

# Urea Assay Kit

Cell Biolabs' Urea Assay Kit is based on the Berthelot reaction. Urea is first degraded into ammonia and carbon dioxide, which further reacts with an alkaline developer to produce a blue-green colored product that can be measured with a standard spectrophotometric plate reader at an optical density between 580-630 nm. Each kit provides sufficient reagents to perform up to 192 assays, including blanks, urea standards and unknown samples.

## Ordering Information

[Order Online »](#)

CBI-STA-382	200 assays
-------------	------------

## Manuals, SDS & CofA

[View Online »](#)

- Sample urea concentrations are determined by comparison with the known urea standards, the standard is linear up to 200 mg/dL urea
- Measures urea levels within urine, serum, plasma, cell lysates, or tissue homogenates
- Provides sufficient reagents to perform up to 192 assays, including blanks, urea standards and unknown samples

## Handling & Storage

**Use/Stability** Upon receipt, prepare aliquots and store the Urea Standard and Urease at -20°C. Store the remaining kit components at 4°C until their expiration dates.

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Colorimetric detection

**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)