

# Urea Assay

A plate based colorimetric assay to measure the concentration of urea in a wide variety of samples. The assay sensitivity is 0.1-0.2 nmol of urea per well.

Urea is a waste product formed in the liver as a result of protein metabolism and is excreted by the kidneys. It is the primary means of nitrogen excretion in mammals. Thus, measuring urea levels can provide information about kidney function or other metabolic conditions. AkrivisBio's Urea Assay is a simple, sensitive way to measure the concentration of urea in a variety of biological samples. In this assay, urea is hydrolyzed by the urease to ammonia, which via a series of reactions results in the formation of hydrogen peroxide. Hydrogen peroxide is then utilized by the peroxidase to oxidize ADHP thereby resulting in the formation of intense color measured at OD 570 nm. The assay sensitivity is 0.1-0.2 nmol of urea per well.

## Ordering Information

[Order Online »](#)

|             |           |
|-------------|-----------|
| AKR-MA-0118 | 100 wells |
|-------------|-----------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application                      Colorimetric detection

Species Reactivity              Species independent

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)