

## Product Information

### Urea Assay Kit

Catalog Number **MAK006**  
Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Urea is the major end product of nitrogen metabolism in most animals and is produced in a series of reactions in the liver called the urea cycle. In the urea cycle, ammonia is converted to urea, which is carried by blood to the kidneys for elimination from the body. High levels of urea in the blood may indicate renal failure. Urea levels may also be elevated in response to treatment with certain drugs such as corticosteroids or in response to decreased kidney filtration due to dehydration or congestive heart failure. Decreased blood urea levels can occur in response to liver disease or malnutrition.

In this assay, Urea concentration is determined by a coupled enzyme reaction, which results in a colorimetric (570 nm) product, proportional to the Urea present.

### Components

The kit is sufficient for 100 assays in 96 well plates.

|   |        |
|---|--------|
| Urea Assay Buffer<br>Catalog Number MAK006A             | 25 mL  |
| Peroxidase Substrate, in DMSO<br>Catalog Number MAK006B | 0.2 mL |
| Enzyme Mix<br>Catalog Number MAK006C                    | 1 vL   |
| Developer<br>Catalog Number MAK006D                     | 1 vL   |
| Converting Enzyme<br>Catalog Number MAK006E             | 1 vL   |
| Urea Standard, 100 mM<br>Catalog Number MAK006F         | 0.1 mL |

### Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate – It is recommended to use clear plates for colorimetric assays.
- Spectrophotometric multiwell plate reader

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Briefly centrifuge vials before opening. To maintain reagent integrity, avoid repeated freeze/thaw cycles. Keep vials tightly capped when not in use to minimize the uptake of  $\text{NH}_3$  from the air.

Urea Assay Buffer – Allow buffer to come to room temperature before use.

Peroxidase Substrate – Thaw at room temperature to melt solution prior to use. Aliquot and store protected from light and moisture at  $-20\text{ }^{\circ}\text{C}$ .

Enzyme Mix, Developer, and Converting Enzyme – Reconstitute each with 220  $\mu\text{L}$  of Urea Assay Buffer. Mix well by pipetting, then aliquot and store, protected from light, at  $-20\text{ }^{\circ}\text{C}$ . Use within 2 months of reconstitution and keep cold while in use.

### Storage/Stability

The kit is shipped on wet ice. Storage at  $-20\text{ }^{\circ}\text{C}$ , protected from light, is recommended.

### Procedure

All samples and standards should be run in duplicate.

### Urea Standards for Colorimetric Detection

Dilute 5  $\mu\text{L}$  of the 100 mM (100 nmole/ $\mu\text{L}$ ) Urea Standard Solution with 995  $\mu\text{L}$  of Urea Assay Buffer to prepare a 0.5 mM (0.5 nmole/ $\mu\text{L}$ ) standard solution. Add 0, 2, 4, 6, 8, and 10  $\mu\text{L}$  of the 0.5 mM Urea standard solution into a 96 well plate, generating 0 (blank), 1, 2, 3, 4, and 5 nmole/well standards. Add Urea Assay Buffer to each well to bring the volume to 50  $\mu\text{L}$ .

**Sample Preparation**

Tissue (20 mg) or cells ( $2 \times 10^6$ ) should be rapidly homogenized in 100  $\mu\text{L}$  of cold Urea Assay buffer. Centrifuge at  $13,000 \times g$  for 10 minutes at  $4^\circ\text{C}$  to remove insoluble material.

Serum and other liquid samples can be directly added to the wells.

Bring samples to a final volume of 50  $\mu\text{L}$  with Urea Assay Buffer.

For unknown samples, it is suggested to test several sample dilutions to ensure the readings are within the linear range of the standard curve.

Ammonium ion,  $\text{NAD}^+/\text{NADP}^+$ , and pyruvate present in the sample can generate background. To control for background, include a blank sample for each sample by omitting the Converting Enzyme in the Reaction Mix.

**Assay Reaction**

1. Set up the Reaction Mixes according to the scheme in Table 1. 50  $\mu\text{L}$  of the appropriate Reaction Mix is required for each reaction (well).

