

# Ammonia Assay II

A plate based colorimetric assay to measure the concentration of ammonia in various sample types.

Ammonia is important both as a source of and a means of eliminating nitrogen in living systems. It is required for the synthesis of amino acids, as well as nucleotides and some lipids. Ammonia is typically formed through amino acid deamination. In the liver, ammonia is converted to the much less toxic urea and in birds and reptiles, into uric acid. AkrivisBio's Ammonia Assay II is a simple, sensitive, reliable assay with a sensitivity below 1 nmole per 100  $\mu$ l well. In the assay, ammonia is converted to indophenol, with a strong absorbance measured at 670 nm.

## Ordering Information

[Order Online »](#)

AKR-MA-0140	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**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)