

Urine Samples - DUR2 Protocol

Standards. Prepare 200 μL 100 mg/dL Premix by mixing 100 μL of the Standard (200 mg/dL) and 100 μL dH₂O. Dilute standards in 1.5-mL centrifuge tubes as described in the Table.

No	100 mg/dL Premix + H ₂ O	Urea (mg/dL)
1	100 μL + 0 μL	100
2	60 μL + 40 μL	60
3	30 μL + 70 μL	30
4	0 μL + 100 μL	0

Urine Samples. Dilute 50-fold in dH₂O (n=50). Add 20 μL of each urine sample to two separate wells in a 96-well plate (each sample requires a sample blank). Then, follow steps 2 & 3 on the kit datasheet for urea detection.

Calculation:

Subtract the blank value (#4) from the standard values and plot the ΔOD against standard concentrations. Determine the slope and calculate the urea concentration of the Samples as follows:

$$[\text{Urea}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{\text{Slope}} \times n \quad (\text{mg/dL})$$

where R_{SAMPLE} and R_{BLANK} are OD readings of the Sample and Sample Blank respectively. n is the sample dilution factor.