**Table 1.**

Reaction Mix

| Reagent              | Sample Blank     | Samples and Standards |
|----------------------|------------------|-----------------------|
| Urea Assay Buffer    | 44 $\mu\text{L}$ | 42 $\mu\text{L}$      |
| Peroxidase Substrate | 2 $\mu\text{L}$  | 2 $\mu\text{L}$       |
| Enzyme Mix           | 2 $\mu\text{L}$  | 2 $\mu\text{L}$       |
| Developer            | 2 $\mu\text{L}$  | 2 $\mu\text{L}$       |
| Converting Enzyme    | –                | 2 $\mu\text{L}$       |

2. Add 50  $\mu\text{L}$  of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting, and incubate the reaction for 60 minutes at  $37^\circ\text{C}$ . Protect the plate from light during the incubation.

3. Measure the absorbance at 570 nm ( $A_{570}$ ).

**Results****Calculations**

The background for the assays is the value obtained for the 0 (blank) Urea Standard. Correct for the background by subtracting the 0 (blank) value from all readings. Background values can be significant and must be subtracted from all readings. Use the values obtained from the appropriate Urea standards to plot a standard curve.

**Note:** A new standard curve must be set up each time the assay is run.

Subtract the blank sample value from the sample reading to obtain the corrected measurement. Using the corrected measurement, the amount of Urea present in the sample may be determined from the standard curve.

**Concentration of Urea**

$$S_a/S_v = C$$

$S_a$  = Amount of Urea in unknown sample (nmole) from standard curve

$S_v$  = Sample volume ( $\mu\text{L}$ ) added into the wells

$C$  = Concentration of Urea in sample

Urea molecular weight: 60.07 g/mole

**Sample Calculation**

Amount of Urea ( $S_a$ ) = 4.84 nmole  
(from standard curve)

Sample volume ( $S_v$ ) = 50  $\mu\text{L}$

Concentration of Urea in sample

$$4.84 \text{ nmole}/50 \mu\text{L} = 0.0968 \text{ nmole}/\mu\text{L}$$

$$0.0968 \text{ nmole}/\mu\text{L} \times 60.07 \text{ ng/nmole} = 5.81 \text{ ng}/\mu\text{L}$$

### Troubleshooting Guide

| Problem  | Possible Cause  | Suggested Solution   |
|--|---|--|
| Assay not working                              | Ice Cold Assay Buffer                                     | Assay Buffer must be at room temperature   |
|  | Omission of step in procedure                             | Refer and follow Technical Bulletin precisely  |
|  | Plate reader at incorrect wavelength                      | Check filter settings of instrument  |
|  | Type of 96 well plate used                                | For colorimetric assays, use clear plates  |
| Samples with erratic readings                  | Samples prepared in different buffer                      | Use the Assay Buffer provided or refer to Technical Bulletin for instructions              |
|  | Cell/Tissue culture samples were incompletely homogenized | Repeat the sample homogenization, increasing the length and extent of homogenization step. |
|  | Samples used after multiple freeze-thaw cycles            | Aliquot and freeze samples if samples will be used multiple times                          |
|  | Presence of interfering substance in the sample           | If possible, dilute sample further   |
|  | Use of old or inappropriately stored samples              | Use fresh samples and store correctly until use  |
| Lower/higher readings in samples and standards | Improperly thawed components                              | Thaw all components completely and mix gently before use                                   |
|  | Use of expired kit or improperly stored reagents          | Check the expiration date and store the components appropriately                           |
|  | Allowing the reagents to sit for extended times on ice    | Prepare fresh Master Reaction Mix before each use  |
|  | Incorrect incubation times or temperatures                | Refer to Technical Bulletin and verify correct incubation times and temperatures           |
|  | Incorrect volumes used                                    | Use calibrated pipettes and aliquot correctly  |
| Non-linear standard curve                      | Use of partially thawed components                        | Thaw and resuspend all components before preparing the reaction mix                        |
|  | Pipetting errors in preparation of standards              | Avoid pipetting small volumes  |
|  | Pipetting errors in the Reaction Mix                      | Prepare a Master Reaction Mix whenever possible  |
|  | Air bubbles formed in well                                | Pipette gently against the wall of the tubes   |
|  | Standard stock is at incorrect concentration              | Refer to the standard dilution instructions in the Technical Bulletin                      |
|  | Calculation errors  | Recheck calculations after referring to Technical Bulletin                                 |
|  | Substituting reagents from older kits/lots                | Use fresh components from the same kit   |
| Unanticipated results                          | Samples measured at incorrect wavelength                  | Check the equipment and filter settings  |
|  | Samples contain interfering substances                    | If possible, dilute sample further   |
|  | Sample readings above/below the linear range              | Concentrate or dilute samples so readings are in the linear range                          |

LS,MAM 10/12-1