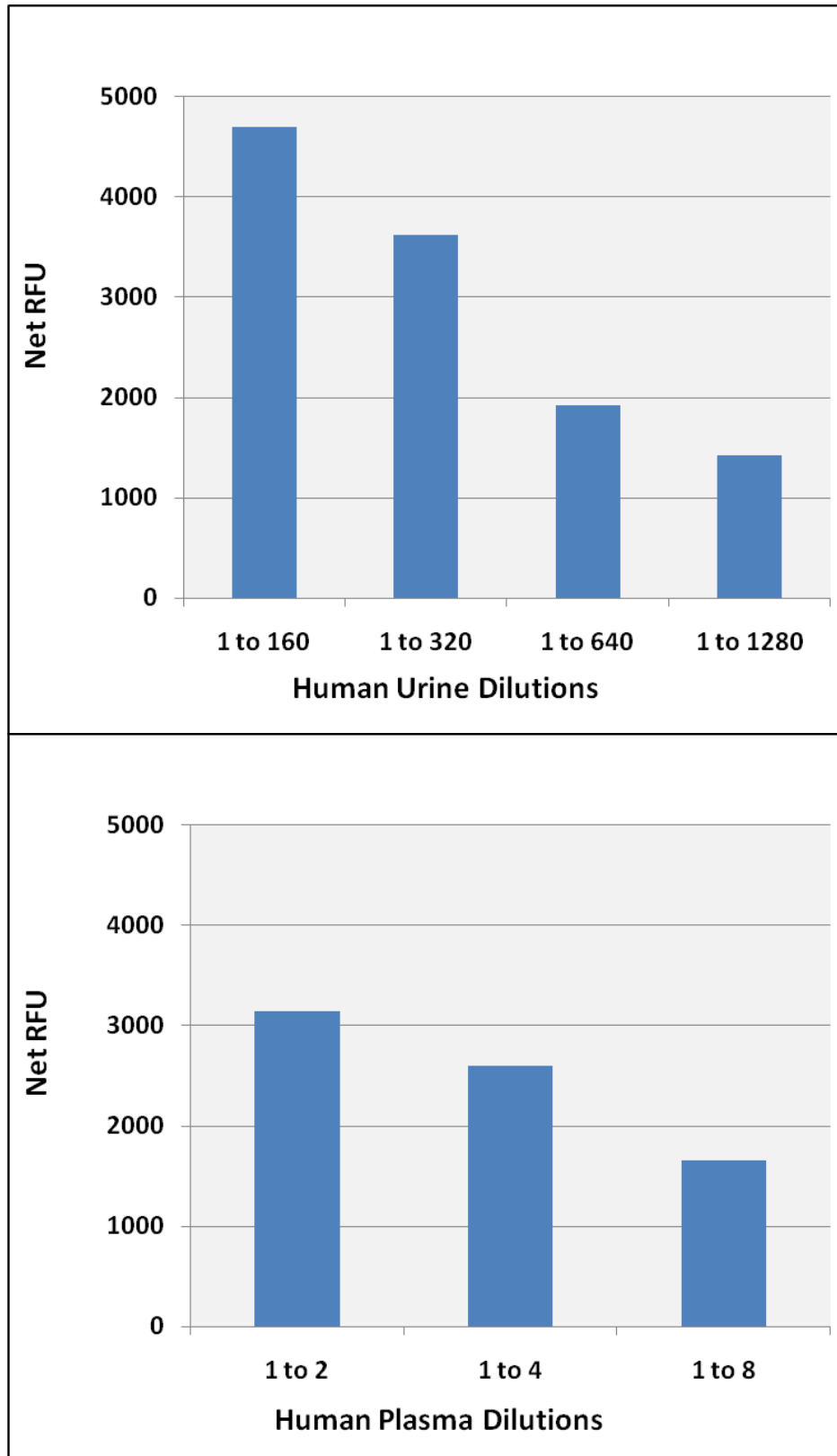


Figure 3: Urine and Plasma Samples. Human urine and plasma samples were tested with the Urea Assay Kit (Fluorometric).

Calculation of Results

1. Determine the average fluorescence values for each sample, control, and standard.
2. Subtract the average zero standard value from itself and all standard and sample values. This is the background correction.
3. Graph the standard curve (see Figure 2).
4. Subtract the sample well fluorescence values without urease (F_{-U}) from the sample well fluorescence values with urease (F_{+U}) to obtain the fluorescence difference. The (F_{-U}) sample value represents the ammonia background concentration within the sample, while the (F_{+U}) sample value is the combined urea and ammonia background concentration within the sample. The fluorescence difference (ΔF) is due to the urea concentration:

$$(\Delta F) = (F_{+U}) - (F_{-U})$$

5. Compare the fluorescence values of each sample to the standard curve to determine and extrapolate the quantity of urea present in the sample. Only use values within the range of the standard curve.

References

1. Friedman and Young. (2000) *Effects of Disease on Clinical Laboratory Tests*, 5th ed. AACC.
2. Sands, J.M. et al. (2009) *Semin Nephrol.* **29(3)**: 178-195.
3. Walker, V. (2009) *Diabetes, Obes and Metab.* **11(9)**: 823-835.
4. Young, D.S. et al. (1972) *Clin. Chem.* **18**: 1041-1303.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858 271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.